Sorvall HT 6 User's Manual



Sorvall HT 6



Carefully read this manual before operating your instrument.

Information contained in this document is the property of Thermo Electron Corporation; it may not be duplicated or distribuited without the owner's authorization.

The validity of the guarantee is subject to the observation of the instructions and precautions described in this document.

MANUAL NUMBER

REV	ECR/ECN	DATE	DESCRIPTION	ВҮ
а	89003530	02/06	Initial release	

Title

Great Britain



WEEE Compliance. This products is required to comply with the European Union's Waste Electrical & Electronic Equipment (WEEE) Directive 2002/96EC. It is marked with the following symbol. Thermo Electron has contracted with one or more recycling/disposal companies in each EU Member State, and this products should be disposed of or recycling through them. Further information on Thermo Electron's compliance with these Directives, the recyclers in your country, and information on Thermo Electron products which may assist the detection of substances subject to the RoHS Directive are available at www.thermo.com/WEEERoHS

Deutschland



WEEE Konformittät. Dieses Produkt muss die EU Waste Electrical & Electronic Equipment (WEEE) Richtlinie 2002/96ECerfüllen. Das Produkt ist durch folgendes Symbol gekennzeichnet. Thermo Electron hat Vereinbarungen getroffen mit Verwertungs-/Entsorgungsanlagen in allen EU-Mitgliederstaaten und dieses Produkt muss durch diese Firmen widerverwetet oder entsorgt werden. Mehr Informationen über die Einhaltung dieser Anweisungen durch Thermo Electron, die Verwerter und Hinweise die Ihnen nützlich sein können, die Thermo Electron Produkte zu identizfizieren, die unter diese RoHS Anweisung fallen, finden Sie unter www.thermo.com/WEEERoHS

Italia



Conformità WEEE. Questo prodotto deve rispondere alla direttiva dell'Unione Europea 2002/96EC in merito ai Rifiuti degli Apparecchi Elettrici ed Elettronici (WEEE). È marcato col seguente simbolo. Thermo Electron ha stipulato contratti con una o diverse società di riciclaggio/smaltimento in ognuno degli Stati Membri Europei. Questo prodotto verrà smaltito o riciclato tramite queste medesime. Ulteriori informazioni sulla conformità di Thermo Electron con queste Direttive, l'elenco delle ditte di riciclaggio nel Vostro paese e informazioni sui prodotti Thermo che possono essere utili alla rilevazione di sostanze soggette alla Direttiva RoHS sono disponibili sul sito www.thermo.com/WEEERoHS





WEEE. Este producto requiere cumplir con la Directiva 2002/96EC WEEE (residuos de quipos electricos y electronicos). Este producto esta marcado con el simbolo WEEE, como se describe. En cada Estado Miembro de la Union Europea (EU), Thermo Electron ha contratado con diversas companias para disponer y/o re-ciclar los residuos de equipos electricos y electronicos. Existe information adictional acerca de la conformidad de Thermo Electron con esta Directiva, incluyendo los nombres de las diversas companias autorizadas para disponer/recyclar nuestros productos. Adicionalmente suministramos informacion acerca de los productos de Thermo Electron que puedes asistir en la deteccion de substancian mensionades en la directiva RoHS, explicado en www.thermo.com/WEEERoHS

France



Conformité WEEE. Ce produit doit être conforme à la directive européenne (2002/96EC) des Déchets d'Equipements Electriques et Electroniques (DEEE). Il est merqué par le symbole suivant. Thermo Electron s'est associé avec une ou plusieurs compagnies de recyclage dans chaque état membre de l'union européenne et ce produit devrait être collecté ou recyclé par celles-ci. Davantage d'informations sur la conformité de Thermo Electron à ces directives, les recycleurs dans votre pays et les informations sur les produits Thermo Electron qui peuvent aider le détection des substances sujettes à la directive RoHS sont disponibles sur www.thermo.com/WEEERoHS

Declaration de conformite	We, Thermo Electron Industries S.A.S., Z.I. d'Aze Bellitourne 53200 CHATEAU-GONTIER, France, declare that the instrument defined here below complies with the following European directives:	
Declaration of compliance	Nous, Thermo Electron Industries S.A.S., Z.I. d'Aze Bellitourne 53200 CHATEAU-GONTIER, France, déclarons que le matériel désigné ci- après est conforme aux directives européennes suivantes:	
Tedesco	Wir, Thermo Electron Industries S.A.S., Z.I. d'Aze Bellitourne 53200, CHATEAU-GONTIER, France, erklären dass das hier beschriebene Gerät sich in Übereinstimmung mit folgenden europäischen Richtlinien befindet:	
Declaración de conformidad	Nostros, Thermo Electron Industries S.A.S., Z.I. d'Aze Bellitourne 53200 CHATEAU-GONTIER, France, declaramos que el instrumento bajo descrito está fabricado según las especificaciones establecidas por las siguientes normas europeas:	
Declaraçao de conformidade	Nós, Thermo Electron Industries S.A.S., Z.I. d'Aze Bellitourne 53200 CHATEAU-GONTIER, France, declaramos que o instrumento aqui apresentado está de acordo com as normas européias abaixo listadas:	
Dichiarazione di conformità	Noi, Thermo Electron Industries S.A.S., Z.I. d'Aze Bellitourne 53200 CHATEAU-GONTIER, France, dichiariamo che lo strumento sotto descritto è conforme alle seguenti Direttive europee:	
	• Machinery: 98 / 37 / CEE	
	• Low voltage: 72 / 23 / CEE - 93 / 68 CEE	
	• EMC: 89 / 336 CEE - 92 / 31 CEE - 93 / 68 CEE	
	• Safety standard: EN 61010-1 - EN 61010-2-020	

Château-Gontier, 28th June 2005
Managing Director Christian Galiana

Guarantee Terms

Thermo guarantees that this unit is free from defects in materials and workmanship when it leaves the factory, and will replace or repair the unit if it proves defective in normal use or during service for a period of ONE YEAR from the delivery.

This guarantee is invalid if the unit is incorrectly used, poorly serviced or neglected, mis-used or accidentally damaged.

There is no explicit guarantee other than as stated above.

For further information, assistance or service

USA/Canada:

Thermo Electron Laboratory Equipment LLC 308 Ridgefield Ct. Asheville, NC 28806 Tel: 800-522-7746 Fax: 828-665-4330 email: info.labequipment@thermo.com

Website: www.thermo.com

Contents

Chapter 1	Presentation of Product	1-1
-	Description	1-1
	Reference Standards	1-1
	Relative Centrifugal Force	1-2
	Drive System	1-3
	Refrigeration System	1-3
	Dimensions and Weight	1-3
	Electrical Specifications	1-4
	Performance	1-4
	Accessories	1-6
	RS485 Communication Specification	1-10
	Mechanical Risks and Mechanical Safety	1-10
	Internal and External Markings of Centrifuge	1-11
	Responsibility	1-11
01		
Chapter 2	Shipment/Iransportation	
	Environmental Conditions for Transport	2-1
	Lifting and Transport Conditions	2-1
Chapter 3	Storage	3-1
Chanter /	Installation	/_1
Chapter 4	Environmental Conditions for Installation and Use	
	Lineadring Instructions	4-1
	Chpacking instructions	
	Electrical Requirements	
		·····4-)
		4-0
	Deter Instellation	4-/ / 7
		4-/
	Power Off Operation	4-10
Chapter 5	Using the product	5-1
	Balancing the Load	5-1
	Controls and Indicators	5-4
	Use	5-4
	Integration	5-13
	Refrigeration System	5-16
	Cautions	5-17
	Rotor and Accessory Precautions	5-18
	Machine Alarms	5-29

Operational Limitations	5-31
Expected Bucket Life	5-31
Corrosion	5-32

Chapter 6	Maintenance	6-1
•	Daily Cleaning	6-1
	Weekly Cleaning	6-1
	Periodic Preventive Maintenance	6-2
	Suggested Template for Annual Centrifuge Safety Inspection	6-2
	Rotor Removal	6-5
	Fuse Replacement	6-6
	Lid Opening and Manual Lid Unlocking	6-6
	1 0 0	

Chapter 7	Disposal of Product7	1-'	1
-----------	----------------------	-----	---

Chapter 1 Presentation of product

Description The HT 6 centrifuge is designed to separate substances of different density by centrifugal force.

Different rotors can be fitted to the drive in order to match the sample containers and the required performance. The maximum speed (and, by consequence, the resultant centrifugal force) is a function of the rotor in use.

A control system permits the speed, acceleration and braking rates, run time and temperature to be monitored and controlled.

The refrigeration is assured by a refrigerant group charged with a CFC-free fluid. The refrigerant fluid circulates around the bowl to achieve maximum efficiency.

The HT 6 is mainly dedicated to the following application fields:

- Large volume industrial processing
- Biochemistry
- Radio immuno assay

It is recommended that the user not process biological fluids known to contain prions. In cases of spillage or container breakage, it is impossible to neutralize the prions without subjecting the equipment to destructive temperature levels.

It is recommended that the operator inspect each sample, before and after centrifugation, to ensure that the sample and its container are in good condition.

This centrifuge must be used in accordance with established laboratory standards in the political entity where the centrifuge is installed.

Reference Standard

This product is supplied with a CE Declaration of Compliance. See the CE declaration for further information.

Relative Centrifugal Force

Suspended solids are separated by the Relative Centrifugal Force (RCF) applied to the sample. The effective force increases with the square of the speed of rotation and the distance from the axis of rotation.

The following formulae permit the calculation of primary parameters and of transformations relating to changes in primary parameters.

Note The value introduced for the radius can be adjusted to allow for position within the tube such as at a boundary. Maximum radii are quoted in the specifications tables (chapter 1.9). ▲

Configuration formule

Legend	
R = radius (in millimetres)	
N = speed (in r.p.m.) ÷ 1000	
RCF = gravitational acceleration 'g'	
M+ = add to memory	_
MR = memory recall	

Note To calculate actual results, press the keys on a pocket calculator in the order shown. ▲

Primary calculations	Key sequence (not valid for CASIO, HP)	
RCF (x g) = $1.118 \text{ R } \text{N}^2$	N x = x 1.118 x R =	
Speed (r.p.m.) = 946 $\sqrt{\frac{\text{RCF}}{2}}$	$RCF \div R = \sqrt{x 946} =$	
Y R		
Radius (mm) = <u>RCF</u>	$N x = x 1.118 = M + RCF \div MR =$	
1.118 N ²		
Transformations	Key sequence	
To determine actual 'g' achieved at a different speed:		
$RCF2 = RCF1 \qquad \left(\frac{N2}{N1}\right)^2$	$N2 \div N1 = x = x RCF1 =$	
To determine actual speed required to achieve a different 'g' at the same	ne radius:	
$N2 = N1 \sqrt{\frac{RCF2}{1}}$	$RCF2 \div RCF1 = \sqrt{x N1} =$	
√ RCF1		
To determine actual speed required to achieve the same 'g' at a different radius:		
N2 = N1 $\sqrt{\frac{R1}{}}$	$R1 \div R2 = \sqrt{x} N1 =$	
$^{ m V}$ R2		

Drive System The rotor is driven by a three phase asynchronous motor. The rotor is situated in the centre of an armour plated centrifugation bowl. A control system ensures the correct drive.

Refrigeration System

Information about the refrigeration system:

- Max pressure: 26 bars
- Type of refrigerant: R404A
- Total weight of refrigerant: 1.5 kg
- The system is CFC free and is without inflammable fluid.

Dimensions and Weight

HT 6	
Height (lid open)	1650/1670 mm
Height (lid closed)	915/935 mm
Width	798 mm
Depth	905 mm
Weight (uncrated)	346 kg
Weight (crated)	380 kg
Rquired surface area	1.2 m ² + safety zone

Electrical Specifications

HT 6			
Mains supply	230 V ±10%, 50 Hz		
	or 208/230 V, 60 Hz		
Fuses	Main pcb	T 5 x 20, 160 mA	26387014
	Supply pcb F1	T 5 x 20, 63 mA	86001807
	Supply pcb F2	T 5 x 20, 250 mA	34000207
Motor	2000 W at 4400 rpm		
Refrigeration	1.5 hp. (1100 W)		
Average power	3500 VA		
Maximum power	4500 VA		

Performance

HT 6		
Max. speed		4,250 rpm
Max. RCF		5,735 x g
Max. capacity		6 x 1000 ml
Speed	Range	500 to 4,250 rpm
	Step	10 rpm
	Accuracy	± 10 rpm
RCF (x g)	Range	80 to 5,735 x g
	Step	1 x g
	Précision	<0.5%
Timer	Range	0 to +99 h 59 min 59 sec + hold position
	Step	1 min / 1 sec
Energy	Max. density	1.2 g/ml
	Max. load on drive shaft	37 kg
	Max. energy	129050 J

Note The maximum rotational energy is not systematically that obtained with the maximum permissible load on the drive shaft.

HT 6		
Integrator	Range	0 to 999 x 10 ¹¹ m.rd ² . sec
	Step	1 m.rd ² . sec
Acceleration rates		10
Braking rates		10
Temperature	Range	-20 to +40 °C
	Step	1 °C
	Accuracy	±1.5 °C
Motor		Induction
Display		High visibility digital display, 240 x 128 pixels
Control		Microprocessor (16 bit)
Memory size		24 programs
Program protection		Individual password

Winshielded rotor RP6 (Cat N° 11178252)							
Bucket	Cat N°	Max speed/RCF	Bucket	Load	Load w	eight	
		with lid	weight (kg)	density	kg/buck	(et*	
Round Cylindrical	11178136	4250 rpm	-	1.375	1.2	1.5	
		5735 x g	-				
Round Sealed	11178135	4250 rpm	3800 rpm	1.375	1.2	1.5	
		5735 x g	4585 x g				
Rectangular	11178138	4250 rpm	3800 rpm	1.54	1.2	1.4	
Open		5392 x g	4310 x g				
Rectangular	11178140	4250 rpm	3800 rpm	1.54	1.2	1.4	
Sealed		5392 x g	4310 x g				

Rotor RHK4B (Cat N° 11178255)						
Bucket	Cat N°	Max speed/RCF	Max speed/RCF	Bucket	Load	Load weight
		with lid	without lid	weight (kg)	density	kg/bucket*
NHK4B	11174023	-	3300 rpm	3.6	1.2	3.7
RIA			2301 x g			

*At bucket maximum speed

Accessories Accessories have been marked with a maximum useable lifetime, that must be respected. For commentary concerning expected lifetime of accessories, please refer to section 5.13.

Only accessories specifically designed for use with the Sorvall HT 6 are authorized for use with this centrifuge.

The accessories used with the HT 6 are:

RP6 windshielded 6 place swing-out rotor:

Designed for multi purpose uses: blood banking (12 quin bags or quad bags), large volume processing (6 x 1000 ml), biochemistry (cylindrical or rectangular plastic inserts), and microtitre carriers.

RHK4B 2 place swing-out rotor:

Designed for RIA, usual RIA counter racks can be directly centrifuged with proper accessories. Maximum RCF is 3300 rpm (2301 x g).

Windshielded Rotor RP6 Rotor Lid (Cat.No. 11178119) and RP6 Rotor Body (Cat. No. 11178252)

Specifications		
Bucket Type	Rectangular	Round
Max. Speed	4,250 rpm	4,250 rpm
Max. RCF	5,392 xg	5,735 xg
Sealed Buckets	Cat. No. 11178140	Cat. No. 11178135
(Set of 6)	includes buckets with sealing cap	includes buckets with sealing cap
	& stainless steel microplate tray	
Unsealed Buckets (Set of 6)	Cat. No. 11178138	Cat. No. 11178136
Buckets Sealing Cap (Each)	Cat. No. 11178216	Cat. No. 25144058
Applications	Microtubes, clinical and blood tubes, conical	Microtubes, clinical and blood tubes, conical
	and flat-bottom tubes, and microplates	and flat-bottom tubes/bottles
Capacity Highlights (per Rotor)	Total Tubes x Volume (mL)	Total Tubes x Volume (mL)
	• 504 x 0.8	• 222 x 3-5
	• 240 x 1.5 Eppendorf tubes	• 180 x 6
	• 360 x 3-5	• 120 x 5-7 vacutainer tubes
	• 240 x 6	• 114 x 10-15 vacutainer tubes
	• 216 x 5-7	• 72 x 15 conical
	• 144 x 10 vacutainer tubes	• 42 x 50
	• 84 x 15 conical	• 30 x 50 conical
	 84 x 15 vacutainer tubes 	• 12 x 100
	• 60 x 25 flat bottom	• 6 x 250 flat bottom
	• 48 x 50 flat bottom	• 6 x 250 conical
	• 36 x 50 conical	• 6 x 500 flat bottom
Capacity Highlights (per Rotor)	Total Tubes x Volume (mL)	Total Tubes x Volume (mL)
	• 12 x 100	• 6 x 1000 flat bottom
	• 12 x 250 flat bottom	
	• 54 x 96-well / 18 x deepwell microplates	

See accessory tables for more details.

Rotor RHK4B (Cat. No. 11178255)

Specifications						
Bucket	Rectangular for Clinical Analyzer Racks & RIA					
Max. Speed						
Max. RCF	2,301 xg					
Bucket Set (Set of 2)	Cat. No. 11174023					
Applications	RIA tubes, clinical rack holders, microtitration plates, microtubes, and other tubes					
Capacity Highlights (per Rotor)	Clinical Racks (place x # of racks)	Total Tubes x Volume (mL)				
	2 x15 LKB racks	• 360 x 1.5 microtubes				
	2 x 3 NUCLEAR racks	• 432 x 3				
	2 x 4 PACKARD racks	• 300 x 6				
	2 x 3 BECKMAN racks	• 270 x 5				
	2 x 5 SMAC-TECHNICON racks	• 228 x 10				
	2 x 7 MICROMEDICS racks	• 180 x 15				
	2 x 6 Microtitration plates	• 50 x 66				
	2 x 8 OLYMPUS racks					

See accessory tables for more details.

RHK4B Accessories in 11174023 Buckets Cat. No. Description Capacity 11178801 2 x 15 LKB 11178804 2 x 3 NUCLEAR racks 11178805 2 x 4 PACKARD racks 11178806 2 x 3 **BECKMAN** racks SMAC-TECHNICON racks 11178807 2 x 5 11178814 2 x 7 **MICROMEDICS** racks 11178808 2 x 6 Microtitration plates 11178820 2 x 8 **OLYMPUS** racks

RHK4B Accessories in 11174023 Buckets and 11178801 Trays						
Cat. No.	Capacity	(3 inserts per bu	cket) x 2 buckets		Ø x Height	
	No. of	Tubes Per	Volume	Total	min/max (mm)	
	Inserts	Insert	(mL)	Tubes		
11179470	6	60	1. 5 microtube	360	11.5 - / 35	
11179476	6	72	3	432	11.5 - 55 / 100	
11179467	6	50	6	300	13.5 - 55 / 100	
11179468	6	45	5	270	13.5 - 65 / 110	
11179469	6	38	10	228	16.5 - 65 / 110	
11179471	6	30	15	180	18 - 65 / 110	
11179472	6	11	50	66	31 - 65 / 110	

RP6 Rotor Accessories for Round Buckets (Cat. No. 11178135, Sealed; and 11178136, Unsealed)

Cat. No.	Color	Capacity (Capacity (1 insert per bucket) x 6 buckets			Ø x Height	Max Radius
		No. of	Tubes Per	Volume	Total	min/max (mm)	RP6 (mm)
		Inserts	Insert	(mL)	Tubes		
11178208	Grey	6	37	3-5	222	12.7 - 50 / 107	266
11178209	Grey	6	37	3-5	222	Decanting	266
11178210	Light blue	6	30	6	180	13.5 - 50 / 107	266

Chapter 1

Presentation of product

11178211	Yellow	6	20	5 - 7 vac.	120	13.5 - 70 / 150	266
11178212	Red	6	19	10 - 15 vac.	114	17 - 70 / 150	266
11178304	White	6	12	15 conical	72	17 - 122 / -	266
11178305	White	6	12	15	72	16 - 75 / 125	266
11178306	Coral	6	12	20	72	22.2 - 75 / 148	271
11178307	Dark red	6	7	50	42	30 - 75 / 150	273
11178308	Blue	6	5	50 conical	30	30 - 115 / -	280
11178309	Blue	6	4	50	24	35 - 75 / 153	283
11178310	Dark Green	6	2	100	12	45 - 75 / 150	281
11178311	Turquoise	6	1	250 flat bottom	6	63 - 75 / 160	266
11178313	Black	6	1	250 conical	6	61 - 127 / -	281
85240076	Black	6	1	375	6	77 - 125 / 146	281
11178312	Black	6	1	500 flat bottom	6	70 - 127 / 171	281
11178063	Black	6	1	500 blood bottle	6	76 - 177	281
11178064	Black	6	1	500 blood bottle	6	83 - 177	283
85240075	Black	6	1	900	6	94 - 177	282
11178203	White	6	1	1000	6	-	282
11178102	White	6	1	1000 mL Nalgene	6	-	282
11178101	White	6	1	1000 mL Sanyo/MSE 6	<u>3</u> -	282	

RP6 Rotor Accessories for Round Buckets (Cat. No. 11178138, Unleased and 11178140, Sealed)

Cat. No.	Color	Capacity (2	inserts per buck	et) x 6 buckets		Ø x Height	Max Radius
		No. of	Tubes Per	Volume	Total	min/max (mm)	RP6 (mm)
		Inserts	Insert	(mL)	Tubes		
11175340	Black	12	42	0.8	504	8.5 - 35 / -	257
11175341	Salmon	12	20	1.5 microtube	240	11 - 35 / -	257
11175342	Grey	12	30	3-5	360	11 - 65 / 107	257
11175343	Beige	12	26	5	312	12.7 - 67 / 107	257
11175357	White	12	7	5 flat bottom	84	18 - 35 / -	257
11175344	Light blue	12	20	6	240	13.5 - 65 / 107	257
11175345	Yellow	12	18	5-7	216	13.5 - 65 / 108	257
11175347	Cream	12	12	5 - 10 flat bottom	144	17.5 - 65 / 112	257
11175346	Red	12	12	10 vac.	144	16.5 - 65 / 112	257
11175358	White	12	7	12 flat bottom	84	18 - 65 / 122	257
11175348	White	12	7	15 conical	84	18 - 90 / 122	259
11175349	Orange	12	7	15 vac.	84	16.5 - 90 / 127	259
11175350	Brown	12	4	15 vac.	48	16.5 - 90 / 133	259
11175351	Coral	12	8	20	96	22.2 - 90 / 116	259
11175352	Light yellow	12	5	25	60	25 - 90 / 120	259
11175375	Light yellow	12	5	25 flat bottom	60	25 - 90 / 120	259
11175376	Light blue	12	3	50 flat bottom	36	30 - 90 / 117	259
11175353	Dark red	12	4	50	48	30.5 - 90 / 122	259
11175354	Light blue	12	3	50 conical	36	30 - 90 / 117	259
11175361	Dark blue	12	2	50	24	35 - 90 / 122	259
11175355	Green	12	1	100	12	45 - 90 / 132	255
11175356	Turquoise	12	1	250 flat bottom	12	63 - 65 / 125	264
				Description			
85540163	-	-	-	Microplate adapter	(54 x 96-well	plates; 18 x deepwell plate	es)
11175359*	Black	12	1	Cytology Carrier for	cytobuckets (11174197 Set of 2 cytobuc	kets)
11175360	-	-		4 sets of 5 decantin	ig rubber pads	(Ø 10.5 - 12.1 mm) for 111	75343 inserts

* Set of 2 inserts

SORVALL HT 6 Centrifuge Specifications

Specifications	
Max Capacity	
Swing-out (RP6)	6000 mL = 6 x 1000 mL
Swing-out (RHK4B)	3300 mL = 66 x 50 mL
Max RCF	
Swing-out (RP6)	5735 xg
Swing-out (RHK4B)	2301 xg
Max Speed	
Swing-out (RP6)	4250 rpm
	Swing-out (RHK4B) 3300 rpm
Temperature range	-20°C to +40°C in 1°C steps with TFT™ system
Temperature condensation	-9°C to +9°C
RCF Range	80 to 5735 xg
Speed Range	500 to 4900 rpm
Timer	1 sec to 99 min 59 sec / 1 min to 99 hr 59 min + Hold
HRF (integral)	100 to 998 x 10E11
Acceleration Rates	10 rates
Braking Rates	10 rates
Control System	Microprocessor; self-diagnostics for servicing
Display	VIDEOset™ graphics (includes run graph); info/Error messages in two languages
Number of Memories	24, individual password
Drive	Direct
Motor	Brushless, induction
Dimensions (H x L x D)	94.5 x 78 x 90 cm (37.2 x 30.7 x 35.4")
Orderimg Information	
Cat. No.	Description
11178191	6L SORVALL HT 6 refrigerated centrifuge 208V/60Hz
11178255	RHK4B swinging bucket RIA rotor
11174023	RHK4B buckets (set of 2)
11178252	RP6 windshielded rotor body (no buckets)
11178119	RP6 windshielded rotor lid (required for 11178252)
11178140	RP6 sealed, rectangular buckets (set of 6)
11178138	RP6 unsealed, rectangular buckets (set of 6)
11178135	RP6 sealed, round buckets (set of 6)
11178136	RP6 unsealed, round buckets (set of 6)

RS485 Specification Communication

Type of transmission line	differential
Maximum number of drivers	32
Maximum number of receivers	31
Maximum cable length	1500 m
Maximum data rate	9.6 K octets/s
Maximum CMV	-7 V
Driver output	1,5-6 V
Driver load	60 Ω

Note For the physical characteristics relating to RS485 communication, please refer to the specifications described in the user manual (N° 89000821, chapter 3). ▲

Mechanical Risks and Mechanical Safety

Mechanical risks under normal use conditions are as follows:

Cause	Effect	Design Safeguards
Bucket that ruptures or escapes	Internal explosion due to impact of	Lid interlock system
from rotor	buckets and rotor on the bowl	 Armored chamber around bowl
		Highly stable chassis
Rupture of rotor or trunnion pin	Internal explosion due to impact of	Lid interlock system
	buckets and rotor on the bowl	 Armored chamber around bowl
		Highly stable chassis
Broken drive shaft or drive system	Internal explosion due to impact of	Lid interlock system
component	buckets and rotor on the bowl	 Armored chamber around bowl
		Highly stable chassis
Rupture or leakage of a sample during	Imbalance that causes severe vibrations,	Automatic imbalance shutdown
centrifugation	and in the worst case, a crash of the	Lid interlock system
	rotor and buckets	 Armored chamber around bowl
		Highly stable chassis
Insufficient dynamic balancing of	Imbalance that causes severe vibrations,	Lid interlock system
samples	and in the worst case, a crash of the	 Armored chamber around bowl
Automatic imbalance shutdown	rotor and buckets	Highly stable chassis
Loss of seal or rupture of the	Pressurized stream of liquid shooting out	The refrigeration system is enclosed by
refrigeration system or one of its	of refrigeration system	the external panels of the centrifuge.
components		There is no risk of explosion in this case.

Internal and External Markings of Centrifuge

Note The following table lists and describes the notices and markings found on the interior or exterior of the machine. ▲

Type of notice or marking	Interior/exterior	Specific Location	Description
Product information label	Exterior	Rear panel	Name of manufacturer
			Reference and serial numbers
			Main characteristics
			of machine
Danger symbol	Exterior	Rear panel	Beware of electric shock
			(precaution: unplug the machine
			prior to performing any service).
Identification sticker	Exterior	Working surface next to bowl	Name of machine
		(visible when lid is open)	Reference and serial
			numbers
Sticker, ISO 9001 certification	Exterior	Front panel	Information concerning Sorvall's
			ISO certification
Environmental protection sticker	Exterior	Front panel	Guaranteed CFC-free: helps to
			preserve the earth's ozone layer
Danger symbol	Exterior	Right panel, next to opening	List of precautions that must be
		used for manual lid unlock	followed to ensure the safety of
		procedure	manual lid opening procedure
Other stickers (for the purpose	Interior	On certain component parts (not	Information and/or danger
of information or warning)		made by Sorvall)	potential related to these
			component parts
	Interior	Compressor	• Type of refrigerant
			• Type of oil

Responsibility

The user of this centrifuge is responsible to respect the recommendations made by the manufacturer, with respect to installation, use, maintenance, storage, hygiene, safety, lifetime limits, etc.

The manufacturer can be held responsible for the non-conformity of the product. For his part, the user can bear a part of the responsibility if he has not verified for himself that the product is in practical conformity, even if it conforms to administrative guidelines. The user is also held responsible if he does something himself to accentuate any possible problems with the product: the responsibility of the manufacturer can be reduced or eliminated taking into account all circumstances when dammages are caused jointly by product problems and by the fault of the victim or by the fault of someone for whom the victim is responsible. It is the responsibility of the user to train (according to the instructions in this manual) everyone who will use or service this equipment. It is also the responsibility of the user to verify the quality of each sample after centrifugation.

Chapter 2 Shipment/Transportation

Enviromental Conditions for Transport

Transport conditions must be in accordance with the following recommendations if the machine is to function properly once installed and used under normal operating conditions.

Note If required to do so, refer to local or national transportation standards. ▲

The machine is transported in an integrated packaging system of carton and protection, fixed to a wooden pallet.

Required environmental conditions for transport are as follows:

- Temperature: $-10^{\circ}C \le \le +50^{\circ}C$
- Humidity : $20\% \leq RH \leq 90\%$
- Light: no particular recommendation (machine is enclosed in opaque packaging)
- Protection from elements: the packaging may not be exposed to precipitation
- Restraints may be neccessary to prevent lateral movement of the package

Stacking is strictly prohibited.

Lifting and Transport

Due to the weight of the machine, all lifting and transporting must be carried out using proper handling equipment (eg : fork lift trolley) that complies with current regulations, and by people having undergone the necessary training.

Note The machine must be supported from underneath. If it has to be transported without its pallet, for example on a staircase, professional handling assistance is required. ▲

Chapter 3 Storage

Whatever the duration of storage, the following conditions apply. Environmental conditions:

- Temperature: $-10^{\circ}C \le \le +50^{\circ}C$
- Humidity: $20\% \leq RH \leq 90\%$
- Light: no particular recommendation (machine is enclosed in opaque packaging)
- Storage in a clean area that is not polluted by industrial activity.

Conditions related to proximity to other stored goods:

- Away from all inflammable or explosive goods
- Away from all corrosives or strong solvents that may be spilled or spread and that may for any reason come into contact with the packaging

Note The machine must be handled using all necessary precautions, using a forklift or pallet jack, by trained and authorized personnel. ▲

Chapter 4 Installation

Note The Sorvall HT 6 centrifuge must be installed by a Sorvall approved technical service provider. The user must fournish the Sorvall technical service provider with a source of electricity (and, when applicable, with a source of water) that is in strict conformity with local codes and regulations. After installation, the integrity of the source(s) should be periodically verified by the user. ▲

Environmental Conditions

This instrument is designed to operate safely under the following conditions:

- Indoor use.
- Temperature: 5° to 40°C.
- Maximum relative humidity of 80% for temperatures up to 22°C.
- Maximum altitude: 2000 m.

Maximum performance is assured across the following ambient temperature range: 15°C to 25°C.

Place the centrifuge and its accessories in a clean, non-corrosive environment to avoid damage.

Install the unit leaving at least 30 cm (12 in.) clearance between its back and any obstruction, and between two adjacent machines, to allow subsequent placement of a vessel to collect condensates.

If the centrifuge is installed next to other equipment or next to a walkway, leave a space of 65 cm between equipment and 85 cm between the equipment and the walkway.

Centrifuges should be placed preferably in a room specifically intended for their use and therefore containing no other apparatus.

Walls and partitions that are within a distance of 1 metre should be strong enough to stop the centrifuge should there be a major accident (mobile partitions, single-width chipboard and plasterboard partitions without reinforcement are not strong enough).

The floor must be completely flat and strong enough to support the weight (346 kg) of the machine without creating or maintaining any

vibrations. Vibrations may be caused by variations during centrifugation. See also chapter 4.4 for positioning the damping feet.

There are no set minimum specifications on the type of floor required as there are many technical factors to be taken into account, such as:

- Composition of the different layers of the floor,
- How the floor is joined to the wall,
- Structures supporting the floor,
- Positions and weights of loads already on the floor,
- Loads already existing on support structures.

The machines must be positioned so that they can be moved easily to enable regular access for maintenance.

This centrifuge should only be installed in a work area that is lighted to the requirements of the local labor code.

The IEC 1010-2-020 standard relative to laboratory centrifuges recommends the user:

• To mark out a clearance envelope 30 cm around the centrifuge or establish special procedures for the non entry of all persons or all dangerous materials into this space during the operation of the centrifuge.

Chapter 4 Installation

Unpacking Instructions



Figure 4.1 Unpacking the HT 6



Read the following instructions carefully before installation. During manipulation, never put your hand underneath the machine.

- Remove staples with a screwdriver.
- Remove both anchor bolts (A) using the 19 mm spanner/wrench supplied in the tool box.
- Remove the front anchor bolt (B).
- Using the prop (C) as a lever:
 - Remove the wooden beam and portable ramp (D) by lifting the machine at the back.
 - Remove the wooden beam and portable ramp (E) by lifting the machine at the right hand side, towards the front.
- Ensure that the ramps (F) are perpendicular to the pallet and in line with the wheels.

2 people are needed to unload the machine: 1 person at the front of the centrifuge to guide and brake the machine as it comes down from the pallet, a second person behind to guide the machine onto the ramps.

Electrical Requirement

The electrical supply circuit must conform to currently applicable regulations. The HT 6 centrifuge is delivered without an electrical plug ; this must be connected by a qualified engineer, trained by SORVALL, according to the following conductor colour code:

• Blue: neutral, brown: phase, yellow/green: ground

Depending on the equipment part number, the electrical supply should be:

- 230 V ±10 % 50 Hz 25 A max.
- 208 / 230 V 60 Hz 25 A max.

Remember that in order to respect the electrical safety standards related to protection of operators against indirect contact, the supply of power to the instrument must be via a power socket fitted with a protection device ensuring automatic cut-off in the case of an insulation fault.

A supply fitted with a cicuit breaker of the correct rating complies with this requirement. It is also recommended that different machines be connected to a circuit breaker that is situated outside the room.



It must be immediately possible to identify the plug of the machine, which must also be easy to reach by the user. \blacktriangle

Concerning the electrical supply to the centrifuge, there must be a circuit breaker outside the room where the centrifuge is installed that can be used to interupt the power supply to the centrifuge in case of emergency.

The IEC 1010-2-020 standard relative to laboratory centrifuges recommends the user:

• To provide an emergency stop switch enabling the mains power to be cut in the case of a malfunction. This switch should be placed at a safe distance from the centrifuge, preferably in a different room from that in which the centrifuge is situated.

Positioning

The 2 front feet

- Put the damping foot (4) under the screw of the foot (3).
- Clean the floor beneath the feet of the machine with a clean, dry cloth (or paper hand towel).
- Remove the protective film from the adhesive face of the pad.
- Centre the pad in line with the foot and stick it to the floor.
- Press firmly on the whole surface of the pad to ensure a good adherence.
- Using the tool (5) set the foot screw in position and turn the nut (2), until the castors are raised 0.5 to 1 cm above the ground.
- Then lock the screw by tightening the nut (1).



Figure 4.2 Positioning the HT 6

The rear foot

- Using the 17 mm box spanner (6) proceed as for the front feet.
 - Set the centrifuge straight on its feet using the spirit level provided with the machine.

Lid Opening - Manual Lid Unlocking

The HT 6 is equipped with a double locking safety device which stops the machine starting if the lid is not completely closed and prevents lid opening before the end of rotation.

When the centrifuge is switched on, wait until display returns to the stand-by mode then pull forwards the latch lever located on the right hand side of the unit : the lid is automatically unlocked and opens.



In the case of a mains power cut, opening of the lid is prevented by the lid locking safety device. It is recommended to wait for the mains to be switched back on so that this safety device enables the lid to be unlocked. \blacktriangle

During rotation, the slightest leak from a sample is enough to create aerosols. Rotation of the rotor can only be detected when the machine is switched on. Opening the lid manually if the rotor is still rotating, would disperse the aerosols in the environment.

Rotating parts are also a risk as these could come in contact with the user or be ejected. There is particularly high risk of injury if:

- The user tries to manually stop the rotor,
- Any object falls inside the centrifugation chamber.

Manual lid opening procedure



If the machine has stoppped because of a mains power cut, the rotor could still be rotating. Wait 10 minutes before opening the lid. In spite of the absence of noise, the rotor could still be rotating when you need to open the lid manually.

Manual lid unlocking must only be done by someone informed of the danger and of the necessary precautions to be undertaken.

Even in the case of a mains power cut, always set the mains switch to the OFF position before carrying out this operation.

- Insert the unlocking device into the hole on the top right of the instrument by piercing the protection sticker.
- While pressing down on the unlocking device, pull the lid lever forward.

• The lid opens upwards automatically.

The protection sticker must be replaced after it has been pierced (cat. n° 85241911: 230 V/50 Hz, cat. n° 85241912: 208 V/60 Hz).

It is recommended that manual lid opening operation be performed only by specially trained individuals who are authorized by the Lab Manager. The person responsible for the laboratory whenever the centrifuge is being used should strictly control the use of the manual lid unlocking tool.



Only use the manual lid opening procendre in case of power failure or failure of the electrical lid lock system, and then only with the main switch in the OFF position. \blacktriangle

Figure 4.4 Manual Lid Opening



Inspection

Before installation, the rotor should be thoroughly inspected for corrosion and cleanliness.

Chemical and stress corrosion will eventually lead to disruption of the rotor with potential severe damage to the centrifuge.

Particles stuck inside the pockets can cause breakage of tubes and lead to major imbalance and / or loss of sample and contamination.



Check that the central hole of the rotor and the drive spindle are clean and undamaged. Remove any particles with a soft cloth.

Refer to chapters 5.9 and 5.10.

Rotor Installation

The rotors and buckets used with the Sorvall HT 6 are very heavy. When handling these accessories, take appropriate precautions to prevent back or other skeleto-muscular injury.

Ensure that the rotor is attached properly to the drive shaft, and that

the windsheild lid is attached properly to the rotor. Any failure to do so may prevent the rotor/buckets from operating properly.



Rotor installation must only be undertaken by qualified and authorized personnel trained by SORVALL. Never lift a rotor by its windshield.

The lid is delivered separately from the rotor.

The rotor packing contains:

- a rotor,
- grease for the trunnions and shaft,
- a user's manual.

The lid packing contains:

- a rotor lid,
- a lid locking device,
- an Allen key,
- a special screw for rotor fastening,
- an instruction sheet, for assembling the lid locking device.
- Check cleanliness of the motor drive shaft and rotor hub.
- Apply a light film of grease to the motor shaft.
- Note the collar position on the shaft and on the rotor; two indicators on the rotor label sticker allow localization of the collar groove.
- Pick up the rotor by two opposite arms and place it on the shaft, ensuring shaft and hub alignment. Position the rotor gently, despite its weight. The rotor slides freely onto the drive shaft.
- Apply a light film of grease to the tip of the trunnions.
- Adjust the lid locking device as mentioned in the instructions.
- Position the rotor lid on rotor, verify that the lid locking device locks on to the rotor.

Windshielded RP 6 Rotors

The packing contains:

- a rotor,
- grease for trunnions and shaft,
- an instruction sheet.
- Check cleanliness of the motor drive shaft and rotor hub.
- Apply a light film of grease to the motor shaft.
- Note collar position on the shaft and the groove on the hub.
- Pick up the rotor by two opposite arms and place it on the drive shaft, ensuring shaft and hub alignment. Position the rotor gently, despite its weight. The rotor slides freely onto the drive shaft.
- Position the washer and the bolt and screw clockwise with the wrench supplied.



Never tighten the rotor bolt without the washer. \blacktriangle

• Apply a light film of grease to the tip of the trunnions.

All Rotors/Winshielded Rotors RP4 - RP6, With LID/RHK4B

Ensure that the accessories are correctly fitted and balanced.

- Check that the rotor is fully locked onto the drive shaft.
- Check that the rotor lid fittings are screwed on tightly.
- Check that there is no water nor other liquid in the base of the windshield.
- Check that all of the buckets are in place and that they swing freely.
- Check that there are no empty transfusion liners in the buckets.

Rotors P40 - P60 - C60 - HK4B Use only rotors and buckets specifically designed for the HT 6 (see chapter 1.8). Please contact us for information on P40, P60, C60 and HK4B rotors.

Power Off Operation

The full power off of the HT 6 centrifuge is obtained by disconnection of the power cord from the mains supply.



Immediately after disconnection and for a duration of not longer than a minute, voltage remains on centrifuge line cord connectors. Do not touch for at least a minute.

Chapter 5 Using the Product



It is essentiel to understand the precautions, risks and limitations of use to achieve the best performance and highest safety with the instrument and its accessories. \blacktriangle

Never centrifuge an empty liner because there is a risk of deformation.

Balancing The Load

When installing tubes into each bucket insert, be certain that the samples are distributed symmetrically. Swing-out rotors should be loaded with a full compliment of buckets to maintain correct balance and safety for the rotor.



Use only centrifuge-rated glass or plastic ware, tubes, vials or bottles. Although the HT 6 can compensate for an imbalance of more than 50 grams, sample containers should be balanced to avoid any long-term effects of vibration and assure good sample separation. If you intend to use old accessories on a new centrifuge, please contact your SORVALL representative for advice.



Never put liquid directly into the buckets.

Refer to the guide on the use of refrigerated centrifuges in blood transfusion.

Controls And Indicators

The mains switch is located on the lower left side of the front panel. The control system includes:

- A dot-matrix graphic LCD screen, able to display numbers, letters, and special characters in addition to animated graphics.
- Eight keys, surrounding 2 sides of the screen, have their functions determined according to the name displayed on the adjacent part of the screen.

Not all keys are necessarily operational at any one time and more than one key can operate the same function when the name of the function is displayed across a wide area of the screen next to these keys. • In addition, a standard numeric keypad (0 to 9 and CE, clear entry) enables parameter values to be keyed in directly when required.



Certain key functions are very common: others are found only on a dedicated screen. They operate as follows:

HT 6		
+	Increases the value of the indicated parameter. Sustaining pressure on the key causes a rapid increase.	
+/-	Changes the arithmetical sign of the current figure (from positive to negative or vice versa).	
-	Decreases the value of the indicated parameter. Sustaining pressure on the key causes a rapid decrease.	
\bigtriangledown	Moves the highlight bar down a list in order to make a selection.	
$\Delta \!$	Moves the s indicator from one parameter to another to allow values to be altered.	
ABOUT INTEGRAL	Shows a graphical representation of the integral, together with values calculated from the previous run	
	(where possible).	
	Enables the acceleration rate to be altered between the 10 choices from 0 (slowest) to 9 (fastest).	
BASIC DISPLAY	Switches the display from the fully detailed run screen to the basic run screen with 3 fundamental parameters.	
BRK	Enables the braking rate to be altered between the 10 choices from 0 (weakest) to 9 (strongest).	
CONFIG	Provides access to fundamental set-up choices, such as language selection.	
DETAILS	Gives access to those parameters associated with the indicated parameter.	
DISPLAY APPLI #	Displays all preselected values of the program whose basic parameters are on screen.	
END MODIF	Returns to the original run screen after temporary values have been entered.	
END PROG	Returns to the stand-by screen, keeping currently selected values in volatile memory ready for use but not in permanent memory for later recall	
ENTER	Beturns to a previous screen, memorizing all changes in parameter values	
ESC	Returns to a previous screen, without memorizing any changes in parameter values	
EXPAND DISPLAY	Switches the display from the basic run screen to the detailed run screen which displays all parameters	
	Selects continuous operation for the time parameter value	
ΜΔΧ	Directly chooses the highest possible speed, or "a" force taking into account the selected rotor and radius	
	(cannot operate without prior rotor selection).	
MODIF & SAVE	Enables parameter values to be changed and saved if the current application has not been already saved.	
MODIF(Y)	Enables parameter values to be changed (Values changed during a run other than PRE COOL will not be memorized).	
MORE INFO	Provides a more detailed explanation of the current parameter, e.g., definition of the integral value.	
NEW APPLI	Gives access to the selection of new parameter values starting from a set of default values.	
NONE	Selects "NO PASSWORD PROTECTION" when memorizing an application.	
PRE COOL	Starts a predetermined precooling cycle (2000 rpm, 4°C, 15 min). During this cycle all other programming facilities	
	are operational.	
RADIUS	Enables the effective radius of a selected rotor to be modified in order to allow for the use of an adaptor or to select	
	a boundary at which a set "g" force can be applied.	
ROTOR	Lists the available rotors from which a selection can be made with their capacity and performance.	
SAVE APPLI	Allows the current set of parameter values (complete or incomplete) to be stored in one of 32 memories, with or	
	without an individual password.	
SCROLL APPLI #	One press displays all parameter values of next used memory. Continued presses will display the values for all	
	memorised applications. Also acts as application (program) selection key.	
START	Starts the run	
STOP	Stops the run, overiding any programmed time or integral.	
Note If the instrument remains switched on, but not used, then 5 minutes after the end of the last run the screensaver operates. The backlight is extinguished although the display can still be read. Press any key to light up the screen then continue with normal operation. ▲

Use Introduction	The VIDEOset graphics based control software is designed to combine power with simplicity. VIDEOset includes on-screen guidance and intuitive programming.
	The screens presented in this manual could differ slightly according to the version of the software.
	Those who use this centrifuge must be well trained according to the instructions in this manual and according to the rules of the institution where the centrifuge is used. It is the responsibility of the purchaser to ensure that users are trained.
	This centrifuge should only be used in a work area that is lighted to the requirements of the local labor code.
Switching On	Immediately after switching on, the screen displays the SORVALL logo, then the stand-by screen is displayed showing the number of the last application program used and its basic parameters.
Precooling	A keypress on CE will cause the display of the value of the actual temperature of the air in the bowl to the left of the set value during the time that the CE key is pressed.
	If the operator wishes to precool the bowl and rotor, he simply selects PRE COOL having installed the rotor, and the HT 6 runs according to the pre-set values which are indicated on the stand-by display.
	The display now indicates PRECOOL ON. The remaining precool time, real temperature and the dynamic rotation symbol are displayed.
	Uniquely, during a PRECOOL run, it is possible to access all programming facilities. The precool run can be halted at any time using the STOP keys

Standard Use

If the program which is currently displayed is the one desired, press a START key to run the program.



Otherwise, another program can be recalled either by selecting the program number directly on the keyboard or by pressing Detail Appli # which first shows all details of the current program, then Scroll Appli # (same key) to scroll through the memorized application programs whose parameters are thus indicated on the screen.

Sustained pressure on the key causes rapid scrolling.

Ot a stallast	3800 RPM APPLI# DETAIL	⊲ ⊙	3800	RPM 24	SCROLL APPLI#	⊲ ⊙
Standby Screen	4 °C PRECOOL: OFF CONFIG		4	°C 28 4534 530 10		
	00 : 15 HR/MN 4'C COOL		00 : 15	8 HB/MN (حات 2 تاخی		
	MODIF NEW START		MODIF N & SAVE AF	IEW PPLI	START	ļ

Once the required program is displayed, simply press a START key to run the program.



During a run, the basic screen displays the 3 basic parameters with both set and live values.

A keypress on Expand Display presents a detailed screen showing all possible run parameters and the values of those selected, both set and live. A keypress on Basic Display returns the basic screen.

Note: If the radius has not been defined by selecting a rotor (and possibly modifying its radius to allow for the use of an adaptor) the screen will not display the RCF.

It is impossible to call up a program or parameter from the keyboard.



Basic Screen During a Rundentailed Screen During a Run

Programming Of Parameters Basic parameters - speed, temperature, time

The values of the 3 basic parameters, usually speed, temperature and time, are quickly introduced after selecting New Applion the stand-by display to show the simple programming screen.

Speed is indicated by a flashing arrowhead \triangleright . Use the keypad to write a new speed value (if required).

Move to temperature using $\Delta \nabla$. Use the keypad to write a new temperature value. If required, use the +/- to change the sign of the value.

Move to time using $\Delta \nabla$. Use the keypad to write a new time value (if required). HOLD is achieved by entering a value of 0 min.

After making the required modification(s), either select End Prog to return to the stand-by screen, ready to run but without storing the values, or select Save Aplli (see 5.4.6) to store the values first.

Note If they had been programmed previously, RCF could replace speed and integral could replace time in the basic parameter menu.

Note In either case, the new values are held in a buffer memory (identified by - -) where they remain until the next time End Prog or New Appli is used. ▲

Advanced parameters

A) Speed related: RCF, acceleration, braking, rotor, radius

To program any of the speed related parameters, select DETAILS from the simple programming screen when the flashing cursor \triangleright indicates SPEED. A screen appears with the speed related parameters and the flashing arrowhead indicates RCF.

A value can be input via the keypad so long as a rotor has already been selected. (No selection of rotor makes it impossible for the software to control by "g" force since this is directly linear with the radius of the base of the sample tube from the centre of the drive axis - see 1.3)

To choose a rotor, select ROTOR and use ∇ to move the cursor down the list of rotors to that which will be used. If an adaptor will be used in the rotor select RADIUS and use the - & + keys to choose the effective radius (see accessory table, 1.9.) and select ENTER. During operation with - &+ the dynamic radius arrow moves and the related RCF is changed in real time. If the user wishes to modify the acceleration and/or braking rates, he selects ACC or BRK and uses the - & + keys to choose the rate(s) indicated on the dynamic graphic screen Select ENTER to return to the speed related parameter screen.

After selection of ROTOR (and maybe RADIUS, ACC, BRK) the operator has the option to select MAX. Without requiring any knowledge on the part of the operator, selection of MAX causes the instrument to program the top speed of the rotor and the related RCF allowing for actual radius.

After all the required parameter changes for this screen have been made, select ENTER to save the changes and return to the simple programming screen (which could indicate control by RCF).

B) Temperature related: Temperature compensation (ΔT) , negative temperature

On the simple programming screen, use $\Delta \nabla$ to move the flashing arrowhead s to indicate the temperature, then select DETAILS.

The flashing arrowhead indicates ΔT (For an explanation of ΔT see 5.8). Use the keypad to input a value and then (if required) the +/- key to change the mathematical sign of the value.

Alternatively, the user has the choice of directly selecting one of the most popular operating temperatures (4, 20 or 37°C) and/or using +/- to introduce a negative value.

If wished use $\Delta \nabla$ to first indicate temperature selection then select the temperature. Select ENTER to return to the simple programming screen.

C) Time related: Integral, hold

On the simple programming screen, use $\Delta \nabla$ to move the flashing arrowhead s to indicate the time then select DETAILS.

The flashing arrowhead indicates integral selection. For more information on integral select About integral and also More info for an accurate definition.

Select ESC/ESC to return to the time / integral screen. If this application has just been used, VIDEOset will indicate the recently totalled integral value.

Selection of ENTER will return the user to the simple programming screen but with control by integral instead of time, ensuring the highest possible run-to-run reproducibility (see 5.7)

	Note this will only be possible if a rotor has already been selected, the integral value being linked to the 'g' force which is calculated from the centrifugation radius of the rotor.
	Alternatively, use $\Delta \nabla$ to indicate time and enter a value or press HOLD for operation with no time limit. Selection of ENTER makes a return to the simple programming screen.
	To use the modified parameter changes without permanently storing them select End Prog.
	During a run controlled by the integrator, the preset value for time will indicate REPROducible, the integral value will count down and the timer will count up the corresponding time.
Saving an Application Pasword Protection	From the simple programming screen select Save Appli (if the flashing curser \triangleright is on temperature , it will be necessary to select $\Delta \nabla$ to make Save Appli reappear). The application save screen appears, proposing the next available empty memory. If required, use \triangleleft and \triangleright to review other memories whose application program parameter values are indicated at the right hand side of the screen.
	Select ENTER to save the application in the indicated file. The user is offered the opportunity to enter an individual password for this application (accepts 0 to 9999 inclusive, via the keypad, then select ENTER) or to select NONE for no password protection to save the application.
	If the user chooses to store his application in a used but unprotected memory, the old values will be replaced by the new values. If he chooses to store his application in a used but protected memory he must first enter the password to gain access, to the memory. In both cases he can password protect the new application.
	Note The Save Appli function will only operate if one or more parameter values have been modified. ▲
Overriding Pasword Protection Erasing A Program	The unit is delivered without any passwords installed. The use of a password prevents easy overwriting of a protected application program.
!	We suggest that you write down the password in case you forget it. \blacktriangle
	In the case of problems there is a special procedure to enable the protection to be overridden. This same procedure is used to clear one or more memories to eliminate unwanted applications.

After Save Appli has been selected, the file screen is displayed.

To eliminate the password protection and erase a program use the \triangleleft and \triangleright keys to move to the required memory (closed file icon) and select CE. Use \triangleleft / \triangleright and CE to clear any other memories. Moving the cursor back over selected memories will then show them to be EMPTY as indicated in the right hand side of the screen. Now either select ESC to return to the standby screen or ENTER to save the modified application in the indicated file memory. The user can decide whether or not to password protect the new protocol which can now be stored in this memory.

To Modify A Program A) From the stand-by screen

With an application program on the stand-by screen, select odif New.

Now alter program parameters exactly as in programming a New Appli.

To SAVE select Save Appli from the simple programming screen and continue as under 5.4.6. A new memory will be proposed. To save in the original memory (erasing original values) use \triangleleft or \triangleright to move the cursor and enter the password if requested.

B) During a run

From the basic or detailed run display, select MODIFY. The three main control parameters (speed or RCF/ temperature/time or integral) will be displayed together with the current set values. Use $\Delta \nabla$ to move the cursor \triangleright to the parameter to be modified and enter the value via the keypad.

Select End Modif to return to the original run screen. Values modified during a run cannot be stored (such a randomly modified run would be impossible to reproduce) but remain in the volatile memory until the end of the run.

Note It is not possible to modify the integral value during a run.

Configuration The HT 6 can be configured to operate in set ways. To set up the instrument, select CONFIG from the stand-by screen.

A) Language programming

Having selected CONFIG, the screen highlights LANGUAGE SELECTION then select ENTER.

Select 1 for ENGLISH or 2 for FRENCH. The stand-by screen will be displayed in the selected language.

B) Selection of refrigeration stand-by

Having selected CONFIG use \bigtriangledown to highlight STAND-BY REFRIGERATION ON/OFF then select ENTER.

Choose 1 for active refrigeration during stand-by and 2 for refrigeration OFF during stand-by.

C) CQCS option (see following page)

D) Timer configuration

Select 1 to program hours/minutes. Select 2 to program minutes/seconds.

E) Service

A special code, reserve

CQCS Option Your centrifuge can be equipped with the CQCS Centrifugation Quality Control System.

This system allows the traceability for centrifuges, the products and the operators and follows the historical use of each instrument. It is thus also considered as a veritable maintenance tool.

With versions 3.0 and later, it is possible to transfer the centrifugation run data from up to 32 centrifuges to a PC.

CQCS ensures traceability of samples by a system of twin bar code readers. This system includes two principal stages.

1st stage : Association Support / Sample

This stage concerns the identification of supports (liners) and samples (blood bags) using a bar code reader connected to the PC (see the BB-CQCS manual).

2nd stage : Association Support / Centrifuge

This stage concerns the identification of the supports (liners) during their insertion into the centrifuge using a bar code reader connected to the centrifuge. This stage is carried out after filling and balancing the supports, ready to load them.

To activate this option, select CONFIG from the stand-by screen. Select CQCS, using V to display it in inverse video and press ENTER.

Enter the centrifuge serial number in the HT 6.

To do so, select the CQCS option in the CONFIG menu, then press START.

On the operator code display, press the top right hand side key to configure the serial number.



Enter the first 4 numbers of the serial number on the digital keyboard and validate with ENTER. These first 4 numbers must stand between 1000 and 9000 non-inclusive (for example 3990).

Then enter the last 4 numbers of the serial number on the digital keyboard and validate with ENTER. These last 4 numbers must stand between 0 and 10000 non-inclusive (for example 2103).

Press ENTER to validate the complete serial number.

The serial number registered in this example is 39902103.



If necessary, enter the Operator Code via the keyboard (the ENTER or ESC to abort).



After the identification of the liners by the PC, identify the liners to be centrifuged by the HT 6 using the bar code reader on the centrifuge. Read the ID of the first support then press ENTER. Continue up to the last support (a maximum of 6 for CQCS V3.0).



Select START to start the run.



For instructions on operating the CQCS application, please refer to the CQCS user manual (cat n° . 89000821).

Integration

Centrifugal acceleration and duration

Setting centrifugation parameters consists of determining the time during which the sample is submitted to the selected centrifugal acceleration.

In practical terms, centrifugal acceleration is set in terms of motor rotation speed (rpm). A given rotation speed corresponds to centrifugal acceleration which is directly proportional to the centrifugation radius, in keeping with the following equation:

"Number of g" Acceleration	=	1.118	Х	R Centrifugation	Х	N ² Speed in thousands
				radius (mm)		of rpm

For example, a given separation may be described as obtainable by centrifuging the sample in a given machine (which determines the accessory, thus the centrifugation radius) for 10 minutes (duration) at $4000 \ge g$ (R.C.F.)

Theoretically, if the centrifuge instantly attains the set point speed and brakes instantly, and if the requested speed attained is constant, reproducibility of results is guaranteed from one centrifugation to the next.

The centrifugation curve is as follows:



Figure 5.2 - Ideal centrifugation integral curve

It is thus obviously equivalent to centrifuge a sample at 4000 g during 10 minutes as to centrifuge it at 2000 g for 20 minutes (or 8000 g for 5 minutes).

In other words, centrifugations yield the same results when the product: DURATION X CENTRIFUGAL ACCELERATION remains constant. This product corresponds to the area of the curve.



Figure 5.3 - Equivalent centrifugation integral curves

Integral of the curve

The integral of the curve is the value of its area. The area represents the total amount of separating force applied to the sample. Under the previous conditions, calculation of this integral is very simple. In reality, the centrifugation curve appears as follows:

Centrifugal acceleration



Figure 5.4 - Actual centrifugation integral curve

The integral of this curve is calculated by adding the elementary areas obtained by calculating the product (elementary time interval x acceleration applied during this time interval).

The curve of two successive centrifugations for which the same speed and duration are applied are not identical. Indeed, according to sample weight, temperature, atmospheric moisture in the chamber, mains

	voltage fluctuations and motor wear, the acceleration rate, as well as constant speed phase, may vary.
	When constant speed durations are very limited or very long, the previously described fluctuations may give rise to 5 to 10% variations in value of the integral. The quality of the resultant separations is sometimes considerably modified, as for example when preparing sub cellular fractions.
	Integration control allows suppression of the effect of the various potential fluctuations. Indeed, rather than programming a given duration of centrifugation, it is possible to set the desired value of the integral for a given speed or centrifugal acceleration level (R.C.F.). Upon attainment of that value, the centrifuge braking phase is started automatically.
	The centrifuge calculator always provides values of the centrifugation integral over time. Thus, the optimization of centrifugation and determination of integrals to be entered is considerably simplified.
Using the Integrator	During a series of runs using the timer as usual, at the run end select MODIFY, $\Delta \nabla$, $\Delta \nabla$ (timer), DETAILS. The instrument will display the value of the integral measured automatically during the run. Note this value together with the sample identities. When the samples from the various runs have been analysed, note those which had the best results and look up the recorded integral value.
	Reprogram the run memory by chosing MODIFY, $\Delta \nabla$, $\Delta \nabla$ (time), DETAILS then write the value of the integral, then save the program by chosing ENTER, SAVE PROG as usual. Now the centrifuge will automatically make small adjustments to the time (slightly longer for high loads, poor balancing etc.). In this way, the identical total separating force will be supplied to all samples.
	The only differences in results will be due to the samples, not to the centrifugation. The use of the integrator is strongly recommended for blood platelet separation.

Refrigeration System

The HT 6 is equipped with a refrigeration system which allows the rotor chamber temperature to be controlled at a lower level than the ambient temperature. The measured value of air temperature in the bowl is displayed and the temperature is controlled by the microprocessor which controls the cooling unit.

The operator can use the temperature compensation setting so that the sample temperature is more accurately controlled according to the user request by measuring the difference between the sample temperature obtained during the run and the set value.

The cooling unit of the HT 6 has either an air cooled condensor or a water cooled condensor.

Air cooled condensor

Accumulation of dust on the condensor fins will seriously reduce the refrigeration power. Every 3 months, disconnect the unit from mains, remove the front panel and clean the condensor with a vacuum cleaner and a brush.

Water cooled condensor

In this configuration, the unit is provided with a water pressostat which controls the input of fresh water. Cold water (10°C and less) as well as warm water (25°C and more) reduce the efficiency of the cooling group.

Determination of the temperature compensation value (ΔT)

To determine the temperature compensation value (ΔT) to be programmed, carry out a run as normal with $\Delta T = 0^{\circ}C$ (no compensation applied). At the end of the run, measure the sample tempereture with a thermometer.

If, for example, the set temperature was 4°C and the measured temperature is 7°C, then program the ΔT as 3°C for future runs. Leave the protocol temperature at 4°C. The centrifuge will control the air at 1°C. The air temperature will be measured independently (1°C) then the value of ΔT will be added (here, 3°C) to give the displayed temperature of the sample (4°C).

In the case where an elevated temperature has been selected and the measured temperature attained at the run end is below the set value (eg. set value 30°C, attained value 28°C) then enter a negative value for ΔT (here, $\Delta T = -2^{\circ}C$).

Caution

1 The following precautions must be observed by the operator:

- Ensure that the accessories are correctly fitted and balanced.
- Check that the rotor is fully locked onto the drive shaft.
- Check that the rotor lid fittings are screwed on tightly.
- Check that there is no water nor other liquid in the base of the windshield.
- Check that all of the buckets are in place and that they swing freely.
- Check that there are no empty transfusion liners in the buckets.
- Do not attempt to override the lid interlock system.
- Do not try to open the lid until display returns to the stand-by mode shown when machine is switched on.
- Use only a properly grounded mains power socket.

Special attention is required to the following:

- Installation of the unit: proper ventilation, levelling of the centrifuge.
- Accessory handling: rotor, buckets.
- Cleaning of the accessories, of the rotor chamber, lubrication of the trunnions.
- Centrifuge ware: shape, material and quality of the vessels must be in accordance with the performance of the centrifuge.
- Load balancing.
- Samples: cleaning of the accessories is particularly necessary when using corrosive products in the samples (saline, acids, bases).
- Ensure that the set point speed is not more than the maximum speed inscribed on either the rotor or the buckets.

Rotor and Accessory Precautions

Corrosion Information

SORVALL rotors made of aluminium alloy are designed to operate at their rated RCF for many years. With careful use they will resist corrosion, lessening the possibility of excessive imbalance, disruption and subsequent damage to the instrument.

The primary conditions for the initation of corrosion exist in every laboratory during daily use of the centrifuge. For this reason it is essential that due care and attention be paid to inspection and cleaning.

Chemical Corrosion

This corrosion is characterised by chemical reactions due to the existence of an electrolyte liquid on the surface of the item. If these substances are allowed to remain on the surface the corrosion will almost certainly occur. This produces first a discoloration of the anodisation then pitting of the metal. Aluminium is a metal which easily reacts with ionic solutions. Anodisation produces a protective layer on the surface of the rotor.

Acidic and alkaline solutions sustaining their pH level will create problems of corrosion in aluminium rotors. Chlorides, present in salts or even in skin contact with the rotor are among the most aggressive and harmful substances commonly found in the laboratory. The chemical products which are the origin of this corrosion do not necessarily originate from broken tubes - for example they could come from:

- Chemical vapours present in the laboratory which are dissolved in the residual humidity, in condensed water (refrigerated centrifuges) present at the base of the rotor pockets.
- Corrosive liquids originating from overfilled uncapped tubes (the liquid overflowing during centrifugation).
- Inserts, adaptors, racks, bottles whose exterior has been soiled by a chemical product or poorly rinsed after decontamination (with bleach, for example).

Note: If the products are very corrosive, simply rinsing is insufficient. Residual traces dissolve little by little in the humidity present in the bottom of the pocket, in the condensed water (such as occurs when putting an ambient temperature insert in a cold rotor). ▲

Beware: of the presence of solid particles beneath tubes, inserts, racks or adaptors. These particles are crushed by the centrifugal force and penetrate the protective, anodised layer of buckets and rotors, thus creating easy pathways for corrosion. ▲

Stress Corrosion

This term relates to the phenomenom of accelerated corrosion due to the effect of centrifugation when a corrosive chemical is in contact with the alloy. From the time when the aluminium alloy has been attacked by chemicals, stress corrosion begins to appear. As it is on a microscopic scale it is even more dangerous than a macroscopic scale since it is invisible.

During centrifugation, chemicals responsible for corrosion are also submitted to the tremendous "g" force which pushes them against the alloy. This close contact facilitates the chemical reaction which occurs much faster than in a static situation. Moreover, centrifugal force is very directional thus corrosion under stress creates, with a very small amount of corrosive product, straight microscopic fissures. Each centrifugation run makes the chemical migrate further and further.

A fissure, although it is microscopic, is a cut in the metal, breaking the cohesion of the material. As one weak link in a chain allows the chain to break, so the microfissures break the chain of resistance of the accessory to centrifugal force.

As accessories are designed with high levels of safety, rupture does not occur as soon as the first microfissures are produced.

Depending upon the location of the fissure, disruption may occur before it reaches the external surface of the accessory. The fissure creates a weakness which makes the accessory less and less resistant to mechanical fatigue.

The corrosion by a small amount of corrosive product does not disrupt the accessory but makes it mechanically weaker and weaker until it disrupts due to both centrifugal effort and number of cycles.

Because stress corrosion is largely invisible, it is essential that rotors are scrutinised regularly paying particular attention to susceptible parts such as the base of the pockets, the outer edges and the base of the rotor. If fissures are suspected, it is forbidden to use the rotor until it has been examined by a specialist.

Cleaning



Users must not use cleaning or decontamination methods different from those recommended without first having verified that these methods do not risk damaging the instrument or its accessories, and obtained prior authorization from the manufacturer.

Ideally, rotors should be washed after every use but at least weekly in warm water containing a few drops of mild non alkaline detergent (domestic washing up liquid is ideal) and EVERY TIME AFTER A

SPILLAGE HAS OCCURRED.

Do not forget to wash the core of the rotor which comes into contact with the drive spindle.

Each pocket of the rotor must be washed thoroughly using a small nylon brush.

DO NOT USE METAL WIRE BRUSHES.

Once the rotor is clean, rinse it in running water, preferably distilled. Dry the rotor with a soft absorbant non-woven cloth or tissue. Drying may be finished off with warm air jet (e.g. a hair dryer).

MAKE CERTAIN THAT THE POCKETS ARE WELL DRIED.

For swing-out rotors, be sure to clean the grease from the trunnions and replace it with a small amount of fresh grease. This will ensure that the buckets swing freely.

The vast majority of apparent imbalance problems arise from the failure of the user to clean and grease the trunnions and not from instrument error.

Pay particular attention that no deposit remains in the bottom of the pocket because the pressure of a flask or tube from above during centrifugation will significantly increase the chances of corrosion.

NEVER LEAVE A DAMP ROTOR ON A METAL SURFACE, PARTICULARLY STAINLESS STEEL because an electrochemical reaction could take place with the aluminium or magnesium in the rotor.

Contamination Hazard Our centrifuges are likely to be used in medical research where hazardous substances, including radioactive chemicals, are frequently found (see also chapter 5.16 : Biological risks).

ALWAYS USE THE APPROPRIATE DECONTAMINATION PROCEDURES WHERE THE ROTOR IS EXPOSED TO THESE CHEMICALS.

Dangerous substance: It is recommended that the user takes all the necessary precautions when centrifuging substances known for their toxicity or radioactivity or when they are contaminated by microbiological pathogens, notably material of Risk Group II as identified in the "Laboratory Biosafety Manual" edited by the World Health Organisation: Centrifugation of Risk 2 microorganisms and materials:

- 1) Centrifuge tubes and specimen containers for use in the centrifuge should be made of thick-walled glass or of plastic and should be inspected for defects before use.
- 2) Tubes and specimen containers should always be securely capped.
- 3) The interior of the bowl should be inspected daily for staining or soiling at the level of the rotor. If this is evident then the centrifugation protocols should be re-evaluated.
- Angle heads should not be used in microbiology except in special, high-speed centrifuges. With ordinary angle heads, some fluid, even from capped tubes, may be ejected because of the geometry of the machine.
- 5) Except in ultracentrifuges and with small prothrombin tubes, a space of at least 2 cm should be left between the level of the fluid and the rim of the centrifuge tube.

Examples of commonly used techniques are outlined below. The information is given as a guide only. It is the responsibility of the owner to use the most suitable procedure. The rotor should always be completely disassembled before being subjected to heat and after external chemical cleaning. Seals, tubes and plastic components should be decontaminated.

Disinfection

Alcohol (70% ethanol or isopropanol) applied for 10 minutes is ideal for bacteria and viruses. Autoclave rotors for 20 minutes at 120°C to destroy micro-organisms. Rotor lids must be disassembled from rotor bodies. 'O' rings, autoclaved separately, should be replaced when deformed.

ANY PART WHICH HAS BEEN SUBJECTED TO TEMPERATURES ABOVE 130°C MUST BE DISCARDED.

Hypochlorite "bleach" used at 0.1% concentration with 10 minutes immersion is effective against bacteria, spores and viruses but, as an oxidising agent, is corrosive to metal alloys and must be thoroughly rinsed off and dried. It should never be used if there is surface damage to the rotor. Note that the black colour on the surface of the rotor will be gradually washed out if the rotor is regularly autoclaved or bleached. This does not necessarily denote a degradation in the anodization.

Formalin (37% formaldehyde in water) in contact for 10 minutes has a similar effect to chlorine bleaches. Rotors should be thoroughly rinsed

under running water for 5 minutes to remove all traces of formalin then dried completely. NOTE: FORMALDEHYDE IS TOXIC.

Gluteraldehyde 2%, sold under many brand names such as Cidex and Glutarex, requires total immersion for 10 minutes to ensure sterility. Thorough rinsing and drying is essential to protect users.

CAUTION : GLUTERALDEHYDE BUILDS UP TO A TOXIC LEVEL IN THE FATTY TISSUES OF THE BODY.

Phenols are very corrosive and should never be used.

Radioactive decontamination

We recommend that all radioactive contamination be referred to your Radioactivity Safety Officer. Rotors may be decontaminated by a mixture of equal volumes of:

a) Distilled water,b) SDS diluted to 10%,c) Ethanol diluted to 70%.

The rotor should then be rinsed with ethanol followed by distilled water and then dried completely. SORVALL makes no claims as to the effectiveness of proprietary brands of decontaminating solutions.

Preventative Maintenance In the interests of prolonging metallic accessory life it is advisable to smear the pockets with a silicone based grease such as SILICONE GREASE 500 from KF or to use a LANOLIN based grease. SORVALL is able to supply cans of Anti Corrosion Spray, catalogue N° 11175399, whose propellant drives out water-based moisture and leaves a protective lanolin film on the surface.

Storage of Rotors It is strongly recommended that rotors not in use should be stored upside down on a non-metallic grid or other support which allows free movement of air. Any moisture contained in the pockets will then be eliminated by gravity and evaporation. If the rotors are to be stored for a long period then we advise that they are dried perfectly and greased with silicone grease.

High speed rotors are frequently preconditioned by storing in a cold room or refrigerator. This tends to advance the onset of corrosion. We recommend, therefore, that such rotors are stored (dismantled) sealed inside plastic bags from which the air has been evacuated.

Chemical Compatibility Table for Rotors, Tubes and Accessories

S = Satisfactory ΒN = Buna N AL = Aluminum D = Discoloration but OK = Cellulose Nitrate CAB = Cellulose acetate butyrate CN P = Pure chemical OKDL = Delrin KΥ = Kynar M = Moderate resistance NO = Noryl NY = Nylon = Polyallomer (= PPCO) PC = Polycarbonate U = Unsatisfactory PA PP PE = Polypropylene X = Explosion risk !!!= Polyethylene 0 = No informationPPCO = Polypropylene Copolymer PS = Polysulfone d = Less resistance if $T > 50^{\circ}C$ SS = Stainless steel TF = Teflon ΤZ = Tefzel ΤI = Titanium t = Unsatisfactory if T $>50^{\circ}$ C

= Viton A

VA

Chemical	Material AL	BN	I CAB	CN	DL	KY	NO	NY	PA PPCC	PC)	PE	PP	PS	SS	TF	ΤZ	TI	VA
Acetaldehyde	S	U	U	U	0	0	0	0	Μ	U	Μ	Μ	0	S	S	Μ	S	U
Acetamide	0	0	0	0	0	0	0	0	S	U	S	S	S	0	0	0	0	0
Acetic Acid (5%)	S	Μ	S	S	Μ	S	S	S	S	S	S	D	S	S	S	S	S	Μ
Acetic Acid (20 %)	S	0	0	0	0	0	0	0	S	Μ	S	S	S	0	0	0	0	0
Acetic Acid (60%)	S	U	U	U	U	S	S	Μ	S	U	Μ	D	S	S	S	S	S	U
Acetic Acid (80 %)	S	0	0	0	0	0	0	0	S	U	St	S	Μ	0	0	0	0	0
Acetic Acid (Glacial)	S	U	U	U	U	S	0	0	S	U	Μ	D	Μ	S	S	S	S	U
Acetic Anhydride	S	0	0	0	0	0	0	0	Sd	U	U	Sd	U	0	0	0	0	0
Acetone	S	U	U	U	Μ	Μ	0	U	S	U	S	Μ	U	S	S	Μ	S	U
Acetonitrile	0	0	0	0	0	0	0	0	Mt	U	S	Mt	U	0	0	0	0	0
Acetylene	S	0	0	0	0	0	0	0	S	0	S	S	U	0	0	0	0	0
Adipic Acid	0	0	0	0	0	0	0	0	S	S	St	S	S	0	0	0	0	0
Alanine	S	0	0	0	0	0	0	0	S	S	U	S	U	0	0	0	0	0
Allyl Alcohol	0	0	U	0	S	0	0	U	0	S	S	S	0	0	S	S	S	0
Aluminum Chloride	0	0	S	S	0	S	0	S	S	S	S	S	0	U	S	S	S	0
Aluminum Fluoride	0	0	0	0	0	S	0	S	S	U	S	S	0	0	S	0	S	0
Aluminum Hydroxide	S	0	0	0	0	0	0	0	S	Mt	Sd	S	S	0	0	0	0	0
Aluminum Nitrate	M	0	0	0	0	0	0	0	S	0	0	S	0	0	0	0	0	0
Aluminum Sulphate	S	0	0	0	0	0	0	0	Sd	0	S	Sd	0	0	0	0	0	0
Amino Acids	S	0	0	0	0	0	0	0	S	S	S	S	S	0	0	0	0	0
Ammonia	S	0	0	0	0	0	0	0	S	U	S	S	Sd	0	0	0	0	0
Ammonium Acetate	0	0	0	0	0	0	0	0	S	S	S	S	0	0	S	S	S	0
Ammonium Carbonate	S	U	S	S	0	S	0	S	S	U	S	S	S	S	S	S	S	0
Ammonium Chloride	M	0	0	0	0	0	0	0	St	0	St	St	0	0	0	0	0	0
Ammonium Hydroxide (10%)	0	S	U	0	0	0	0	S	D	U	S	D	S	S	S	S	S	S
Ammonium Hydroxide (conc)	0	U	U	0	0	0	0	S	D	U	S	D	0	S	S	S	S	U
Ammonium Oxalate	0	0	0	0	0	0	0	0	Sd	S	S	Sd	S	0	0	0	0	0
Ammonium Phosphate	U	0	0	0	0	0	0	0	S	Μ	S	S	0	0	0	0	0	0
Ammonium Sulphate	S	S	0	0	U	0	0	S	S	S	S	S	0	S	S	S	S	0
Ammonium Sulphide	0	0	0	0	0	S	0	0	S	U	0	S	0	0	S	0	0	0
n-Amyl Acetate	S	0	0	0	0	0	0	0	Sd	U	Sd	Sd	U	0	0	0	0	0
Amyl Alcohol	S	Μ	U	0	S	0	0	S	S	S	S	S	0	0	S	S	S	Μ
Amyl Chloride	S	0	0	0	0	0	0	0	U	U	U	U	U	0	0	0	0	0
Aniline	S	0	0	0	0	S	0	0	U	0	S	Μ	0	0	S	S	S	0

	Material	AL	BN	CAB	CN	DL	KY	NO	NY	PA	PC	PE	PP	PS	SS	TF	ΤZ	TI	VA	
	Chemical									PPCO)									
	Aqua Regia	U	U	U	0	U	0	0	0	U	U	U	U	0	0	S	S	S	Μ	
	Barium Chloride	U	0	0	0	0	0	0	0	S	0	S	S	0	0	0	0	0	0	
	Barium Hydroxide	U	0	0	0	0	0	0	0	S	0	S	S	0	0	0	0	0	0	-
	Barium Sulphate	S	0	0	0	0	0	0	0	S	0	St	S	0	0	0	0	0	0	-
	Benzaldehyde	S	0	0	0	0	0	0	0	Sd	Mt	S	Sd	Μ	0	0	0	0	0	
	Benzene	S	U	Р	0	Μ	0	0	S	U	U	U	U	U	S	S	S	S	S	
	Benzoic Acid, Sat	St	0	0	0	0	0	0	0	Sd	Sd	S	Sd	Μ	0	0	0	0	0	-
	Benzyl Acetate	0	0	0	0	0	0	0	0	Sd	Mt	S	Sd	U	0	0	0	0	0	-
	Benzyl Alcohol	S	U	U	0	Μ	0	0	U	U	U	U	U	0	0	S	S	S	S	-
	Boric Acid	U	0	0	0	0	0	0	0	S	U	Sd	S	U	0	0	0	0	0	-
	Bromine	U	0	0	0	0	0	0	0	U	Mt	U	U	U	0	0	0	0	0	-
	Bromobenzene	U	0	0	0	0	0	0	0	U	U	U	U	U	0	0	0	0	0	-
	Bromoform	U	0	0	0	0	0	0	0	U	U	U	U	U	0	0	0	0	0	-
	Butadiene	S	0	0	0	0	0	0	0	U	U	U	U	U	0	0	0	0	0	-
	Butane	S	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	-
	n-Butanol	S	S	U	0	S	0	0	U	0	Μ	S	S	Μ	0	S	S	S	S	-
	n-Butyl Acetate	S	0	0	0	0	0	0	0	S	U	S	S	U	0	0	0	0	0	-
	Butylene	S	0	0	0	0	0	0	0	0	0	S	0	0	0	0	0	0	0	-
	Butyl Chloride	0	0	0	0	0	0	0	0	U	U	U	U	U	0	0	0	0	0	-
	Butyric Acid	S	0	0	0	0	0	0	0	U	Mt	U	U	S	0	0	0	0	0	-
	Caesium Acetate	Μ	0	0	0	0	0	0	0	S	S	S	S	S	0	S	S	0	0	-
	Caesium Bromide	Μ	0	0	0	0	0	0	0	S	S	S	S	S	0	S	S	0	0	-
	Caesium Chloride	Μ	0	S	S	0	0	0	0	S	S	S	S	S	S	S	S	S	0	-
	Caesium Formiate	Μ	0	0	0	0	0	0	0	S	S	S	S	S	0	S	S	0	0	-
	Caesium lodide	Μ	0	0	0	0	0	0	0	S	S	S	S	S	0	S	S	0	0	-
	Caesium Sulphate	Μ	0	0	0	0	0	0	0	S	S	S	S	S	0	S	S	0	0	-
	Caesium Trifluoracetate	Μ	0	Μ	U	0	0	0	0	S	S	S	S	S	Μ	S	0	0	0	-
	Calcium Carbonate	U	0	0	0	0	0	0	0	S	0	St	S	0	0	0	0	0	0	-
	Calcium Chloride	Μ	S	S	0	S	0	0	S	S	Μ	0	D	S	S	S	0	S	S	-
	Calcium Hypochlorite	Μ	U	0	0	Μ	S	0	S	S	Μ	S	S	S	U	S	0	S	S	-
	Calcium Sulphate	Μ	0	0	0	0	0	0	0	S	0	St	S	0	0	0	0	0	0	-
	Carbazole	0	0	0	0	0	0	0	0	S	U	S	S	U	0	0	0	0	0	-
	Carbon Sulphide	S	0	0	0	0	0	0	0	U	U	U	U	U	0	0	0	0	0	-
	Carbon Tetrachloride	Х	U	S	S	Μ	S	0	S	U	U	U	U	S	Μ	S	S	S	S	-
	Cedarwood Essence	0	0	0	0	0	0	0	0	U	St	U	U	Μ	0	0	0	0	0	-
	Chlorine, dry	Μ	0	0	0	0	0	0	0	St	Sd	St	St	0	0	0	0	0	0	-
	Chlorine, moist	U	0	0	0	0	0	0	0	Mt	Sd	St	Mt	0	0	0	0	0	0	-
-	Chloroacetic Acid	U	0	0	0	0	0	0	0	Sd	Mt	S	Sd	U	0	0	0	0	0	-
	p-Chloroacetophenone	0	0	0	0	0	0	0	0	S	U	S	S	U	0	0	0	0	0	-
	Chlorobenzene	0	0	U	U	0	S	0	0	U	U	U	U	0	0	S	0	S	0	-
-	Chloroform	Х	U	Μ	S	Μ	S	0	U	U	U	U	U	U	S	S	Μ	S	S	-
-	Chlorosulphonic	Μ	0	0	0	0	0	0	0	Μ	0	Μ	Μ	0	0	0	0	0	0	-
	Chromic Acid (5 %)	U	0	0	0	0	0	0	0	S	Μ	S	S	U	0	0	0	0	0	-
-	Chromic Acid (10%)	Μ	U	U	U	U	S	S	0	S	Μ	S	S	U	U	S	S	S	S	-
-	Chromic Acid (50%)	U	U	U	S	U	S	0	0	D	U	S	S	U	U	S	S	Μ	S	-
-	Cinnamon Essence	0	0	0	0	0	0	0	0	U	St	U	U	Μ	0	0	0	0	0	-
_	Citric Acid (10%)	S	S	S	0	Μ	S	S	Μ	S	S	S	S	S	S	S	S	S	S	-
-	Copper Nitrate	U	0	0	0	0	0	0	0	S	0	S	S	0	0	0	0	0	0	-

	Material AL	BN	CAB	CN	DL	KY	NO	NY	PA	PC	PE	PP	PS	SS	TF	ΤZ	TI	VA
Chemical									PPCO)								
Copper Sulphate	U	0	0	0	0	0	0	0	S	0	St	S	0	0	0	0	0	0
Croesol	S	0	0	0	0	S	0	U	S	U	S	S	0	0	S	Μ	S	0
Cyclohexane	S	0	0	0	0	0	0	0	Mt	Sd	Mt	Mt	U	0	0	0	0	0
Cyclohexanol	S	0	U	0	0	0	0	S	S	Μ	S	S	0	0	S	0	S	0
Cyclohexanone	0	0	0	0	0	0	0	0	Mt	U	U	U	U	U	0	0	0	0
Cyclopentane	0	0	0	0	0	0	0	0	Mt	U	U	U	U	U	0	0	0	0
Decane	0	0	0	0	0	0	0	0	Mt	Mt	Mt	Mt	Sd	0	0	0	0	0
Dextran Sulphate	М	0	0	0	0	0	0	0	S	S	S	S	S	0	S	S	0	0
Diacetone	S	0	U	0	0	0	0	0	S	0	S	S	0	0	S	0	S	0
Diacetone Alcohol	S	0	0	0	0	0	0	0	S	0	S	S	0	0	0	0	0	0
o-Dichlorobenzene	0	0	0	0	0	0	0	0	Mt	U	Mt	Mt	U	0	0	0	0	0
p-Dichlorobenzene	0	0	0	0	0	0	0	0	Mt	U	Mt	Mt	U	0	0	0	0	0
Dichloroethane	0	U	U	U	S	0	S	S	U	U	U	U	0	0	S	S	S	S
Dichlorophenol	0	0	0	0	0	0	0	0	U	U	U	U	U	0	0	0	0	0
Diethylamine	S	0	0	0	0	0	0	0	St	U	U	St	Sd	0	0	0	0	0
Diethyl Benzene	0	0	0	0	0	0	0	0	U	Mt	U	U	U	0	0	0	0	0
Diethylene Glycol	S	S	S	S	S	S	0	U	S	S	S	S	S	0	S	S	S	S
Diethylene Glycol Ethyl Ether	0	0	0	0	0	0	0	0	S	Mt	S	S	Μ	0	0	0	0	0
Diethyl Ether	S	0	0	0	0	0	0	0	U	U	0	U	0	0	0	0	0	0
Diethyl Ketone	S	0	U	U	Μ	0	0	U	U	U	Μ	Μ	0	0	S	Μ	S	0
Dimethylacetamide	0	0	0	0	0	0	0	0	S	U	St	S	U	0	0	0	0	0
Dimethylformamide	S	0	0	0	0	0	0	0	S	U	S	S	0	0	S	Μ	S	0
Dimethylsulphoxide	S	0	0	0	0	0	0	0	S	U	0	S	0	S	S	Μ	S	0
Dioxane	S	U	U	0	Μ	S	0	0	Μ	U	Μ	Μ	0	0	S	S	S	U
Diphenyloxide	S	0	0	0	0	0	0	0	U	0	0	U	0	0	0	0	0	0
Dipropylene Glycol	0	0	0	0	0	0	0	0	S	Sd	S	S	S	0	0	0	0	0
Distilled Water	S	S	S	S	S	0	S	S	S	S	S	S	S	S	S	S	S	S
Ethanol (50%)	S	S	S	S	Μ	S	S	U	S	Μ	S	S	S	S	S	S	S	S
Ethanol (95%)	S	S	U	U	Μ	S	S	U	S	U	S	S	S	S	S	S	S	S
Ethyl Acetate	М	U	U	U	Μ	S	0	U	Μ	U	S	U	U	0	S	S	S	U
Ethyl Benzene	0	0	0	0	0	0	0	0	Mt	U	St	Mt	U	0	0	0	0	0
Ethyl Benzoate	0	0	0	0	0	0	0	0	Sd	Μ	S	Sd	U	0	0	0	0	0
Ethyl Butyrate	0	0	0	0	0	0	0	0	St	U	St	St	U	0	0	0	0	0
Ethyl Chloride	S	0	0	0	0	0	0	0	St	U	St	St	U	0	0	0	0	0
Ethylene Chloride	S	0	0	0	0	0	0	0	St	U	St	St	U	0	0	0	0	0
Ethylene Glycol	S	S	S	S	S	S	0	U	S	S	S	S	S	0	S	S	S	S
Ethylene Oxide	0	0	0	0	0	0	0	0	Μ	Mt	Μ	Μ	S	0	0	0	0	0
Ethyl Ether	S	0	U	U	0	0	0	0	Μ	U	Μ	М	0	0	S	Μ	S	0
Ethyl Lactate	0	0	0	0	0	0	0	0	0	S	St	S	S	Μ	0	0	0	0
Ethyl Malonate	0	0	0	0	0	0	0	0	0	S	Mt	S	S	Μ	0	0	0	0
Fatty Acids	S	0	0	0	0	0	0	0	Sd	Sd	Sd	Sd	S	0	0	0	0	0
Ferric Chloride	U	S	0	0	Μ	S	S	S	S	0	S	S	0	U	S	S	S	S
Ferric Nitrate	М	0	0	0	0	0	0	0	S	0	St	S	Μ	0	0	0	0	0
Ferric Sulphate	S	0	0	0	0	0	0	0	S	0	S	S	0	0	0	0	0	0
Ficoll Paque	Μ	0	0	0	0	0	0	0	S	S	S	S	S	0	S	S	0	0
Fluorine	S	0	0	0	0	0	0	0	Mt	Sd	St	Mt	U	0	0	0	0	0
Fluorhydric Acid (10%)	U	U	М	Μ	U	S	0	S	S	Μ	S	S	S	U	S	S	U	0
Eluorhydric Acid (50%)		U	U	U	IJ	S	0	0	S	IJ	S	S	M	U	S	S	U	M

Material	AL	BN	CAB	CN	DL	KY	NO	NY	PA	PC	PE	PP	PS	SS	TF	ΤZ	TI	VA
Chemical									PPCC									
Formaldehyde (20%)	S	0	0	0	0	0	0	0	S	Sd	S	S	Sd	0	0	0	0	0
Formaldehyde (40%)	M	Μ	0	S	0	S	S	S	S	S	S	D	S	S	S	S	S	S
Formaldehyde (50%)	S	0	0	0	0	0	0	0	S	Sd	S	S	Sd	0	0	0	0	0
Formic Acid (100%)	S	Μ	U	0	U	S	S	U	S	Μ	S	S	0	U	S	S	S	U
Freon TF	U	0	0	0	0	0	0	0	Sd	Sd	Sd	Sd	Sd	0	0	0	0	0
Fuel Oil	0	0	0	0	0	0	0	0	Sd	Sd	Mt	Sd	Sd	0	0	0	0	0
Glucose	S	0	0	0	0	0	0	0	S	0	S	S	0	0	0	0	0	0
Glutaraldehyde	0	0	0	0	0	0	0	0	Sd	Sd	S	Sd	Sd	0	0	0	0	0
Glycerine	S	0	0	0	0	0	0	0	S	S	S	S	S	0	0	0	0	0
Glycerol	S	0	0	S	0	S	S	0	S	S	S	S	S	S	S	S	S	0
Heptane	S	0	0	0	0	0	0	0	Μ	0	S	Μ	0	0	0	0	0	0
Hexane	S	0	0	0	0	0	0	0	S	0	Μ	S	0	0	0	0	0	0
Hydrazine	0	0	0	0	0	0	0	0	U	U	U	U	U	0	0	0	0	0
Hydrochloric Acid (5%)	U	0	0	0	0	0	0	0	S	S	S	S	S	0	0	0	0	0
Hydrochloric Acid (37 %)	U	0	0	0	0	0	0	0	Sd	U	S	Sd	S	0	0	0	0	0
Hydrochloric Acid (50%)	U	U	U	U	U	S	S	0	Μ	U	S	Μ	0	U	S	S	S	М
Hydrochloric Acid (conc).	U	U	U	0	U	0	0	0	S	U	0	S	0	U	S	S	S	0
Hydrofluoric Acid (10%)	U	U	Μ	Μ	U	S	0	S	S	Μ	S	S	S	U	S	S	U	0
Hydrofluoric Acid (50%)	U	U	U	U	U	S	0	0	S	U	S	S	Μ	U	S	S	U	М
Hydrogen Peroxide (3%)	S	Μ	S	S	S	0	S	S	S	S	S	D	S	S	S	S	S	S
Hydrogen Peroxide (100%)	S	U	S	S	U	0	S	0	S	S	S	D	S	S	S	S	U	М
lodine, Crystals	S	0	0	0	0	0	0	0	Mt	U	U	Mt	U	0	0	0	0	0
Isobutyl Alcohol	0	Μ	U	0	S	0	0	U	S	S	S	S	0	0	S	S	S	S
Isopropyl Alcohol	U	Μ	U	U	S	0	S	U	S	Μ	S	S	Μ	0	S	S	S	S
Isopropylbenzene	0	0	0	0	0	0	0	0	Mt	U	Mt	Mt	U	0	0	0	0	0
Kerosene	S	0	0	0	0	0	0	0	Sd	S	Mt	Sd	St	0	0	0	0	0
Lactic Acid (20%)	0	S	0	0	0	0	S	0	S	S	S	S	S	S	S	S	S	S
Lactic Acid (100%)	0	S	0	0	0	0	0	0	S	S	S	S	0	S	S	S	S	S
Lead Acetate (ag.)	U	0	0	0	0	0	0	0	S	S	St	S	0	0	0	0	0	0
Lemon Essence	U	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0
Lime, (conc).	Μ	0	0	0	0	0	0	0	S	U	S	S	S	0	0	0	0	0
Magnesium Chloride	M	0	0	0	0	0	0	0	S	S	S	S	S	0	S	S	0	0
Magnesium Hydroxide	U	0	U	0	0	S	S	0	S	U	S	S	0	0	S	0	S	0
Magnesium Nitrate	Μ	0	0	0	0	0	0	0	S	0	Sd	S	0	0	0	0	0	0
Magnesium Sulphate	S	0	0	0	0	0	0	0	S	0	St	S	0	0	0	0	0	0
Manganese Salts	Μ	0	S	0	0	0	0	0	S	0	S	S	0	0	S	0	S	0
Mercury	0	0	0	0	0	0	0	0	S	U	S	S	S	0	0	0	0	0
Methanol (100%)	S	S	U	U	Μ	S	0	U	S	Μ	S	D	S	S	S	S	S	U
Methyl Acetate	S	0	0	0	0	0	0	0	Μ	U	М	М	U	0	0	0	0	0
Methyl Alcohol (100%)	S	S	U	U	Μ	S	0	U	S	М	S	D	S	S	S	S	S	U
Methyl Butyl Ketone	0	0	0	0	0	0	0	0	U	0	0	0	0	0	0	0	0	0
Methyl Ethyl Ketone	S	U	U	U	M	M	0	U	S	U	S	S	U	0	S	M	S	U
Methyl Isobutyl Ketone	St	0	0	0	0	0	0	0	S		S	S	<u> </u>	0	0	0	0	0
Methyl Isopropyl Ketone	0	0	0	0	0	0	0	0		1	S			0	0	0	0	
Methylene Chloride	<u>х</u>		1		S	S	0	1	1	1	M	1	1	S	S	S	S	
Mineral Oil	0	0	0	0	0	0	0	0	Mt	<u> </u>	Mt	Mt	Mt	0	0	0	0	0
Nickel Chloride	<u></u>			<u> </u>	0		0		S		St	S	0	0	0		<u> </u>	0
Nickel Salts	M	2	<u>S</u>	0	0	0	0	5	S	S	S	S	0	S	<u>S</u>	2	5	S
	1 1 1	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0

Ma	terial AL	BN	CAB	CN	DL	KY	NO	NY	PA	PC	PE	PP	PS	SS	TF	ΤZ	TI	VA
Chemical									PPCO)								
Nickel Sulphate	U	0	0	0	0	0	0	0	S	S	St	S	0	0	0	0	0	0
Nitric Acid (10%)	Μ	U	S	S	U	S	S	Μ	D	S	S	S	S	D	S	S	S	S
Nitric Acid (20%)	U	0	0	0	0	0	0	0	S	Sd	S	S	St	0	0	0	0	0
Nitric Acid (50%)	М	U	Μ	Μ	U	S	S	Μ	D	Μ	Μ	Μ	0	D	S	S	S	S
Nitric Acid (95%)	М	U	U	0	U	0	0	U	Μ	U	U	Μ	U	S	S	S	S	S
Nitric Acid (conc.)	St	0	0	0	0	0	0	0	U	U	Mt	U	U	0	0	0	0	0
Nitrobenzene	0	0	0	0	0	0	0	0	U	U	U	U	U	0	0	0	0	0
Octane	0	0	0	0	0	0	0	0	S	Sd	S	S	Sd	0	0	0	0	0
Octyl Alcohol	S	0	0	0	0	0	0	0	0	0	S	0	0	0	0	0	0	0
Oleic Acid	S	U	S	S	S	S	0	S	S	S	S	S	S	S	S	S	S	Μ
Orange Essence	S	0	0	0	0	0	0	0	Mt	Μ	Mt	Mt	Μ	0	0	0	0	0
Oxalic Acid	Μ	Μ	S	S	0	S	0	S	S	S	S	S	S	S	S	S	Μ	S
Oxygenated Water (20%)	S	0	0	0	0	0	0	0	S	S	S	S	S	0	0	0	0	0
Oxygenated Water (50%)	S	0	0	0	0	0	0	0	S	S	S	S	S	0	0	0	0	0
Oxygenated Water (90%)	S	0	0	0	0	0	0	0	S	S	S	S	S	0	0	0	0	0
Ozone	0	0	0	0	0	0	0	0	Sd	Sd	S	Sd	S	0	0	0	0	0
Paraffin	S	0	0	0	0	0	0	0	S	0	S	S	0	0	0	0	0	0
Pentane	S	0	0	0	0	0	0	0	U	0	U	U	0	0	0	0	0	0
Perchlorethylene	S	0	0	0	0	0	0	0	U	U	U	U	U	0	0	0	0	0
Perchloric Acid	Х	0	0	0	0	0	0	0	St	U	St	St	U	0	0	0	0	0
Perchloric Acid (10%)	U	0	0	0	U	S	0	0	S	U	Μ	Μ	0	U	S	S	S	S
Perchloric Acid (70%)	Х	0	0	0	0	0	0	0	Μ	U	Μ	Μ	U	0	S	S	0	0
Petrol	S	0	0	0	0	0	0	0	St	Μ	St	St	Μ	0	0	0	0	0
Phenol (5%)	S	U	0	0	U	S	0	U	Μ	U	S	Μ	U	S	S	S	U	S
Phenol (50%)	U	0	0	0	0	0	0	0	U	U	U	U	U	0	S	Μ	0	0
Phenol, crystals	U	0	0	0	0	0	0	0	St	U	St	St	U	0	0	0	0	0
Phenol, liquid	U	0	0	0	0	0	0	0	U	U	U	U	U	0	0	0	0	0
Phenyl Ethyl Alcohol	0	0	U	0	0	0	0	S	S	0	S	S	0	0	S	0	S	0
Phosphoric Acid (10%)	0	Μ	S	S	U	S	S	0	S	S	S	S	S	S	S	S	0	S
Phosphoric Acid (conc.)	0	U	Μ	Μ	U	S	0	0	S	U	S	S	S	Μ	S	S	Μ	S
Picric Acid	S	0	0	0	0	0	0	0	U	U	U	U	U	0	0	0	0	0
Pine Oil	0	0	0	0	0	0	0	0	Sd	St	St	Sd	Μ	0	0	0	0	0
Potash, conc.	U	0	0	0	0	0	0	0	S	U	S	S	S	0	0	0	0	0
Potassium Bromide	U	0	0	0	0	0	0	0	S	S	S	S	0	0	0	0	0	0
Potassium Carbonate	Μ	0	S	S	0	S	S	S	S	U	S	S	0	S	S	S	S	0
Potassium Chlorate	М	0	S	S	0	S	S	S	S	S	S	S	0	S	S	0	S	0
Potassium Chloride	U	0	0	0	0	0	0	0	S	S	S	S	S	0	S	S	0	0
Potassium Hydroxide (5%)	U	Μ	S	Μ	U	0	0	S	S	U	S	S	S	S	S	S	Μ	S
Potassium Hydroxide (conc.)	U	Μ	U	U	U	0	0	0	S	U	S	S	0	S	S	S	U	Μ
Potassium Nitrate	S	0	0	0	0	0	0	0	S	0	S	S	0	0	0	0	0	0
Potassium Permanganate	S	0	0	0	0	0	0	0	S	S	S	S	D	0	S	S	0	0
Propane Gas	S	0	0	0	0	0	0	0	U	St	U	U	Μ	0	0	0	0	0
Propionic Acid	0	0	0	0	0	0	0	0	Sd	U	Μ	Sd	S	0	0	0	0	0
Propyl Alcohol	S	0	0	0	0	0	0	0	S	0	St	S	0	0	0	0	0	0
Propylene Glycol	S	0	0	0	0	0	0	0	S	Sd	S	S	S	0	0	0	0	0
Propylene Oxide	0	0	0	0	0	0	0	0	S	St	S	S	S	0	0	0	0	0
Pyridine	U	0	0	0	0	0	0	0	Μ	U	S	Μ	0	0	0	0	0	0
Resorcinol, Sat'd., Sol	0	0	0	0	0	0	0	0	S	Sd	S	S	U	0	0	0	0	0

Material	AL	BN	CAB	CN	DL	KY	NO	NY	PA	PC	PE	PP	PS	SS	TF	ΤZ	TI	VA
Chemical									PPCC)								
Rubidium Bromide	Μ	0	0	0	0	0	0	0	S	S	0	S	0	0	0	0	0	0
Saccharose	U	0	0	0	0	0	0	0	S	S	S	S	S	0	0	0	0	0
Salycylic Acid, Sat	0	0	0	0	0	0	0	0	S	Sd	S	S	S	0	0	0	0	0
Serum	S	0	0	0	0	0	0	0	S	S	S	S	S	0	S	S	0	0
Silver Acetate	0	0	0	0	0	0	0	0	S	Sd	S	S	S	0	0	0	0	0
Silver Nitrate	U	0	0	0	0	0	0	0	S	S	S	S	S	0	S	S	0	0
Sodium Acetate	S	0	0	0	0	0	0	0	S	Sd	S	S	S	0	0	0	0	0
Sodium Bisulphate	Μ	0	S	S	0	S	S	S	S	S	S	S	0	S	S	0	S	0
Sodium Borate	Μ	0	0	0	0	0	0	0	S	S	Sd	S	0	0	0	0	0	0
Sodium Bromide	U	0	0	0	0	0	0	0	S	S	0	S	0	0	0	0	0	0
Sodium Carbonate (2%)	Μ	S	S	S	S	S	S	S	S	S	S	D	0	S	S	S	S	S
Sodium Chloride (10%)	S	S	S	S	S	0	0	S	S	S	S	S	S	S	S	S	Μ	S
Sodium Chloride (Sat'd.)	S	S	0	0	S	0	0	S	S	0	S	S	0	S	S	S	S	S
Sodium Hydroxide (>1%)	U	Μ	S	S	U	S	0	S	S	U	S	S	S	S	S	S	S	S
Sodium Hydroxide (10%)	U	Μ	U	U	U	S	0	S	S	U	S	S	S	S	S	S	S	S
Sodium Hydroxide (conc.)	U	Μ	U	U	U	0	0	0	Μ	U	S	Μ	0	S	S	S	Μ	U
Sodium Hypochlorite (5%)	Μ	Μ	S	S	U	S	S	S	D	S	S	S	S	Μ	S	S	S	S
Sodium Iodide	Μ	0	0	0	0	0	0	0	S	S	0	S	0	0	0	0	0	0
Sodium Nitrate	S	0	0	0	0	0	0	0	S	0	Sd	S	0	0	0	0	0	0
Sodium Sulfate	S	0	0	0	0	0	0	0	Sd	0	Sd	Sd	0	0	0	0	0	0
Sodium Sulphide	S	S	S	0	0	S	0	S	S	U	S	S	0	S	S	S	Μ	S
Stearic Acid	S	0	0	0	0	0	0	0	S	Sd	S	S	S	0	0	0	0	0
Sulphuric Acid (10%)	Μ	U	S	S	U	S	S	S	S	М	S	S	S	U	S	S	S	S
Sulphuric Acid (20%)	U	0	0	0	0	0	0	0	Sd	Sd	S	Sd	S	0	0	0	0	0
Sulphuric Acid (50%)	U	U	U	U	U	S	S	U	S	S	S	S	S	U	S	S	M	S
Sulphuric Acid (conc.)	U		U	U	U	S	0	U	D		M	D		M	S	S	U	S
Sulphuric Anhydride, dry or moist	S	0	0	0	0	0	0	0	S	Sd	S	S	S	0	0	0	0	0
Tetrachloroethane	M	0	0	0	0	0	0	0	M	0	0	M	0	0	0	0	0	
Tetrachlorethylene	0	0	0	0	0	0	0	0		0	S	<u> </u>	0	0	0	0	0	
Tannic Acid	M	0	0	0	0	0	0	0	S	0	bZ	S	0	0	0	0	0	0
Tartaric Acid	M	0	0	0	0	0	0	0	S	bZ	S	S	S	0	0	0	0	0
Tetrahydrofuran	S	0	0	0	0	0	0	0	11	11				0	S	S	0	0
Thionyl Chloride	0	0	0	0	0	0	0	0		11				0	0	0	0	0
Toluene	S		<u>Р</u>	S	M	S	0							S	S	S	M	
Trichlorethylene	S	11	0	0	0	S	0	11	11					11	S	S	M	S
Trichloroacetic Acid		0	0	0	0	0	0	0	Mt	Mt	Mt	Mt	ک	0	0	0	0	0
1.2.4 - Trichlorobenzen	0	0	0	0	0	0	0	0	11	11	11	11		0	0	0	0	0
Trichloroethane	0 2		0	0	M	0	0	0					M	0	0	0	0	<u> </u>
Triethylamine	0	0	0	0	0	0	0	0	11	0	0			0	0	0	0	0
	0	0	0	0	0	0	0	0	0	64 62	0	0	0	0	0	0	0	0
Tris Ruffer (neutral)	0	0	0	0	0	0	0	0	ى د	Su C	ა ი	ა ი	ა ი	0	0	0	0	0
Trisedium Decembero	0	0	0	0	0 M	0	0	0	ۍ د	0	ა 	ۍ د	0	0	<u>с</u>	ۍ د	0	<u> </u>
	0	0	<u> </u>	0		0	0	<u> </u>	<u>د</u>	0	ა 	ა 	0	0	ა ი	<u>о</u>	<u>ہ</u>	<u> </u>
Inton X-100	<u> </u>	0	0	0	0	0	0	0	2	2	2	2	3	0	<u> </u>		0	0
	0	0	0	0	0	0	0	0	50	JIVI	ა[50 50	U	0	0	0	0	0
	U	0	0	0	0	0	0	0	50	50	ət	50		0	0	0	0	<u> </u>
	IVI	0	5	5	5	0	0	5	5	5	5	5	5	5	5	5	5	0
	0	0	5	0	5	0	0	5	5	5	5	5	U	0	5	5	5	0
Vinylidene chloride	0	0	0	0	0	0	0	0	U	U	U	U	U	0	0	0	0	U

	Material	AL	BN	CAB	CN	DL	KY	NO	NY	PA	PC	PE	PP	PS	SS	TF	ΤZ	TI	VA
Chemical										PPCO)								
Xylene		S	U	Р	0	Μ	S	0	U	U	U	U	U	U	S	S	S	S	S
Zinc Chloride		Μ	S	S	0	0	0	0	S	S	S	S	S	S	Μ	S	S	S	S
Zinc Hydrosulphite		U	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0
Zinc Sulphate		U	0	0	0	0	0	0	0	Sd	0	S	Sd	0	0	0	0	0	0

* This table is intended as a guide only because of the difficulty in quantifying, cross-checking and monitoring the results under different conditions of temperature, pressure and purity relating to the solvents and samples dissolved therein. It is strongly recommended that you carry out your own trials, particularly before protracted work periods

Machine Alarms

N°	Message	Descritpion	Action
1	TACHOMETER FAILURE (error code 0)	Speed measured at value other than zero when the	Wait for the motor to stop and the lid to unlock,
		centrifuge is powered up.	before trying to restart the machine
2	TACHOMETER FAILURE (error code 1)	No signal from the tachometer	Contact Sorvall technical service
3	TACHOMETER FAILURE (error code 2)	Variation of tachometer signal is too sharp.	Contact Sorvall technical service
4	TACHOMETER FAILURE (error code 3)	Inconsistent tachometer signal	Contact Sorvall technical service
5	POWER FAILURE	Motor slip abnomally high	Contact Sorvall technical service
6	LID UNLOCKED	Attempt to start the centrifuge while the lid is open	Close the lid completely and try again to start
			the centrifuge. If the problem persists, contact
			Sorvall technical service
7	LID UNLOCKED	Warning, this is a major problem: The lid has been	Wait for the rotor to come to a complete stop. If
		opened while the rotor is in motion.	problem persists, contact Sorvall technical
			service.
8	IMBALANCE	Buckets/samples are not balanced.	Wait until the rotor has stopped and the lid has
			been unlocked. Rebalance the samples, then try
			to start the run again. If problem persists,
			contact Sorvall technical service.
9	MOTOR OVERTEMPERATURE	Motor temperature is too high	Contact Sorvall technical service
10	CHAMBER OVERTEMPERATURE	Bowl temperature is higher than 43°C	May be because the ambient temperature around
			the centrifuge is too high. Verify that the
			centrifuge is installed far enough away from heat
			sources and that there is enough space around
			the machine to allow proper ventilation. If
			problem persists, contact Sorvall technical
			service.
	COMMUNICATION FAILURE	(only if the centrifuge is in CQCS mode) The	Verify that the supervision application is active
		centrifuge has not received a request from the	on the supervisor PC.
		supervisor PC for over 2 minutes.	

Safety

At any time when the instrument is able to detect and inform the operator of a fault condition the screen will display a special message from among the following.

Imbalance: Important vibrations which can damage the accessories, the unit and the samples have been detected. The cause is the bad balancing of the inserts. As soon as the motor stops, open the centrifuge and

redistribute the samples to produce an equal weight on diametrically opposite sides of the rotor and, for swing-out rotors, on each side of the trunnion for each bucket.

Using the grease supplied ensure that the trunnions are regularly lubricated removing the old grease first with a clean tissue.

Note: Never mix buckets from different sets on the same rotor. If the message persists, call an engineer. ▲

LID unlocked: Displayed if the lid was not correctly shut. The instrument will inform you to close it correctly in order that it can start the requested run.

Tachometer failure: Displayed when an unrealistic speed situation is detected. \blacktriangle

For example, if the measured speeds in two sequential measurement steps are very different, or if a power cut occured or if the unit was switched off momentarily during a run, in which case speed would be detected immediately upon switch-on. If speed is readable (certainly true in the later case) its real value will be displayed and in every case a TIME OUT safety lock will be applied, preventing the user from opening the lid for a set period after the speed reaches zero, or at least effects this lock-out period in the case of no detected speed.

Motor overtemperature: When using the centrifuge in bad conditions (e.g. : motor failing) a sensor detects an overtemperature in the motor and the run is automatically stopped. ▲

Chamber overtemperature: When using the centrifuge in bad conditions or if the refrigeration system is failing, the chamber temperature can seriously increase depending on the speed level and the heat can damage the samples. At 45°C the run is automatically stopped. ▲

Other safety features are also included in the centrifuge. There is no corresponding message but these features ensure safe use of your centrifuge:

- overspeed safety
- current limitation
- armour plating

Operational	Caution: The HT 6 Centrifuge is not explosing proof. \blacktriangle
Limitations	Using this centrifuge with explosive samples is entirely at the user's own risk. The use of inflammable or explosive materials or materials which give a violent chemical reaction is strictly forbidden.
Expected bucket life Mechanical strength	Centrifuge buckets are manufactured from aluminium alloys of high mechanical strength and elasticity. During centrifugation their own mass and most importantly the load that they contain submit them to a stress load varying from some hundreds of kilograms to several tonnes. This stress varies according to the total mass and increases considerably with speed: it doubles with a speed increase of only 41%.
	It is therefore essential to rigorously respect the limits of speed and load indicated in the instructions and even on the accessories themselves. Exceeding these limits causes weakening followed by a total rupture of the metal.
Mechanical fatigue	The phenomenon
	This normal physical phenomenon appears when a mechanical force is applied and relaxed repeatedly. The accumulation of alternate application and releasing of force (as in the number of runs) translates after a given number of runs into a very rapid reduction in mechanical strength. The item cans no longer resist the same force as at the start of its life. If the force is applied repetitively with same intensity, it will rupture.
	It is this phenomenon of excessive fatigue which is used when repeatedly bending and unbending a piece of wire or plastic in order to break-it.
	Each centrifugation run is a fatigue run for the buckets.
	Bucket lifetime
	Rotors and buckets are marked with maximum lifetime ratings (in years and max number of cycles. In no case should these accessories ever be used past the lifetime for which they are rated. Ignoring the ratings exposes the user to serious danger
	Remarks: The graphs show that "soft spin" runs such as for obtaining PRP and for platelets from the buffy coat (up to 1500 x g or about 2200 rpm) do not incur any fatigue. So the number of runs carried out at low speed is not counted in the total for the calculation of the lifetime of the buckets.

Corrosion

The above graphs have been established from tests carried out on samples of metal in perfect condition and over a period of many hours.

As the useable lifetime of the buckets is several years, it is important to take into account the wear due to corrosion, to which the buckets may have been submitted, independently of the fatigue phenomenon.



Figure 5.5 - P6 Liners (11178112) In KR4.22



2100 runs/year = 7 runs/day, 300 days/year 1800 runs/year = 6 runs/day, 300 days/year 1500 runs/year = 5 runs/day, 300 days/year

Figure 5.6 - P4 LINERS (11178074) IN KR4.22

Electrical risks	High voltage is present behind the panels of the HT 6. It is not necessary for the user to intervene inside the machine. Only qualified techniciens are authorized to do this. Do not remove these panels unless the centrifuge has been first disconnected from the power outlet.
Improper use	• Never spin a rotor or its accessories at more than the maximum speed, indicated in chapter 3.
	• Do not attempt to override the lid interlock assembly.
	• Use only accessories designed for use in the HT 6.
	• Load the rotor only in the recommended way (see 5.1 and 5.2) and do not spin higher than the speed indicated on the rotor and accessories. As the centrifuge starts to spin, an improperly loaded rotor could cause enough force to damage the drive shaft and the rotor chamber.
Biological risks	By the action of the turbulence created in the bowl a centrifuge is an aerosol generator. In the case of products presenting a biological risk you must take suitable precautions to prevent or reduce this risk and also train operators in the use of this instrument and its accessories in the context of their applications.
	The following information should be given to all users including those persons using the equipment infrequently:
	• A ventilated centrifuge (without built-in refrigeration) permanently exhausts into the room air which has passed through its bowl while a refrigerated (or thermostated) centrifuge contains a relatively sealed bowl. Thus the risks from the spread of aerosols are minimised in a refrigerated centrifuge and significant in a ventilated centrifuge.
	• Aerosols are produced naturally when using uncapped tubes or bottles. They are also produced when a container deforms or breaks provoking a smaller or greater sample leak. We draw your attention to the use of tubes used in angle rotors. Even if the liquid level does not touch the cap when in a static state, it could do so during centrifugation and leak. You should therefore follow the



instructions of the tube supplier concerning the fill level and

sealing of the tube.

Once a biological risk is known or suspected, samples should be placed in a sealed container. Should such containers not be available, sealed accessories should be used such as buckets or angle rotors with sealed lids. Sealed rectangular buckets (11178140) are certified by CAMR -Porton Down for their containment efficiency. Unsealed buckets (11178138) are sealed if fitted with lid (11178216).

- Duration of aerosol presence : when aerosols are created in a centrifuge bowl they persist in the air for a period of 3-5 minutes after the rotor stops moving. Indeed, the action of opening the lid itself provokes the dispersion of aerosols in the environment.
- When an elevated risk of aerosols or of breakage is perceived the accessories should be handled using gloves and opened in an environment ensuring the protection of the operator (safety cabinet, glove chamber, wearing a mask ...), even if they are apparently sealed.
- Those who use or service this equipment must wear appropriate personal protective gear given the type of solutions being centrifuged.

Dangerous substances:

It is recommended tat the user takes all the necessary precautions when centrifuging substances known for their toxicity or radioactivity or when they are contaminated by microbiological pathogens (notably material of Risk Group II as identified in the "Biological Safety Manual for Laboratories" edited by the World Health Organisation). See also chapter 5.10.3.

Chapter 6 Maintenance



Caution: All cleaning should be done with the unit disconnected from mains.

Users must not use cleaning or decontamination methods different from those recommended without first having verified that these methods do not risk damaging the instrument or its accessories, and obtaining prior authorization from the manufacturer. Users must use the adaquate protection needed for the cleaning product used in each case.

Please note that no maintenance performed by the user requires the user to open the panels to expose the internal components of the centrifuge.

All service (other than routine service that can be performed by the user) must be performed by qualified technicians trained by Sorvall.

Daily Cleaning

Daily cleaning is not required, except in the event of glass breakage or a large amount of spillage in the rotor chamber or fairing. If you wish to clean the control panel:

- Turn off the centrifuge,
- Clean with a moist sponge and a soap product and rinse with a moist sponge. Dry with a lint-free cloth. To disinfect, if required, use 70% alcohol.

Weekly Cleaning

- a) In the event of spillage between the chamber and the windshield, clean the rotor chamber with a detergent and water only. First, apply a mild water/detergent solution to a clean rag. Then use the rag to remove all traces of the spill from the bowl of the centrifuge and from the outside of the windshield if necessary. Next, take a clean rag and apply the mild water/detergent solution to it. Reclean the surfaces of the bowl and the windshield with this rag. Next, take a dry rag and remove all traces of the solution from the surfaces.
 - b) The rotors and buckets should be washed with warm water (at < 50°C) and a mild soap or detergent (for example : a household detergent for handwashing).

Rinse the rotor thoroughly with (preferably distilled) water. Dry the rotor or bucket, and preferably with hot compressed air store upside down.

c) The HT 6 centrifuge is equipped with a drainage hose, attached to the rotor chamber to enable rinsing the chamber when required.

To rinse and drain, operate the unit at 3000 rpm, at 20°C temperature for 1 minute.

The following preventive maintenance procedures are required for the HT 6 centrifuge:

- Ensure that the positioning instructions, as described in chapter 2.5, are always carried out (horizontability, anti-skid pads, castor positioning).
- Check the condition of all accessories (rotors, buckets, inserts, sealed lids) see chapter 6.
- Check earth continuity protection (when undertaking any intervention inside the machine).

Stability

- Is the centrifuge level?
- Is the centrifuge stable?
- Is the centrifuge sitting firmly on the floor?
- Is the rotor firmly attached to the shaft?
- Is the seal between the bowl and motor attached?

Protective Elements

- Is the tachometer port plug in place?
- Are the lid lock latches in place?
- Are the lid lock eyes in place?
- Is the imbalance detector in place?
- Is the motor fixed firmly in place?

Periodic Preventive Maintenance

Suggested Template for Annual Centrifuge Safety Inspection

Visual verification of equipment

- Are the lid springs and hinges in place?
- Is the lid lock electrical system in place?
- When the lid is locked, is the bowl sealed?
- Are the external panels firmly attached?

Condition of materials

- Rotor
- Buckets
- Lid/bowl seal
- Bowl/motor seal
- Motor struts
- Lid latches
- Drive shaft
- Plastic bucket adapters
- Feet
- Dust on the condenser filter?
- Foreign objects in the bowl?
- Foreign objects in the winshield?
- Water in the bowl drain?

Condition of electrical connections

- Control panel attachments
- Attachments to power source, condensor, locks, switches, etc.
- Grounding and power cord

Functional verification of equipment

Presence and function of safety systems

- Lid lock switch
- Lid must not open while rotor is moving
- Machine must not start while lid is open
- Lid open detection system

Abnormalities during use?

- Noises
- Vibrations
- Temperature effects
- Abrupt impacts

Function of automatic/manual shutdown systems

- Imbalance shutdown
- Lid open shutdown
- Stop button shutdown
- Overtemperature shutdown
- Circuit breaker shutdown

Verification of adjustments

Condition of springs

- Lid lock springs
- Lid support springs

Recognition of abnormal play in mechanical systems

- Lid opening process
- Motor/bearings

Condition of pieces susceptible to wear

- Bearings
- Carriers, adapters, linersa
- Buckets
- Rotors

Adjustment of 'end of run' systems

- Lid lock mechanisms
- Lid open detector

Verification of condition of indicators

Condition of measurement systems

- Tachometer
- Other measurement systems (temperature, etc.)

Condition of displays

- Read-outs
- Indicator lights

Rotor Removal



Changing the type of rotor on the HT 6 must be carried out by qualified and authorised personnel, given the adjustment of overspeed control necessary for the type of rotor installed. Please consult your SORVALL supplier for any operations of this nature.

- Using the wrench supplied, remove the rotor mounting screw.
- Use the hub-puller supplied : place its screw above the top of the motor shaft.
- Tighten the puller screw clockwise until the rotor is pulled out completely (use wrench supplied if necessary).
• To reinstall the rotor, refer to section 4.7.

This operation must only be done by trained and qualified staff.

The HT 6 and buckets used with the Sorvall HT 6 are very heavy. When handling these accessories, take appropriate precautions to prevent back or other skeleto-muscular injury.

Fuse Replacement Changing of fuses must not be carried out by the operator but by a qualified engineer who will diagnose the fault before replacing the fuses.

Fuses:

Main pcb	T 5 x 20, 160 mA	26387014
Supply pcb F1	T 5 x 20, 63 mA	86001807
Supply pcb F2	T 5 x 20, 250 mA	34000207

LID Opening - Manual LID Unlocking

See chapter 4.5.

Chapter 7 Disposal of Product

The following disposal recommendations are valid for the machine (or its accessories) at the end of its useful life, or the parts that must be disposed of because they are worn out or degraded during use.

Prior to disposal, the machine and accessories must be decontaminated. In this respect the user is required to submit a decontamination certificate (Sorvall document number J84S, available upon request), completed and signed, to the company or persons who are in charge of disposal.

The elements of the product that may become noxious during the life of the product are the following:

- Electric batteries and accumulators situated in the front panel of the machine among the other electrical and electronic components,
- Oil in the compressor of the refrigeration system,
- Refrigerant.

Disposal of these elements must be in accordance with applicable regulations and recommendations defined by the relevant regulatory bodies. In addition, these elements must be disposed of differently based on the different noxious substances they contain (possibly by a specialized disposal company).

Certificate of Decontamination and Cleaning



Read carefully the instructions below before sending an instrument, or parts of it, to Thermo Service Dept. or to any Authorized Technical Assistance Service.

Mr / Mrs (name)	Establishment
Department	
Address	Post / Zip code
City	State / Country

Declares the cleaning and decontamination of the following:

Product type	Serial N°	
Rotor	Serial N°	
Rotor	Serial N°	
Accessory - Description	Serial N°	
Accessory - Description	Serial N°	

Nature of contamination

	•••••	 	 ••••••	••••••	••••••

Decontamination Procedure used

Decontamination certified by:

Mr / Mrs	Institution
Date	Signature

When an instrument, or parts of it, requires servicing by Thermo SAT or CSAT personnel, the following procedure must be accomplished to ensure personnel safety:

- Clean the instrument and proceed to its decontamination from any kind of dangerous products.
- Compile this Decontamination Certificate with all the information required.

• Attach this Certificate to the instrument (or part) before sending it to Thermo Service or other authorized technical service.

SAT and CSAT personnel will not accept to work on instruments deprived of this Decontamination Certificate. If an instrument is received at our Service facilities and, in our opinion, is a radioactive or biological hazard, the item will be refused and resent to the Customer. Disposition costs will be borne by the sender. Instructions for decontamination and cleaning are explained in the User Manual. Additional certificates are available from your local technical or Customer Service representative. In the event these certificates are not available, a written statement certifying that the instrument or part has been properly decontaminated and outlining the procedures used will be acceptable.

For your convenience, use a copy of this page