



Thermo Fisher Scientific

TV-860

Instruction Manual

53292-4

March 2010

© 2009 Thermo Fisher Scientific Inc. All rights reserved.

ULTRACRIMP® is either a registered trademark or a trademark of Thermo Fisher Scientific.

All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.

Thermo Fisher Scientific Inc. provides this document to its customers with a product purchase to use in the product operation. This document is copyright protected and any reproduction of the whole or any part of this document is strictly prohibited, except with the written authorization of Thermo Fisher Scientific Inc.

The contents of this document are subject to change without notice. All technical information in this document is for reference purposes only. System configurations and specifications in this document supersede all previous information received by the purchaser.

Thermo Fisher Scientific Inc. makes no representations that this document is complete, accurate or error-free and assumes no responsibility and will not be liable for any errors, omissions, damage or loss that might result from any use of this document, even if the information in the document is followed properly.

This document is not part of any sales contract between Thermo Fisher Scientific Inc. and a purchaser. This document shall in no way govern or modify any Terms and Conditions of Sale, which Terms and Conditions of Sale shall govern all conflicting information between the two documents.

Release history: 53292-4 printed in March 2010.

For Research Use Only. Not for use in diagnostic procedures.

Table of contents

	Thermo Scientific TV-860 Titanium Vertical Ultraspeed Centrifuge Rotor	iii
	Important Safety Information	iv
Chapter 1	DESCRIPTION	1-1
	Rotor Description	1-2
	Rotor Application	1-2
	Rotor Specifications	1-2
	Accessories	1-3
Chapter 2	RUN PREPARATION	2-1
	Prerun Safety Checks	2-2
	Rotor Holding Fixture	2-2
	Rotor Precool	2-3
	Chemical Compatibility	2-3
	Sample Preparation	2-3
Chapter 3	SPECIAL OPERATING CONSIDERATIONS	3-1
	Compartment Loads in Excess of Design Mass	3-2
	Precautions to Prevent Precipitation of Cesium Chloride	3-2
	Critical Speed	3-7
Chapter 4	OPERATION	4-1
	Rotor Loading and Sealing	4-2
	Rotor Balancing	4-5
	Rotor Installation	4-5
	Acceleration and Deceleration	4-6
	Tube Removal	4-6
	Tube Trimmer Accessory (for use with Optional Tubes)	4-7
	Rotor Log Book	4-7
Chapter 5	OPERATION	5-1
	Relative Centrifugal Force (RCF) Determination	5-2
	Calculation of Sedimentation Time in Aqueous (Non-gradient) Solutions	5-5
	Calculation of Sedimentation Time in Gradient Solutions	5-6
Chapter 6	DESCRIPTION	6-1
	Corrosion	6-2
	Cleaning	6-2
	Storage	6-3
	Inspection	6-3
	Overspeed Decal Replacement	6-3
	Service Decontamination Policy	6-4

Table of contents

Chemical Compatibility ChartA-1
Contact InformationB-1

This manual is a guide for the use of

Thermo Scientific TV-860 Titanium Vertical Ultraspeed Centrifuge Rotor

Data herein has been verified and is believed adequate for the intended use of the rotor. Because failure to follow the recommendations set forth in this manual could produce personal injury or property damage, always follow the recommendations set forth herein. Thermo Fisher Scientific does not guarantee results and assumes no obligation for the performance of rotors or other products that are not used in accordance with the instructions provided. This publication is not a license to operate under, nor a recommendation to infringe upon, any process patents.

Publications prior to the Issue Date of this manual may contain data in apparent conflict with that provided herein. Please consider all data in this manual to be the most current.

NOTES, CAUTIONS, and WARNINGS within the text of this manual are used to emphasize important and critical instructions.

WARNING informs the operator of a hazard or unsafe practice that could result in personal injury, affect the operator's health, or contaminate the environment.

CAUTION informs the operator of an unsafe practice that could result in damage of equipment.

NOTE highlights essential information.



CAUTION and WARNING are accompanied by a hazard symbol and appear near the information they correspond to.

© 1997, 19990, 2010 by Thermo Fisher Scientific

Important Safety Information

Certain potentially dangerous conditions are inherent to the use of all centrifuge rotors. To ensure safe operation of this rotor, anyone using it should be aware of all safe practices and take all precautions described below and throughout this manual.

WARNING

When using radioactive, toxic, or pathogenic materials, beware of all characteristics of the materials and the hazards associated with them in the event leakage occurs during centrifugation. If leakage does occur, neither the centrifuge nor the rotor can protect you from particles dispersed in the air. To protect yourself, we recommend additional precautions be taken to prevent exposure to these materials, for example, use of controlled ventilation or isolation areas.

Always be aware of the possibility of contamination when using radioactive, toxic, or pathogenic materials. Take all necessary precautions and use appropriate decontamination procedures if exposure occurs.

Never use any material capable of producing flammable or explosive vapors.

Never exceed the maximum rated speed of the installed rotor, to do so can cause rotor failure.

Always reduce (derate) rotor speed as instructed in this manual whenever:

- the rotor speed/temperature combination used exceeds the solubility of the gradient material and causes it to precipitate.
- the compartment load exceeds the maximum allowable compartment load specified. See Chapter 3, paragraph 3-1.

Failure to reduce rotor speed under these conditions can cause rotor failure.



CAUTION

Do not expose aluminum rotor components to: strong acids, bases, or laboratory detergents; liquid chlorine bleach; or salts (chlorides) of heavy metals such as cesium, lead, silver, or mercury. Use of these materials with aluminum can cause a chemical reaction that initiates corrosion.



Do not operate or precool the rotor at the critical speed, as this will have a detrimental effect on centrifuge component life. See Chapter , paragraph 3-3.

Do not operate the rotor unless it is symmetrically balanced as described in this manual. Operating the rotor out of balance can cause damage to the centrifuge drive assembly.

Always maintain the rotor in the recommended manner. The rotor and all accessories must be clean and inspected prior to each run; do not use rotors showing signs of corrosion or cracking. See Chapter 6, Care and Maintenance.

Do not autoclave or expose any aluminum rotor parts to temperatures in excess of 100°C.

DESCRIPTION

This manual contains information required to operate and maintain the Thermo Scientific TV-860 Vertical Ultraspeed Centrifuge Rotor. If you require additional information regarding operation or maintenance, please contact Thermo Fisher Scientific for assistance. In the United States, call Thermo Fisher Scientific toll-free 1-800-522-7746; outside the United States, contact the nearest Thermo Fisher Scientific office (see back cover) or your local representative for Thermo Scientific products. Thermo Scientific product information is available on our internet web site at <http://www.thermo.com/centrifuge>.

Contents

- “Rotor Description” on page 1-2
- “Rotor Application” on page 1-2
- “Rotor Specifications” on page 1-2
- “Accessories” on page 1-3

Rotor Description

The TV-860 Rotor assembly consists of eight sealing assemblies and an eight-compartment rotor body that is machined from a titanium forging for strength and corrosion resistance. The rotor has a black coating which provides surface protection and aids temperature regulation. The tube compartments are bored parallel to the rotor axis and, occasionally, are counterbored to achieve precision balance. A disc with alternative black and reflective segments is attached to the bottom of the rotor; this disc will prevent the rotor from being operated at a speed above the recommended maximum speed of the rotor.

Rotor Application

The TV-860 Rotor is used in an ultracentrifuge, primarily for density gradient separations; the design of the vertical rotor allows these separations to be done in shorter times than the conventional fixed-angle or swinging-bucket rotor. Specifically, the TV-860 can be used for many isopycnic and rate zonal applications. Proper operation of the TV-860 Rotor requires use of the automatic slow acceleration and deceleration controls that are standard on all Thermo Fisher Scientific Ultracentrifuges. Without the use of this feature, the gradient will orient and reorient too fast—the slow acceleration and slow deceleration prevents disturbance of the gradient.

Rotor Specifications

Table 1-1. Rotor Specifications

Rotor Type	Vertical
Maximum Speed *	60 000 rpm **
Relative Centrifugal Force (RCF) at Maximum Speed:	
at r_{maximum} , 8.47 cm	340 596
at r_{average} , 7.195 cm	289 325
at r_{minimum} , 5.92 cm	238 055
Number of Tube Compartments	8
ULTRACRIMP® Tube Size	25 mm x 97 mm
Tube Compartment Diameter	2.54 cm (1 in)
Tube Compartment Length (Nominal)	8.89 cm (3.5 in)
Tube Volume (Nominal)	35 ml
Total Rotor Capacity (Nominal)	280 ml
Critical Speed	700 rpm
Maximum Compartment Mass (includes tube, contents, sealing assembly and rotor cap)	82.5 g
Rotor Mass (Weight)	11.4 kg (24.5 lb)
K Factor at Maximum Speed	25.2

*Speed in revolutions per minute (rpm) is related to angular velocity, w , according to the following:

$$\omega = (\text{rpm}) \left(\frac{2\pi}{60} \right) = (\text{rpm})(0.10472)$$

Where $w = \text{rad/s}$. All further references in this manual to speed will be designated as rpm.

**With tubes filled with a homogeneous solution having an average density of 1.7 g/ml or less.

Accessories

a. Accessories Supplied

The accessories that are supplied with the TV-860 Rotor, Catalog Number 53290 (rotor with complete accessories), are listed in Table 1-2. Those items indicated by an asterisk (*) are not supplied with the TV-860 Rotor, Catalog Number 53289, which is the rotor with basic accessories only.

Note In order to seal the ULTRACRIMP® Tubes supplied with the rotor (Catalog No. 53290), you must have an ULTRACRIMP® Sealing Tool (Catalog No. 03920) and a Tool Kit (Catalog No. 12771)

To order or for information, in the United States, call Thermo Fisher Scientific toll-free 1-800-522-7746; outside the United States, contact the nearest Thermo Fisher Scientific office (see back cover) or your local representative for Thermo Fisher Scientific products. Be sure to provide a description of the part, the catalog number, plus the rotor model and serial number. Thermo Fisher Scientific product information is available on our internet web site at <http://www.thermo.com/centrifuge>.

Table 1-2. Accessories Supplied

Quantity	Catalog Number	Description
8	03989*	ULTRACRIMP® Tubes, Polyallomer, 35 ml (25/pkg)**
8	03994*	Rotor Cap
1	03935*	Tube Rack, 4-place (1 pair)
1	12826*	ULTRACRIMP® Tube Removal Tool
1	51346	Overspeed Decal, 60 000 rpm (extra)
1	51942	Rotor Stand
1	61556	Lubricant
1	52384	Ultraspeed Centrifuge Rotor Log Book
1	53292	Instruction Manual

*Not supplied with Cat. No. 54289 (rotor with basic accessories only).

**Each package of tubes is supplied with an equal number of tube caps and tube plugs.

b. Tool Kit, Catalog No. 12771

This kit is required for use with the TV-860 Rotor when used with the ULTRACRIMP® Tube Sealing System. Table 1-3 lists the items supplied in the Tool Kit. If you already have these tools, it is not necessary to order the kit. Any of the tools listed can also be ordered individually; to do so, specify the catalog number given in the table.

Table 1-3. Tool Kit (Cat. No. 12771) for use with ULTRACRIMP® Tube Sealing System

Catalog Number	Description
52580	Torque Wrench, 3/8-inch drive
54519	Rotor Body Cap Tool
52540	Rotor Holding Fixture (with mounting screws)

c. Optional Sealing Assembly

If you prefer, you can use the TV-860 with the optional sealing assembly which consists of a rotor body cap and a tube plug. When the body cap is tightened, the tube plug is pushed into the tube to form the seal. To use this style sealing assembly, you must purchase the TV-860 Rotor with basic accessories only, Catalog No. 52389; a Tool Kit, Catalog No. 52599; the tubes; and the sealing assembly components listed in Table 1-4. Table 1-5 lists the items supplied in the Tool Kit; any of these tools can also be purchased individually.

Table 1-4. Optional Sealing Assembly Components

Catalog Number	Description
03141	Tube, Polyallomer, 36 ml (package of 25)
52556	Rotor Body Cap, Red (each); 8 required
52557	Tube Plug (package of 8)
52555	Tube Plug O-ring (package of 16)
52596	Titanium Seal Ring, Blue (package of 8)
09073	Tube Trimmer (optional)

Table 1-5. Tool Kit (Cat. No. 52599) for use with Optional Sealing Assembly

Quantity	Catalog Number	Description
1	52448	Tube Plug Removal Tool
1	52577	Torque Wrench Adapter
1	52580	Torque Wrench
1	52540	Rotor Holding Fixture (with mounting screws)

RUN PREPARATION

Contents

- “Prerun Safety Checks” on page 2-2
- “Rotor Holding Fixture” on page 2-2
- “Rotor Precool” on page 2-3
- “Chemical Compatibility” on page 2-3
- “Sample Preparation” on page 2-3

Prerun Safety Checks

To ensure safe performance of the rotor, the following items should be checked before every run:

- a. read the Safety Information Page in front of this manual.
- b. make sure each tube compartment and rotor cap is clean and that there is no sign of corrosion, and that the centrifuge chamber and the drive spindle are clean.
- c. make sure the rotor itself is clean and shows no sign of corrosion or cracking.
- d. check the centrifuge chamber and drive spindle to be sure they are clean and free of scratches and burrs.
- e. verify that the proper overspeed decal is firmly attached to the bottom of the rotor; the decal should have 15 black segments (refer to paragraph 6-5 for Overspeed Decal Replacement procedure).
- f. check the chemical compatibility of all materials used (see Appendix A).
- g. be sure the rotor caps are torqued to the recommended value (paragraph 4-1).
- h. be sure the proper environment has been selected for operation; for example, controlled ventilation or isolation, if required.
- i. check the top speed capability of the tube being used; observe the CAUTION below.



WARNING If using a Thermo Fisher Scientific tube (or bottle) assembly other than the one specified, refer to the Rotors, Tubes, Bottles and Adapters Catalog for the maximum speed. When using a tube or bottle assembly other than those supplied by Thermo Fisher Scientific, be sure to check its top speed capability. When in doubt, do a test run for the desired application. Exceeding the top speed capability of the tube or bottle assembly can result in tube breakage.

Rotor Holding Fixture

The rotor holding fixture (Cat. No. 52540) is used to hold the rotor during the rotor loading and unloading procedure to prevent disturbance of the gradient.

Fasten the fixture to a workbench with the two mounting screws supplied as shown in figure 2-1.

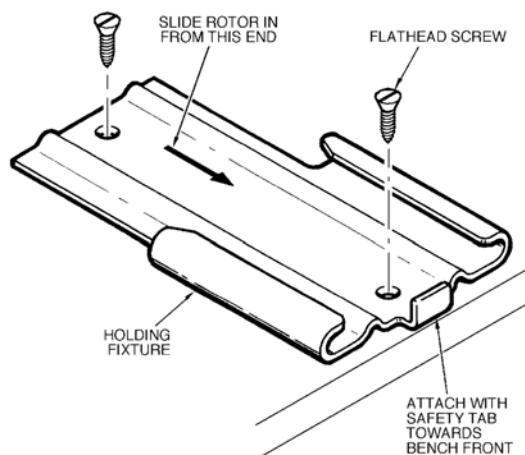


Figure 2-1. Rotor Holding Fixture Installation

Rotor Precool

If samples are routinely processed around 4°C or below, the rotor should be stored in a refrigerator or a cold room. If this is not possible, the rotor can easily be precooled in a Thermo Fisher Scientific Ultraspeed Centrifuge. Refer to the Ultraspeed Centrifuge Instruction Manual for precooling directions. Be careful not to precool the rotor at its critical speed (read paragraph 3-3).

Chemical Compatibility

The critical components of the TV-860 Rotor that are apt to come in contact with solution are: the rotor body (titanium), the tubes (polyallomer), and all sealing assembly components. ULTRACRIMP® sealing assembly components include: rotor caps and tube caps (aluminum) and the tube plugs (Buna N). Optional sealing assembly components include: tube plugs (Delrin®) and rotor body caps (aluminum). The chemical compatibility of rotor elements and accessory materials is given in the Appendix A.

Because no organized chemical resistance data exists for materials under the stress of centrifugation, this data is intended to be used only as a guide to the selection of tube materials. When in doubt, we recommend pretesting of sample lots.

Sample Preparation

a. ULTRACRIMP® Tubes, Catalog No. 03989

Prepare the ULTRACRIMP® tubes for use by following the tube filling procedure in the ULTRACRIMP® Tube Sealing System Instruction Manual.

b. Optional Tubes, Catalog No. 03141

When using optional tubes, Catalog Number 03141, each tube must contain a total of 34 ml of solution; the recommended capacity for each tube is: 32.0 ml of gradient, 1.0 ml of sample and 1.0 ml of overlay (e.g. mineral oil).



CAUTION To prevent the tube from collapsing or leaking during centrifugation, each tube must always contain 34 ml of solution. Also, use a new tube for each operation — do not reuse these tubes.

SPECIAL OPERATING CONSIDERATIONS

Contents

- “Compartment Loads in Excess of Design Mass” on page 3-2
- “Precautions to Prevent Precipitation of Cesium Chloride” on page 3-2
- “Critical Speed” on page 3-7

Compartment Loads in Excess of Design Mass

A recommended design mass has been established for each ultraspeed centrifuge rotor, representing the maximum mass that each tube compartment can contain at top speed operation. To prevent rotor failure, the total contents of each compartment, including the specimen, rotor cap, tube cap, tube plug, and the tube, should not exceed the recommended figure unless the rotor speed is reduced proportionately. (If using the optional sealing assembly, include the weight of all components.)

Strict adherence to the maximum allowable compartment mass or reduced speed is required to prevent rotor failure. Observe the WARNING on the Safety Information Page in front of this manual.

The design mass for each tube compartment of the TV-860 Rotor is 82.5 g at 60 000 rpm. This figure is based on the use of an ULTRACRIMP[®] tube filled with a liquid at 1.7 specific gravity. If the density is greater than 1.7 g/ml, use the following formula to determine the reduced speed:

$$\text{Reduced Speed} = 60000 \sqrt{\frac{1.7}{\text{Average Fluid Density (g/ml)}}}$$

Refer to figures 3-1, 3-2, and 3-3 for derating factors when using cesium chloride solutions.

Precautions to Prevent Precipitation of Cesium Chloride

a. Reducing Speed to Prevent Precipitation

Observe the WARNING on reducing rotor speed found on the Safety Information Page in front of this manual.

Maximum speed must be reduced for an average fluid density greater than 1.7 g/ml (square-root reduction) to prevent excessive hydraulic pressure in a rotor. Although the standard formula (see paragraph 3-1) pertains to sucrose and similar gradient materials, it will not prevent precipitation of heavy crystals when material such as cesium chloride (CsCl) is used in an ultraspeed centrifuge.

When solid, crystalline CsCl forms, it places a density of 4 g/ml along the outside edge of each tube compartment. This density is dangerously high and can cause the rotor to fail, with subsequent sample loss. Therefore, cesium chloride solutions must be run at a reduced speed to avoid this precipitation. The allowable speed is determined by the average density of the CsCl solution and the run temperature. Figure 3-1 should be used to determine the maximum operating speed. Curves are given for specific average densities at 5°C and 25°C that will prevent precipitation. Also, figure 3-1 includes a curve that shows the standard speed (square-root) reduction to avoid excessive hydraulic pressure only. The standard, square-root reduction formula cannot be used when precipitation of CsCl must be considered. For example, standard speed reduction would allow you to run a solution with a density of 1.6g/ml at 60 000 rpm; however, you can see that a cesium chloride solution will precipitate at this speed. The graph shows that the highest speed you can run a cesium chloride solution having a density of 1.6 g/ml is 55 000 rpm at 25°C or 45 000 rpm at 5°C.

Note The solubility limit of cesium chloride in an aqueous solution is 1.86 g/ml at 25°C and 1.81 g/ml at 5°C.

b. The Gradient Shape

The curves in figures 3-2 and 3-3 show the gradient shape at equilibrium, 5°C and 25°C respectively, for tubes filled with a CsCl solution. Figure 3-2 illustrates the shape of the gradient curve at various speeds. For example, to find the gradient profile of a tube filled with a 1.6 g/ml solution at 45000 rpm at 5 °C, locate the 45 000 rpm curve. The gradient profile is 1.43 g/ml at the meniscus to 1.81 g/ml at the bottom of the tube. Use figure 3-3 to find the shape of the gradient (1.6 g/ml) that corresponds to 55 000 rpm at 25°C. Interpolate to find gradient shapes at speeds other than those illustrated.

Next, use figures 3-2 and 3-3 to determine the position of a band of particles in the tube when run at 5°C and 25°C respectively. Using the example given above (i.e., 45 000 rpm at 5°C), use figure 3-2 to locate the position of a band of particles having a density of 1.52g/ml. Find 1.52 on the side scale (DENSITY , g/ml) and follow the line across until it intersects the curve at 45000 rpm. At this point, follow the line down the graph until it intersects the bottom scale (RADIUS, cm) in this example the line corresponds to a radius of 6.53cm. To find this band of particles after reorientation, continue to follow the line to the VOLUME (ml) of Fluid Above Band scale; you can see that the band of particles after reorientation is at 7.5. This means that there will be 7.5 ml of gradient above the band and 26.5ml below the band; therefore, measuring from the top of the meniscus, the particles will be located at 78% of the fluid column.

Use figure 3-3 to locate a band of particles having the same density (i.e., 1.52 g/ml) run at 25°C. In this example, the band of particles can be located at a radius of 6.55 cm. Continue to follow the line to the VOLUME (ml) of Fluid Above the Band scale to locate the band of particles after reorientation which is at 9.2. This means there is 9.2ml above the band and 24.8 ml below the band. Again, measuring from the top of the meniscus, the particles will be located at 73% of the fluid column.

3 SPECIAL OPERATING CONSIDERATIONS

Precautions to Prevent Precipitation of Cesium Chloride

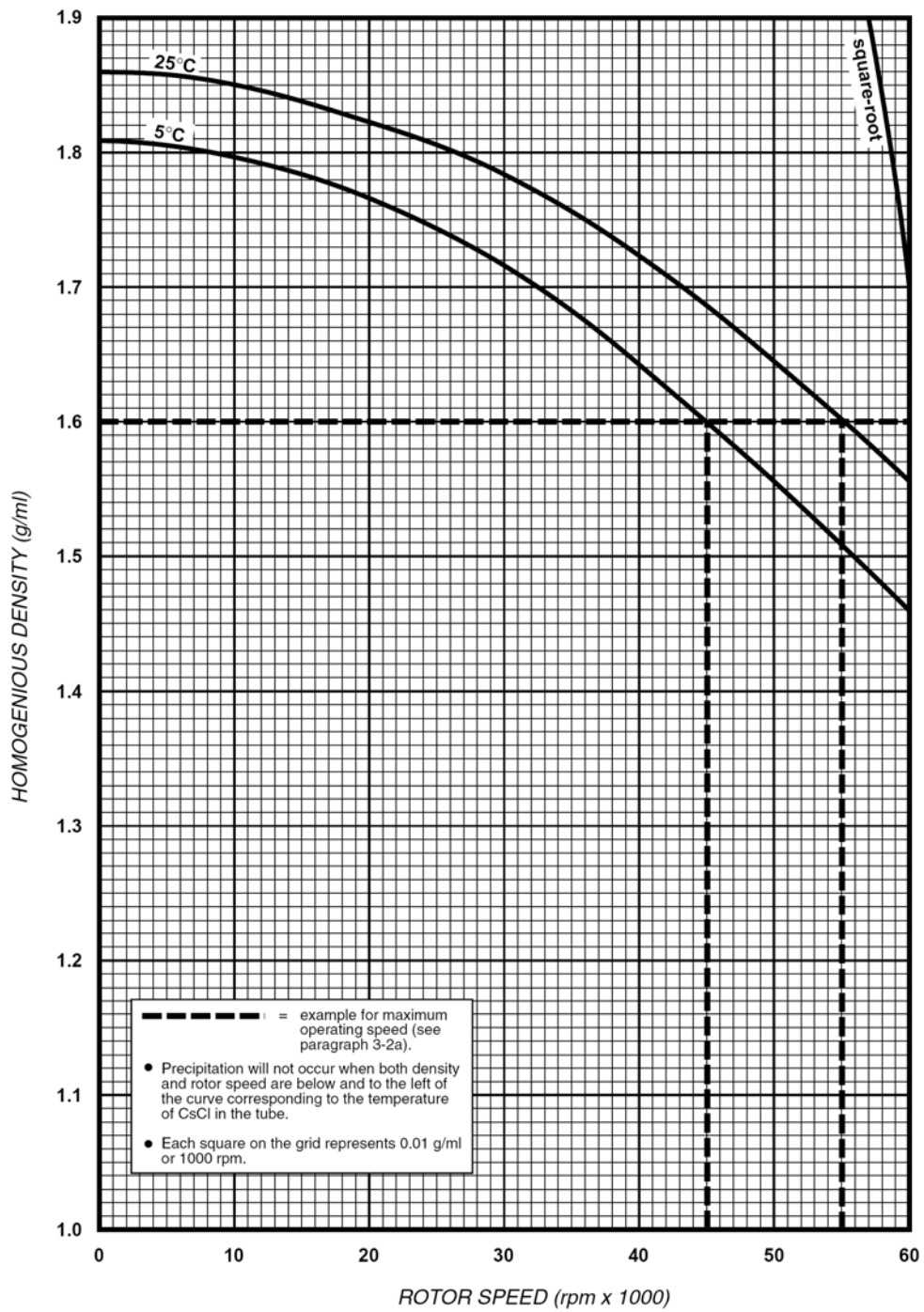


Figure 3-1. CsCl Precipitation Curves for the TV-860 Rotor

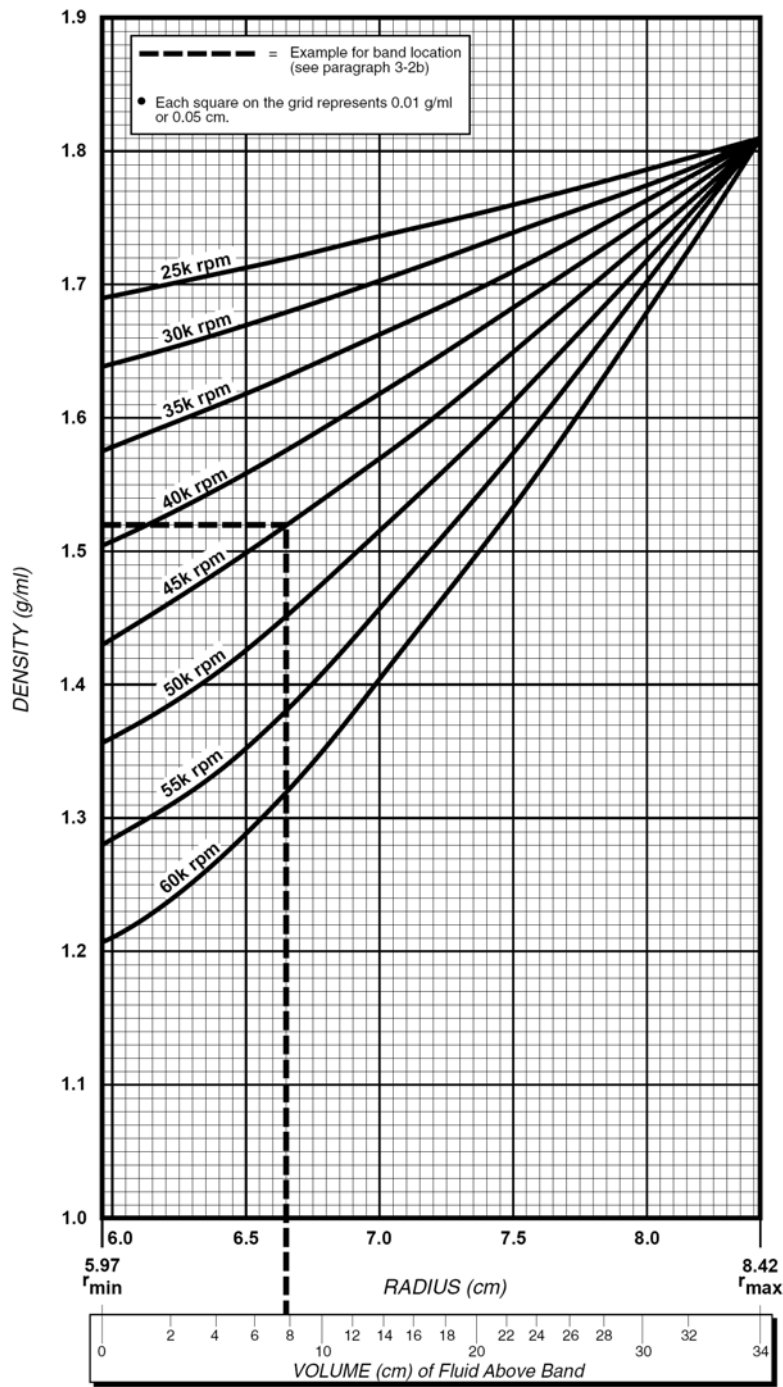


Figure 3-2. CsCl Density Gradient Shapes at Equilibrium (5°C)

3 SPECIAL OPERATING CONSIDERATIONS

Precautions to Prevent Precipitation of Cesium Chloride

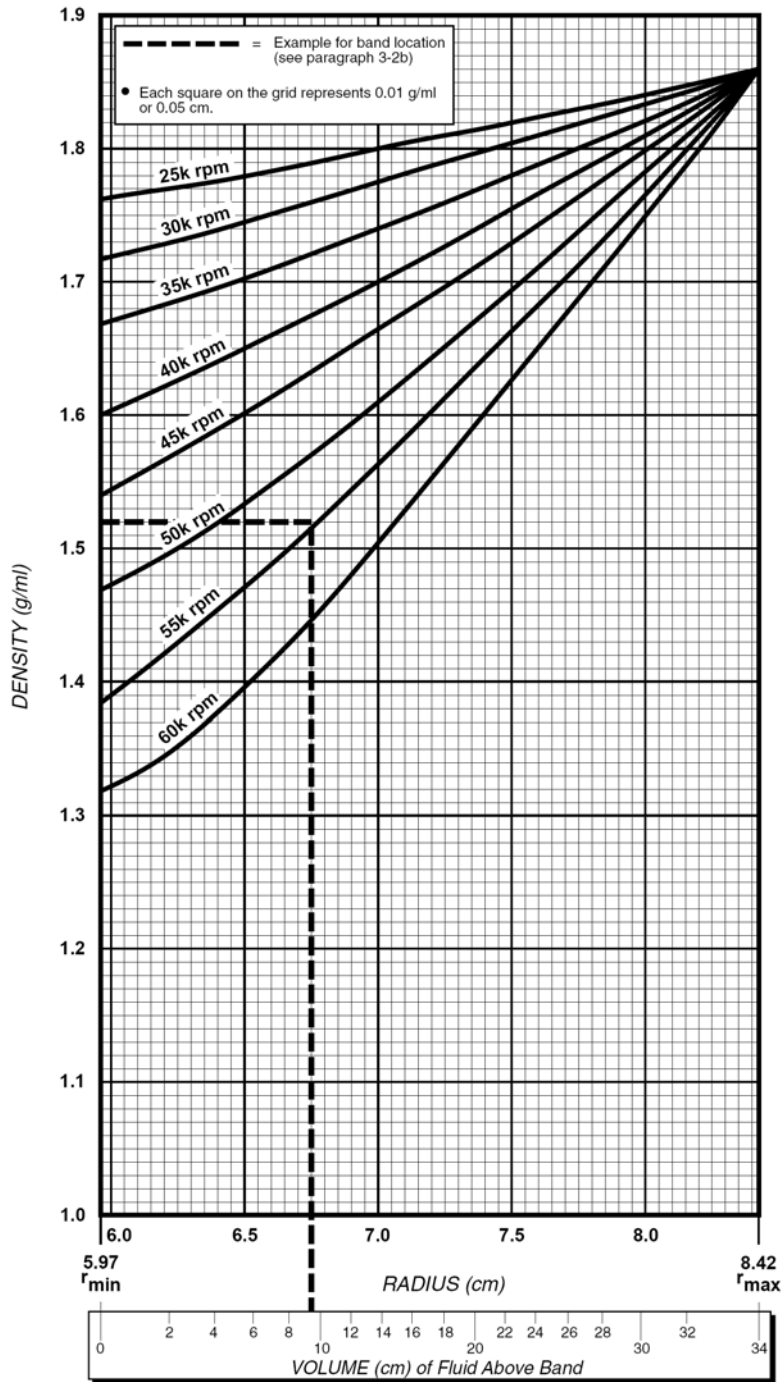


Figure 3-3. CsCl Density Gradient Shapes at Equilibrium (25°C)

Critical Speed

The critical speed is that speed at which any rotor imbalance will produce a driving frequency equal to the resonant frequency of the rotating system (for example, the rotor and the centrifuge drive). At this speed, the rotor may produce large amplitude vibrations which can be felt in the centrifuge frame. Mass imbalance will contribute to increased vibration intensity at the critical speed. Avoid operating the rotor at the critical speed, which is 700 rpm for the TV-860 Rotor. Operation at the critical speed will have a detrimental effect on centrifuge component life. Observe the CAUTION on the Safety Information Page in the front of this manual.

OPERATION

Contents

- “Rotor Loading and Sealing” on page 4-2
- “Rotor Balancing” on page 4-5
- “Rotor Installation” on page 4-5
- “Acceleration and Deceleration” on page 4-6
- “Tube Removal” on page 4-6
- “Tube Trimmer Accessory (for use with Optional Tubes)” on page 4-7
- “Rotor Log Book” on page 4-7



WARNING The stresses that this rotor withstands during centrifugation eventually cause the rotor body to weaken, increasing the possibility of rotor failure. For your safety, we recommend that this rotor be withdrawn from service after it has been used for 5 000 runs or 10 000 hours, whichever comes first. Failure to do so can cause rotor failure with subsequent sample loss and damage to the rotor and/or centrifuge. If the material being processed is hazardous, take all necessary precautions to prevent exposure should rotor failure occur.

Rotor Loading and Sealing

To be sure that each tube and rotor compartment is properly sealed for operation, these instructions must be carefully followed. If necessary, precool the rotor, tube plugs and rotor body caps to the same temperature. (See the Ultraspeed Centrifuge Instruction Manual for precooling instructions.)

Prepare samples according to the instructions in Chapter 2, paragraph 2-5 then load and seal the rotor as follows:

a. Using ULTRACRIMP® Tubes and Sealing System

1. Place the rotor in the rotor holding fixture (Cat. No. 52540). Gently put filled tubes into the tube compartments, balancing the rotor as explained in paragraph 4-2.
2. Apply a light coat of lubricant (Cat. No. 61556) to each rotor cap (Cat. No. 03994).
3. Place a rotor cap in each compartment that contains a filled tube.



CAUTION Do not put a rotor cap in an empty tube compartment. Rotor caps should only be installed in compartments that contain a filled tube.

4. Hand tighten each rotor cap.
5. Using the torque wrench (Cat. No. 52580) with the rotor body cap tool (Cat. No. 54519), tighten each rotor cap to 180 in lbs (20.3 Nm).



CAUTION Do not undertighten or overtighten the rotor caps. Always use a torque value not less than 180 in lbs (20.3 Nm) or greater than 200 in lbs (22.6 Nm).

b. Using Optional Tubes and Sealing Assembly (refer to figure 4-1)

Precool the rotor, seal rings, tube plugs and body caps to the same temperature if necessary.

Prepare sample as explained in paragraph 2-5 then load and seal the rotor as follows:



CAUTION Before every run inspect the counterbore area where the seal ring fits into the tube compartment be sure there are no nicks or scratches in the area and that the area is clean.

1. Place the rotor in the rotor holding fixture with either compartment #1 or #5 facing you.

2. Install a seal ring (Cat. No. 52596), chamfer side up, into the counterbore of each rotor compartment that will contain a tube (see figure 4-1).



CAUTION The chamfer side of each seal ring must be facing upward, if not the tube may stick in the rotor compartment.

3. Gently put the filled tubes into the rotor tube compartments (balancing the rotor as described in paragraph 4-2).
4. Lubricate each tube plug O-ring (Cat. No. 52555) with a light film of vacuum grease (Cat. No. 65937).
5. Insert a lubricated O-ring in the groove of each tube plug (Cat. No. 52557). Place a tube plug (with O-ring) into the top of each tube in the rotor.
6. Apply a light coat of lubricant (Cat. No. 61556) to the threads of the rotor body caps, then place a body cap into each tube compartment that contains a filled tube. Finger tighten each body cap.



CAUTION Do not put a body cap in an empty tube compartment. Body caps should only be installed in compartments that contain a filled tube.

4 OPERATION

Rotor Loading and Sealing

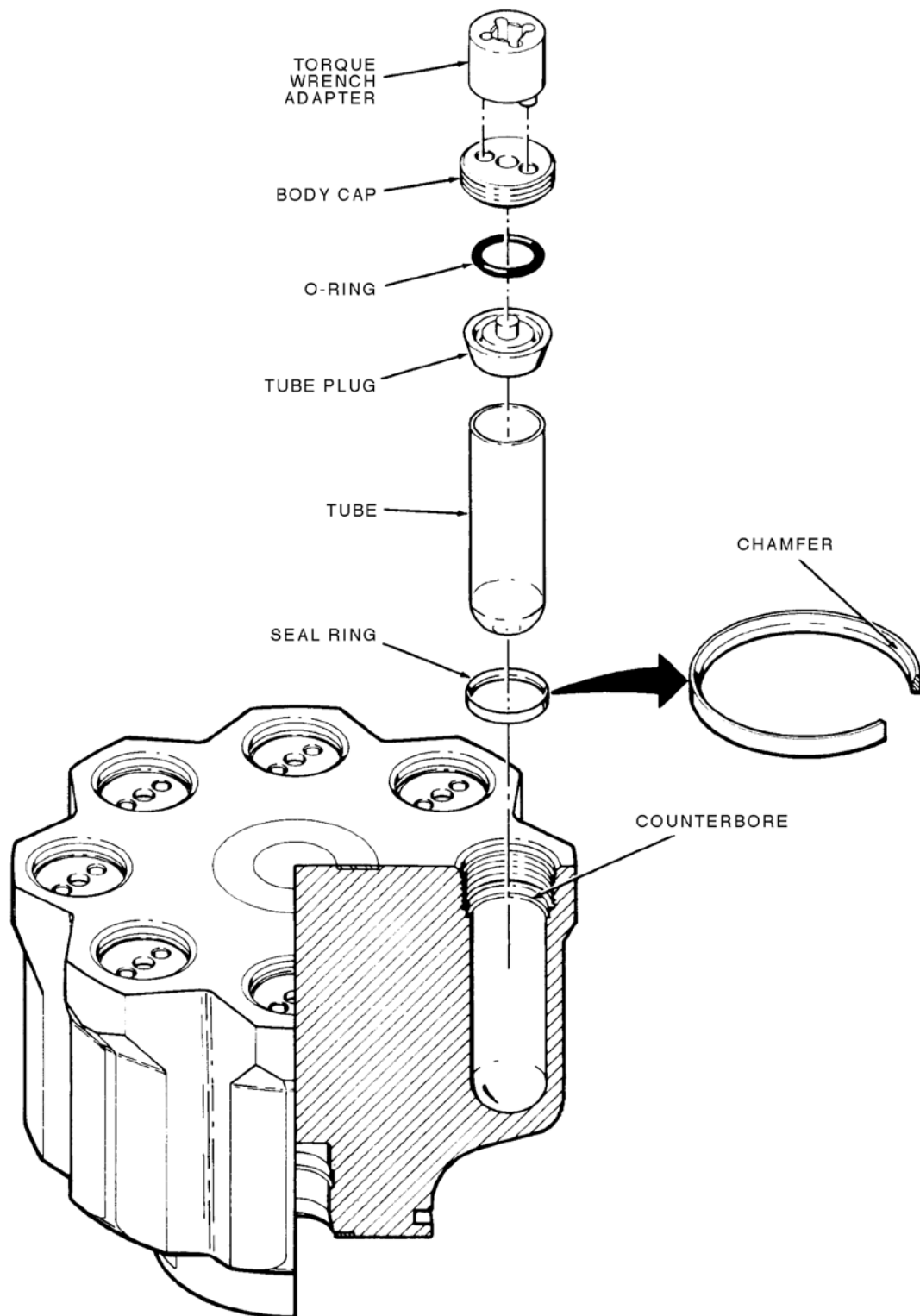


Figure 4-1. Optional Rotor Sealing Assembly

7. Preset torque wrench (Cat. No. 52580) to 220 in lb. (24.9 Nm) according to the instructions supplied. Attach the torque wrench to the torque wrench adapter (Cat. No. 52577) and tighten each body cap to 220 in lb. When the torque being applied equals the preset torque value, the wrench will automatically release (approximately 10° arc). This release, or break away, can be distinctly felt and heard by the operator and signals that the operation is complete. When pressure is released, the wrench resets itself and is ready for use.



CAUTION Do not undertighten or overtighten the body caps. Always use a torque value between 180 in lb. (20.3 Nm) and 240 in lb. (27.1 Nm).

Rotor Balancing

Always balance the rotor according to the following criteria:

- a. observe CAUTION on the Safety Information Page in the front of this manual.
- b. pairs of tubes containing a sample of identical specific gravity must be balanced to within 0.5 gram and placed in opposing compartments.
- c. the rotor can be operated at any speed up to its maximum speed with groups of two, four, six or eight samples, as long as each sample is paired to another of identical specific gravity and the tubes are placed in opposite rotor compartments as shown in figure 4-2.



CAUTION When the rotor is operated with two, four or six samples placed in opposing tube compartments, the remaining tube compartments must be empty. Do not put rotor body caps in empty compartments.

- d. if an uneven number of samples is to be run, the odd sample must be counterbalanced with a tube that contains a solution of the same specific gravity.

Rotor Installation

Note Whenever you carry the rotor after it has been loaded, hold it vertically with both hands underneath it; the rotor must always be moved slowly and carefully so the gradient will not be disturbed.

To install the rotor in the centrifuge, remove it from the rotor holding fixture and carefully lower it onto the drive adapter in the centrifuge rotor chamber. The rotor will snap in place on the drive adapter. To make sure the rotor is fully seated, pull up on it gently if there is a small amount of resistance, it is properly seated.

The rotor is ready for operation per form the centrifuge run according to the instructions for the ultracentrifuge.

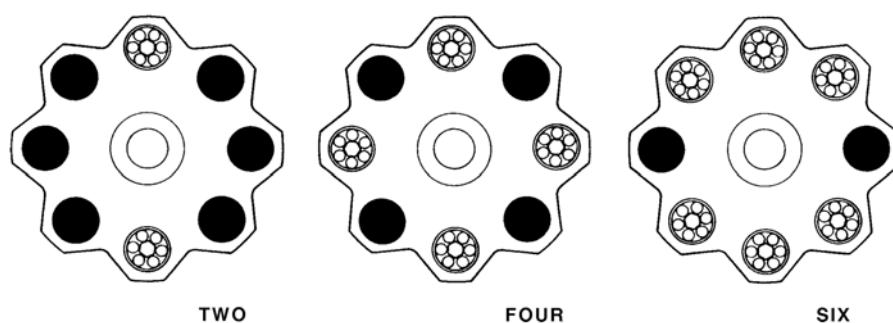


Figure 4-2. Rotor Balancing

Acceleration and Deceleration

When using the TV-860 Rotor for density gradient separations, the automatic rate controller or the REOGRAD mode must be used for slow acceleration and slow deceleration. For instructions on the use of these controls, refer to the ultracentrifuge instruction manual.

Tube Removal

To remove the tubes from the rotor:

a. ULTRACRIMP® Tubes

1. Place the rotor in the rotor holding fixture.
2. Remove the rotor caps from the rotor using the torque wrench with the body cap tool.
3. Remove tubes from the rotor. Because the tubes expand slightly during centrifugation, use the ULTRACRIMP® Tube Removal Tool (Cat. No. 12826), needle-nose pliers, or a hemostat to grasp the tube and gently pull it from the rotor.

Note You may notice that the inner side wall of a tube may appear slightly flattened after centrifugation. This is a normal effect of centrifugation and should not be cause for any concern.

b. Optional Tubes, Catalog No. 03141

1. Place the rotor in the rotor holding fixture.
2. Remove the rotor body caps using the torque wrench and torque wrench adapter.
3. Push the tube plug tool onto the stem of the tube plug and gently rock the plug until it pops loose, then remove the plug.



WARNING Be extremely careful when removing the tube plugs, particularly if the samples are radioactive or pathogenic. If liquid is trapped between the plug and the tube wall, it may be forced out by an air bubble when the plug is loosened. This is one reason we recommend the use of an overlay, such as mineral oil, on the sample.

4. Grasp the edge of each tube with a hemostat or forceps and slowly pull the tube out of the rotor. If you have difficulty getting hold of the edge of the tube, insert a small tool such as an 18 gauge needle or a spatula between the wall of the rotor tube compartment and the tube and pry the tube until it can be gripped with the hemostat or forceps.
5. Remove seal rings.

Clean all sealing assembly components according to the procedure given in paragraph 6-2.

Tube Trimmer Accessory (for use with Optional Tubes)

The rim of the optional tubes (Cat. No. 03141) will be deformed by the sealing assembly during operation. This may cause the tubes to leak when used with certain density gradient fractionators. To correct this problem, a tube trimmer is supplied with the rotor. After the tube has been removed from the rotor, insert it into the trimmer. The deformed portion of the tube will extend above the top of the trimmer. Use a razor blade and very carefully cut it away from the rest of the tube.

Note The tube trimmer is not required when using the TV-860 Rotor with ULTRACRIMP[®] tubes (Cat. No. 03989).

Rotor Log Book

An Ultraspeed Centrifuge Rotor Log Book is supplied with the TV-860 so you can easily record all data necessary to meet the warranty stipulation that if a defective rotor is returned to Thermo Fisher Scientific it is accompanied by an up-to-date history of the rotor.

Each time the TV-860 Rotor is used, record the run in the log book as shown in Figure 4-3, Sample Rotor Log Sheet..

Thermo Scientific Centrifuge and Rotor Log Book					RUN TIME (List by Rotor Used)								This log is for use with one centrifuge ONLY:	
Date	Operator	Rev. Count @ Run Start	TEMP	SPEED	Rotor AH-629 S/N 8731384		Rotor TV-865 S/N 9130129		Rotor T-1270 S/N 8931255		Rotor T-880 S/N 9030040		Model: ULTRA 80	Ser. No.: 9102448
					HRS	MIN	HRS	MIN	HRS	MIN	HRS	MIN	Remarks*	
09/04/91	J. JONES	00410290	4	57.0			05	30						PLASMID PREP.
09/05/91	B. SMITH	00429100	4	21.0	26	00								SUCROSE GRADIENT
09/07/91	J. JONES	00461860	21	70.0					18	00				LIPOPROTEIN SEP.

Figure 4-3. Sample Centrifuge/Rotor Log Sheet

OPERATION

This chapter contains the information necessary to prepare the TFT-45.6 Rotor for operation and includes important safety information.

Contents

- “Relative Centrifugal Force (RCF) Determination” on page 5-2
- “Calculation of Sedimentation Time in Aqueous (Non-gradient) Solutions” on page 5-5
- “Calculation of Sedimentation Time in Gradient Solutions” on page 5-6

Relative Centrifugal Force (RCF) Determination

Relative centrifugal force (RCF) refers to the force during centrifugation that moves the particulate outward from the center of rotation. This force is proportional to the radial distance and the square of the rotor speed. The RCF value, also known as g force, is determined by the following formula:

$$\text{RCF} = 11.17(r) \left(\frac{\text{rpm}}{1000} \right)^2$$

when r = the radius in centimeters from the centerline of the rotor to the point in the tube where RCF value is required

and rpm = the rotor speed in revolutions per minute

Figure 5-1 shows the minimum, average, and maximum radii of the TV-860. Table 5-1 gives the RCF value at each radius at speeds from 20 000 rpm to 60 000 rpm (in increments of 500 rpm). The RCF value at any other speed can be calculated by using the given formula.

Note The radii values given are the actual rotor specifications. These values do not take the thickness of the tube into consideration.

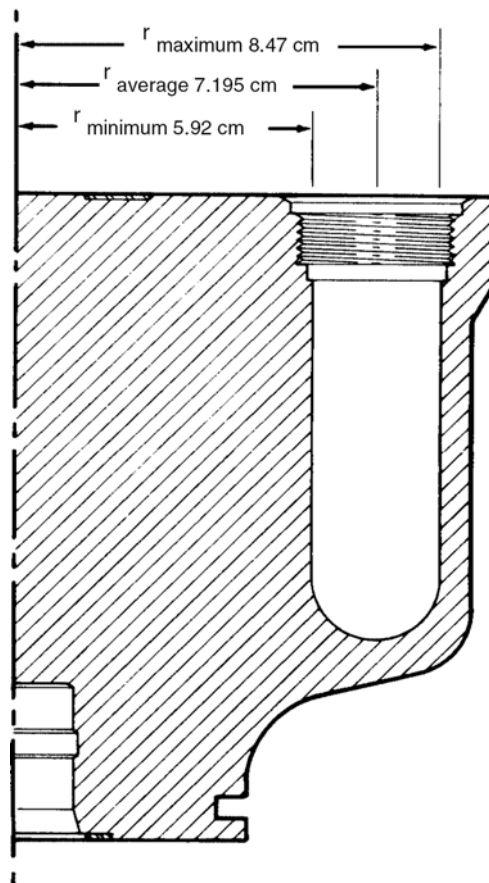


Figure 5-4. TV-860 Rotor Radii

Table 5-1. TV-860 Rotor: RCF Values and K Factors *

Speed (rpm)	RCF			K Factors
	r _{max} 8.47 cm	r _{avg} 7.195cm	r _{min} 5.92cm	
20 000	37 844	32 147	26 451	227
20 500	39 760	33 775	27 790	216
21 000	41 723	35 442	29 162	205
21 500	43 733	37 150	30 567	196
22 000	45 791	38 898	32 005	187
22 500	47 896	40 686	33 476	179
23 000	50 049	42 515	34 981	171
23 500	52 248	44 383	36 518	164
24 000	54 495	46 292	38 089	157
24 500	56 790	48 241	39 692	151
25 000	59 131	50 230	41 329	145
25 500	61 520	52 259	42 999	139
26 000	63 956	54 329	44 701	134
26 500	66 440	56 439	46 437	129
27 000	68 971	58 588	48 206	124
27 500	71 549	60 778	50 008	120
28 000	74 174	63 009	51 843	116
28 500	76 847	65 279	53 711	112
29 000	79 567	67 590	55 612	108
29 500	82 334	69 940	57 546	104
30 000	85 149	72 331	59 514	101
30 500	88 011	74 762	61 514	97.4
31 000	90 920	77 234	63 547	94.3
31 500	93 877	79 745	65 614	91.3
32 000	96 881	82 297	67 713	88.5
32 500	99 932	84 889	69 846	85.8
33 000	103 030	87 521	72 012	83.2
33 500	106 176	90 193	74 210	80.8
34 000	109 369	92 906	76 442	78.4
34 500	112 609	95 658	78 707	76.1
35 000	115 897	98 451	81 005	74.0
35 500	119 232	101 284	83 336	71.9
36 000	122 614	104 157	85 700	69.9

5 OPERATION

Relative Centrifugal Force (RCF) Determination

Table 5-1.TV-860 Rotor: RCF Values and K Factors*

Speed (rpm)	RCF			K Factors
	r _{max.} 8.47 cm	r _{avg.} 7.195cm	r _{min.} 5.92cm	
36 500	126 044	107 070	88 097	68.0
37 000	129 521	110 024	90 527	66.2
37 500	133 045	113 018	92 990	64.4
38 000	136 617	115 052	95 487	62.8
38 500	140 236	119 126	98 016	61.1
39 000	143 902	122 240	100 578	59.6
39 500	147 615	125 394	103 174	58.1
40 000	151 376	128 589	105 802	56.6
40 500	155 184	131 824	108 464	55.2
41 000	159 039	135 099	111 158	53.9
41 500	162 942	138 414	113 886	52.6
42 000	166 892	141 769	116 647	51.4
42 500	170 889	145 165	119 441	50.2
43 000	174 934	148 601	122 268	49.0
43 500	179 026	152 077	125 128	47.9
44 000	183 185	155 593	128 021	46.8
44 500	187 351	159 149	130 947	45.8
45 000	191 585	162 746	133 906	44.8
45 500	195 866	166 382	136 898	43.8
46 000	200 195	170 059	139 923	42.8
46 500	204 570	173 776	142 982	41.9
47 000	208 993	177 533	146 073	41.0
47 500	213 464	181 331	149 198	40.2
48 000	217 981	185 168	152 355	39.3
48 500	222 546	189 046	155 546	38.5
49 000	227 158	192 964	158 769	37.7
49 500	231 818	196 922	162 026	37.0
50 000	236 525	200 920	165 316	36.2
50 500	241 279	204 959	168 639	35.5
51 000	246 080	209 038	171 995	34.8
51 500	250 929	213 156	175 384	34.2
52 000	255 825	217 315	178 806	33.5
52 500	260 769	221 515	182 261	32.9
53 000	265 759	225 754	185 749	32.3

Table 5-1. TV-860 Rotor: RCF Values and K Factors*

Speed (rpm)	RCF			K Factors
	r _{max} 8.47 cm	r _{avg} 7.195cm	r _{min} 5.92cm	
53 500	270 979	230 034	189 270	31.7
54 000	275 882	234 354	192 825	31.1
54 500	281 015	238 713	196 412	30.5
55 000	286 195	243 114	200 032	30.0
55 500	291 422	247 554	203 686	29.4
56 000	296 697	252 035	207 372	28.9
56 500	302 018	256 555	211 092	28.4
57 000	307 388	261 116	214 845	27.9
57 500	312 804	265 717	218 630	27.4
58 000	318 268	270 358	222 449	26.9
58 500	323 779	275 040	226 301	26.5
59 000	329 337	279 762	230 186	26.0
59 500	334 943	284 523	234 104	25.6
60 000	340 596	289 325	238 055	25.2

* These values do not take the thickness of the tube into consideration.

Calculation of Sedimentation Time in Aqueous (Non-gradient) Solutions

The time required to sediment a particle in water at 20°C through the maximum rotor path length (that is, the distance between r_{minimum} and r_{maximum}) can be calculated using the equation:

$$t = \frac{K}{S_{20,w}}$$

where:

t = sedimentation time in hours

K = the clearing factor for the rotor (defined on the next page)

S_{20, w} = the sedimentation coefficient for the particle of interest in water at 20°C as expressed in Svedbergs¹

¹ The sedimentation coefficient (S) in seconds, for a particle in a centrifugal field is defined by the equation $S = (dx/dt) [1/(\omega^2 x)]$; where dx/dt = sedimentation velocity of the particle in cm/s; ω = rotor speed in rad/s; and x = the distance of the particle from the axis of rotation in centimeters. Conventionally, experimentally determined values of sedimentation coefficients are multiplied by 10¹³ to convert them to Svedberg units (S), so a particle with an experimentally determined sedimentation coefficient of 10⁻¹¹ seconds is usually referred to in the literature as a "100 S particle." Since the value determined for the sedimentation coefficient is dependent on the density and viscosity of the solution in which centrifugation is performed, values are usually reported for the standard conditions of infinite dilution in water at 20°C, and designated S_{20, w}.

5 OPERATION

Calculation of Sedimentation Time in Gradient Solutions

The clearing (or K) factor is defined by the equation:

$$K = (253000) \left[\ln \left(\frac{r_{\text{maximum}}}{r_{\text{minimum}}} \right) \right] \div \left(\frac{\text{rotor speed}}{1000} \right)^2$$

Where r_{maximum} and r_{minimum} are the maximum and minimum rotor radii, respectively, and rotor speed is expressed in rpm.

K factors for the TV-860 Rotor at speeds from 20 000 rpm to 60 000 rpm (in increments of 500 rpm) are listed in Table 5-1.

For example: the TV-860 Rotor has a K factor of 25.2 at the maximum permitted speed (60 000 rpm). If the particles to be sedimented have a sedimentation coefficient of 10 S, the estimated run time required at maximum speed for the outer row will be:

$$t = \frac{25.2}{10S} \text{ 2.52 hours} = 2 \text{ hours, 31 minutes}$$

Note that the calculation assumes particles in water at 20°C. If the suspending medium is denser or more viscous than water, the sedimentation time will be greater.

Calculation of Sedimentation Time in Gradient Solutions

The time required to sediment a particle through a density gradient can be calculated using the following formula:

$$t = \frac{K^1}{S_{20,w}}$$

where:

t = sedimentation time in hours

K^1 = the clearing factor for the rotor (the value of K^1 is dependent on the gradient being used, the temperature of the gradient, and the density of the particle being sedimented).

$S_{20,w}$ = the sedimentation coefficient for the particle of interest in water at 20°C as expressed in Svedbergs¹

Table 5-2 gives K^1 factors for the TV-860 Rotor when operated at maximum speed (i.e., 60000 rpm) with particles ranging in density from 1.1 g/cm³ to 1.5 g/cm³. In this case, the K^1 factors are based on the use of a 5% -20% (w/w) linear sucrose density gradient at 4°C.

Table 5-2. K^1 Factors for TV-860 (at maximum speed)

Particle Density (g/cm ³)	K^1 Factor (at 60000 rpm)
1.1	151
1.2	80
1.3	70
1.4	66
1.5	64

DESCRIPTION

Observe all WARNINGS and CAUTIONS found on the Safety Information Page in the front of this manual.

Contents

- “Corrosion” on page 6-2
- “Cleaning” on page 6-2
- “Storage” on page 6-3
- “Inspection” on page 6-3
- “Overspeed Decal Replacement” on page 6-3
- “Service Decontamination Policy” on page 6-4

Corrosion

The TV-860 rotor body is made from titanium, making it more resistant to corrosion than an aluminum rotor; however, it should be maintained and kept clean the same as an aluminum rotor. Proper care will lessen the chances of rotor failure and significantly prolong the useful life of the rotor.

Corrosion commonly refers to chemical reactions at the surface (that is, rusting or pitting) recognized by the growing areas of visible deterioration. On the other hand, stress corrosion attacks the inside of the metal as well; barely detectable surface cracks grow inward, weakening the part without visible warning. Stress corrosion applies to most commonly used alloys, even the corrosion-resistant alloys have been found susceptible.

Stress corrosion is thought to be initiated by certain combinations of stress and chemical reaction. The most common chemical causing harmful effects is chloride, whether in a solution such as ammonium salts or as a subtle form as hand perspiration. If the rotor is not kept clean and chemicals remain on the rotor body, corrosion will result. Also, any moisture left on the rotor for an extended period of time can initiate corrosion. Therefore, it is important to keep the rotor clean and thoroughly dry after use.

In general, conditions for corrosion are present in all rotor applications; proper care and maintenance will minimize its effects.

Cleaning

These procedures are for general cleaning purposes only. If the rotor or any components are exposed to a contaminant, they must be decontaminated first, then washed.

a. Washing

Wash the rotor body, and sealing assembly components with warm water and mild soap or detergent at least once a week or, ideally, after each use. It is particularly important that the rotor is washed immediately after any spills have occurred. Most laboratory chemicals can be removed with a lukewarm, 1% solution of a mild, non-alkaline detergent such as a mild dishwashing liquid. Rinse and dry thoroughly before storing.

Do not use strong laboratory detergents to clean the rotor surface. Use a bristle brush to loosen encrusted materials only if necessary; be careful not to scratch the rotor surface.

b. Decontamination

Ethylene oxide, a 2% glutaraldehyde solution, or ultraviolet radiation are the recommended methods of sterilizing; however, the titanium rotor body of the TV-860 can be autoclaved at temperatures up to 121°C. Do not autoclave the ULTRACRIMP[®] rotor caps or the optional sealing assembly components. If these parts are subjected to temperatures above 100°C, they should not be used.



CAUTION Do not autoclave any sealing assembly components. If any of these parts are subjected to a temperature above 100°C, they should not be used.

For general radioactive decontamination, use a solution of equal parts of 70% ethanol, 10% SDS, and water. Follow this with ethanol rinses, then deionized water rinses, and dry with a soft absorbent cloth. Dispose of all wash solutions in proper radioactive waste containers.



CAUTION Most commercially available radioactive decontamination solutions are not compatible with titanium or aluminum.

Storage

After the rotor has been cleaned and dried, it should be stored upside-down without rotor body sealing caps; this will keep the tube compartments dry.

Inspection

Periodically, inspect the rotor body and sealing assembly components for signs of: stress, including cracks, tears, or abrasions; wear (particularly the threads of the rotor body caps); corrosion or deformation. If such problems are found, contact Thermo Fisher Scientific for information on factory inspection or replacement.

Overspeed Decal Replacement

Before replacing the decal, be sure the rotor is dry and at room temperature; if it is not, the new decal will not adhere properly.

To replace the decal (Cat. No. 51346):

1. Remove the existing decal from the bottom of the rotor being careful not to scratch the rotor surface.
2. Clean the adhesive from the rotor surface using either acetone or 3M General Adhesive Remover #8984.
3. Wipe the surface dry with a clean, soft cloth.

Note Check that the new decal has the correct number of black segments the TV-860 Rotor decal should have 15 black segments.

4. Peel the paper backing off the new decal. Fit the decal into the recess in the bottom of the rotor. Be sure the decal is properly centered, then press the decal firmly into place.

Service Decontamination Policy



WARNING Because of the characteristics of the samples likely to be processed in this centrifuge, biological or radioactive contamination may occur. Always be aware of this possibility, and take normal precautions. Use appropriate decontamination procedures should exposure occur.

Observe the WARNING on hazardous materials found on the Safety Information Page in the front of this manual.

If a centrifuge or rotor that has been used with radioactive or pathogenic material requires servicing by Thermo Fisher Scientific personnel, either at the customers laboratory or at a Thermo Fisher Scientific facility, comply with the following procedure to ensure the safety of all personnel:

1. Clean the equipment to be serviced of all encrusted material and decontaminate it (see Care and Maintenance section of centrifuge or rotor instruction manual) prior to servicing by the Thermo Fisher Scientific representative or returning it to the Thermo Fisher Scientific facility. There must be no radioactivity detectable by survey equipment.

The Thermo Fisher Scientific Product Guide contains descriptions of commonly used decontamination methods and a chart showing method compatibility with various materials. The Care and Maintenance Section of the centrifuge or rotor instruction manual contains specific guidance about cleaning and decontamination methods appropriate for the product it describes.

Clean and decontaminate your centrifuge or rotor as follows:

For ultracentrifuges:

- a. Remove rotor from the rotor chamber.
- b. Decontaminate door and rotor chamber using an appropriate method.

For rotors:

Remove tubes, bottles, and adapters from the rotor and decontaminate rotor using an appropriate method. If tubes or rotor caps are stuck in the rotor, or the rotor lid is stuck, notify Thermo Fisher Scientific representative; be prepared with the name and nature of the sample so the Thermo Fisher Scientific Chemical Hazards Officer can decide whether to authorize the rotor's return to a Thermo Fisher Scientific facility.

Do not leave a loaded rotor locked inside a centrifuge that requires servicing. If, with a loaded rotor installed in the chamber, a centrifuge malfunction makes it so that the chamber door will not open by normal means, follow the Emergency Sample Recovery procedure found in your centrifuge operating instructions manual to gain access to the rotor.

2. Complete and attach Decontamination Information Certificate (in the back of your rotor or instrument manual) to the centrifuge or rotor before servicing.

Decontamination Information Certificates are included with this book. Additional certificates are available from the local Thermo Fisher Scientific Representative or Field Service Engineer. In the event these certificates are not available, a signed, written statement certifying that the unit has been properly decontaminated, identifying what the contaminants were and outlining the decontamination procedures used will be acceptable.

Note The Field Service Engineer will note on the Customer Service Repair Report if decontamination was required and, if so, what the contaminant was and what procedure was used. If no decontamination was required, it will be so stated.

If a centrifuge or rotor to be serviced does not have a Decontamination Information Certificate attached and, in Thermo Fisher Scientific's opinion presents a potential radioactive or biological hazard, the Thermo Fisher Scientific representative will not service the equipment until proper decontamination and certification is complete.

If the centrifuge or rotor must be returned to a Thermo Fisher Scientific facility:

1. Contact your Thermo Fisher Scientific representative to obtain an Equipment Return Decontamination Form; be prepared with the name and serial number of the centrifuge or rotor and the repairs required.
2. Complete the Equipment Return Decontamination Form and returned it to Thermo Fisher Scientific. Upon receipt of a completed form, a Returned Material Authorization Number (RMA Number) will be issued to you.
3. With the RMA Number clearly marked on the outside of packaging, send the items to the address obtained from your Kendro representative.



CAUTION Do not ship or transport a centrifuge with a rotor installed on the drive spindle. If a centrifuge chamber door cannot be opened using conventional methods, refer to the Emergency Sample Recovery (mechanical override) instructions that are provided in your centrifuge operating manual.

Note United States federal regulations require that parts and instruments must be decontaminated before being transported. Outside the United States, check local regulations.

If equipment is received at Thermo Fisher Scientific facilities without a valid RMA Number on the outside of the shipping container and a completed Equipment Return Decontamination Form on file, the equipment will be treated as a potential contamination hazard, and will not be serviced until decontamination certification has been completed. The sender will be contacted for instructions regarding disposition of the equipment in question; all disposition costs will be borne by the sender. If contaminated equipment is received at Thermo Fisher Scientific facilities, both the carrier and appropriate authorities shall be notified.

Chemical Compatibility Chart

CHEMICAL	MATERIAL																											
	ALUMINUM	ANODIC COATING for ALUMINUM	BUNA N	CELLULOSE ACETATE BUTYRATE	POLYURETHANE ROTOR PAINT	COMPOSITE Carbon Fiber/Epoxy	DELRIN®	ETHYLENE PROPYLENE	GLASS	NEOPRENE	NORYL®	NYLON	PET, POLYCLEAR®, CLEARCRIMP®, CCCLEARCRIMP®	POLYALLOMER	POLYCARBONATE	POLYESTER, GLASS THERMOSET	POLYETHERIMIDE	POLYRTHYLENE	POLYPROPYLENE	POLYSULFONE	POLYVINYL CHLORIDE	RULON A®, TEFLON®	SILICONE RUBBER	STAINLESS STEEL	TITANIUM	TYGON®	VITON®	
2-mercaptoethanol	S	S	U	-	S	M	S	-	S	U	S	S	U	S	S	-	S	S	S	S	U	S	S	S	S	S	S	S
Acetaldehyde	S	-	U	U	-	-	-	M	-	U	-	-	-	M	U	U	U	M	M	-	M	S	U	-	S	-	U	
Acetone	M	S	U	U	S	U	M	S	S	U	U	S	U	S	U	U	U	S	S	U	U	S	M	M	S	U	U	
Acetonitrile	S	S	U	-	S	M	S	-	S	S	U	S	U	M	U	U	-	S	M	U	U	S	S	S	S	U	U	
Alconox®	U	U	S	-	S	S	S	-	S	S	S	S	S	S	M	S	S	S	S	S	S	S	S	S	S	S	U	
Allyl Alcohol	-	-	-	U	-	-	S	-	-	-	-	S	-	S	S	M	S	S	S	S	-	M	S	-	-	S	-	-
Aluminum Chloride	U	U	S	S	S	S	U	S	S	S	S	M	S	S	S	S	-	S	S	S	S	S	M	U	U	S	S	
Formic Acid (100%)	-	S	M	U	-	-	U	-	-	-	-	U	-	S	M	U	U	S	S	-	U	S	-	U	S	-	U	
Ammonium Acetate	S	S	U	-	S	S	S	-	S	S	S	S	S	S	S	U	-	S	S	S	S	S	S	S	S	S	S	S
Ammonium Carbonate	M	S	U	S	S	S	S	S	S	S	S	S	S	S	U	U	-	S	S	S	S	S	S	M	S	S	S	
Ammonium Hydroxide (10%)	U	U	S	U	S	S	M	S	S	S	S	S	-	S	U	M	S	S	S	S	S	S	S	S	S	S	M	S
Ammonium Hydroxide (28%)	U	U	S	U	S	U	M	S	S	S	S	S	U	S	U	M	S	S	S	S	S	S	S	S	S	S	M	S
Ammonium Hydroxide (conc.)	U	U	U	U	S	U	M	S	-	S	-	S	U	S	U	U	S	S	S	-	M	S	S	S	S	-	U	
Ammonium Phosphate	U	-	S	-	S	S	S	S	S	S	S	S	-	S	S	M	-	S	S	S	S	S	S	M	S	S	S	S
Ammonium Sulfate	U	M	S	-	S	S	U	S	S	S	S	S	S	S	S	S	-	S	S	S	S	S	S	U	S	S	U	
Amyl Alcohol	S	-	M	U	-	-	S	S	-	M	-	S	-	M	S	S	S	S	M	-	-	-	U	-	S	-	M	
Aniline	S	S	U	U	S	U	S	M	S	U	U	U	U	U	U	U	-	S	M	U	U	S	S	S	S	U	S	
Sodium Hydroxide (<1%)	U	-	M	S	S	S	-	-	S	M	S	S	-	S	M	M	S	S	S	S	S	S	M	S	S	-	U	
Sodium Hydroxide (10%)	U	-	M	U	-	-	U	-	M	M	S	S	U	S	U	U	S	S	S	S	S	S	M	S	S	-	U	
Barium Salts	M	U	S	-	S	S	S	S	S	S	S	S	S	S	M	-	S	S	S	S	S	S	M	S	S	S	S	
Benzene	S	S	U	U	S	U	M	U	S	U	U	S	U	U	U	M	U	M	U	U	U	S	U	U	S	U	S	
Benzyl Alcohol	S	-	U	U	-	-	M	M	-	M	-	S	U	U	U	U	U	U	U	-	M	S	M	-	S	-	S	
Boric Acid	U	S	S	M	S	S	U	S	S	S	S	S	S	S	S	S	U	S	S	S	S	S	S	S	S	S	S	S

A Chemical Compatibility Chart

CHEMICAL	MATERIAL																										
	ALUMINUM	ANODIC COATING for ALUMINUM	BUNA N	CELLULOSE ACETATE BUTYRATE	POLYURETHANE ROTOR PAINT	COMPOSITE Carbon Fiber/Epoxy	DELRIN®	ETHYLENE PROPYLENE	GLASS	NEOPRENE	NORYL®	NYLON	PET®, POLYCLEAR®, CLEARCRIMP®, CCCCLEARCRIMP®	POLYALLUMER	POLYCARBONATE	POLYESTER, GLASS THERMOSET	POLYETHERIMIDE	POLYETHYLENE	POLYPROPYLENE	POLYSULFONE	POLYVINYL CHLORIDE	RULON A®, TEFLON®	SILICONE RUBBER	STAINLESS STEEL	TITANIUM	TYGON®	VITON®
Cesium Acetate	M	-	S	-	S	S	S	-	S	S	S	S	-	S	S	-	-	S	S	S	S	S	S	M	S	S	S
Cesium Bromide	M	S	S	-	S	S	S	-	S	S	S	S	S	S	S	-	-	S	S	S	S	S	S	M	S	S	S
Cesium Chloride	M	S	S	U	S	S	S	-	S	S	S	S	S	S	S	-	-	S	S	S	S	S	S	M	S	S	S
Cesium Formate	M	S	S	-	S	S	S	-	S	S	S	S	S	S	S	-	-	S	S	S	S	S	S	M	S	S	S
Cesium Iodide	M	S	S	-	S	S	S	-	S	S	S	S	S	S	S	-	-	S	S	S	S	S	S	M	S	S	S
Cesium Sulfate	M	S	S	-	S	S	S	-	S	S	S	S	S	S	S	-	-	S	S	S	S	S	S	M	S	S	S
Chloroform	U	U	U	U	S	S	M	U	S	U	U	M	U	M	U	U	U	M	M	U	U	S	U	U	U	M	S
Chromic Acid (10%)	U	-	U	U	S	U	U	-	S	S	S	U	S	S	M	U	M	S	S	U	M	S	M	U	S	S	S
Chromic Acid (50%)	U	-	U	U	-	U	U	-	-	-	S	U	U	S	M	U	M	S	S	U	M	S	-	U	M	-	S
Cresol Mixture	S	S	U	-	-	-	S	-	S	U	U	U	U	U	U	-	-	U	U	-	U	S	S	S	S	U	S
Cyclohexane	S	S	S	-	S	S	S	U	S	U	S	S	U	U	U	M	S	M	U	M	M	S	U	M	M	U	S
Deoxycholate	S	S	S	-	S	S	S	-	S	S	S	S	S	S	S	-	-	S	S	S	S	S	S	S	S	S	S
Distilled Water	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S
Dextran	M	S	S	S	S	S	S	-	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	M	S	S	S
Diethyl Ether	S	S	U	U	S	S	S	U	S	U	U	S	U	U	U	U	U	U	U	U	U	U	S	S	S	M	U
Diethyl Ketone	S	-	U	U	-	-	M	-	S	U	-	S	-	M	U	U	U	M	M	-	U	S	-	-	S	U	U
Diethylpyrocarbonate	S	S	U	-	S	S	S	-	S	S	U	S	U	S	U	-	-	S	S	S	M	S	S	S	S	S	S
Dimethylsulfoxide	S	S	U	U	S	S	S	-	S	U	S	S	U	S	U	U	-	S	S	U	U	S	S	S	S	U	U
Dioxane	M	S	U	U	S	S	M	M	S	U	U	S	U	M	U	U	-	M	M	M	U	S	S	S	S	U	U
Ferric Chloride	U	U	S	-	-	-	M	S	-	M	-	S	-	S	-	-	-	S	S	-	-	-	M	U	S	-	S
Acetic Acid (Glacial)	S	S	U	U	S	S	U	M	S	U	S	U	U	U	U	U	M	S	U	M	U	S	U	U	S	-	U
Acetic Acid (5%)	S	S	M	S	S	S	M	S	S	S	S	S	M	S	S	S	S	S	S	S	M	S	S	M	S	S	M
Acetic Acid (60%)	S	S	U	U	S	S	U	-	S	M	S	U	U	M	U	S	M	S	M	S	M	S	M	U	S	M	U
Ethyl Acetate	M	M	U	U	S	S	M	M	S	S	U	S	U	M	U	U	-	S	S	U	U	S	M	M	S	U	U
Ethyl Alcohol (50%)	S	S	S	S	S	S	M	S	S	S	S	S	U	S	U	S	S	S	S	S	S	S	S	M	S	M	U
Ethyl Alcohol (95%)	S	S	S	U	S	S	M	S	S	S	S	S	U	S	U	-	S	S	S	M	S	S	S	U	S	M	U
Ethylene Dichloride	S	-	U	U	-	-	S	M	-	U	U	S	U	U	U	U	U	U	U	-	U	S	U	-	S	-	S
Ethylene Glycol	S	S	S	S	S	S	S	S	S	S	S	S	-	S	U	S	S	S	S	S	S	S	S	M	S	M	S
Ethylene Oxide Vapor	S	-	U	-	-	U	-	-	S	U	-	S	-	S	M	-	-	S	S	S	U	S	U	S	S	S	U
Ficoll-Hypaque®	M	S	S	-	S	S	S	-	S	S	S	S	-	S	S	-	S	S	S	S	S	S	S	M	S	S	S

CHEMICAL	MATERIAL	ALUMINUM	ANODIC COATING for ALUMINUM	BUNA N	CELLULOSE ACETATE BUTYRATE	POLYURETHANE ROTOR PAINT	COMPOSITE Carbon Fiber/Epoxy	DELRIN®	ETHYLENE PROPYLENE	GLASS	NEOPRENE	NORYL®	NYLON	PET®, POLYCLEAR®, CLEARCRIMP®, CIRCLECRIMP®	POLYALLUMER	POLYCARBONATE	POLYESTER, GLASS THERMOSET	POLYETHERIMIDE	POLYETHYLENE	POLYPROPYLENE	POLYSULFONE	POLYVINYL CHLORIDE	RULON A®, TEFLON®	SILICONE RUBBER	STAINLESS STEEL	TITANIUM	TYGON®	VITON®
Hydrofluoric Acid (10%)	U	U	U	M	-	-	U	-	-	U	U	S	-	S	M	U	S	S	S	S	M	S	U	U	U	-	-	
Hydrofluoric Acid (50%)	U	U	U	U	-	-	U	-	-	U	U	U	U	S	U	U	U	S	S	M	M	S	U	U	U	-	M	
Hydrochloric Acid (conc.)	U	U	U	U	-	U	U	M	-	U	M	U	U	M	U	U	U	-	S	-	U	S	U	U	U	-	-	
Formaldehyde (40%)	M	M	M	S	S	S	S	M	S	S	S	S	M	S	S	S	U	S	S	M	S	S	S	M	S	M	U	
Glutaraldehyde	S	S	S	S	-	-	S	-	S	S	S	S	S	S	S	-	-	S	S	S	-	-	S	S	S	-	-	
Glycerol	M	S	S	-	S	S	S	S	S	S	S	S	S	S	S	S	-	S	S	S	S	S	S	S	S	S	S	
Guanidine Hydrochloride	U	U	S	-	S	S	S	-	S	S	S	S	S	S	S	-	-	S	S	S	S	S	S	U	S	S	S	
Haemo-Sol®	S	S	S	-	-	-	S	-	S	S	S	S	S	S	S	-	-	S	S	S	S	S	S	S	S	S	S	
Hexane	S	S	S	-	S	S	S	-	S	S	U	S	U	M	U	S	S	U	S	S	M	S	U	S	S	U	S	
Isobutyl Alcohol	-	-	M	U	-	-	S	S	-	U	-	S	U	S	S	M	S	S	S	-	S	S	S	-	S	-	S	
Isopropyl Alcohol	M	M	M	U	S	S	S	S	S	U	S	S	U	S	U	M	S	S	S	S	S	S	S	M	M	M	S	
Iodoacetic Acid	S	S	M	-	S	S	S	-	S	M	S	S	M	S	S	-	M	S	S	S	S	S	M	S	S	M	M	
Potassium Bromide	U	S	S	-	S	S	S	-	S	S	S	S	S	S	S	S	S	S	S	-	S	S	S	M	S	S	S	
Potassium Carbonate	M	U	S	S	S	S	S	-	S	S	S	S	S	S	U	S	S	S	S	S	S	S	S	S	S	S	S	
Potassium Chloride	U	S	S	-	S	S	S	S	S	S	S	S	S	S	S	-	S	S	S	S	S	S	S	U	S	S	S	
Potassium Hydroxide (5%)	U	U	S	S	S	S	M	-	S	S	S	S	-	S	U	S	S	S	S	S	S	S	M	U	M	S	U	
Potassium Hydroxide (conc.)	U	U	M	U	-	-	M	-	M	S	S	-	U	M	U	U	U	S	M	-	M	U	-	U	U	-	U	
Potassium Permanganate	S	S	S	-	S	S	S	-	S	S	S	U	S	S	S	M	-	S	M	S	U	S	S	M	S	U	S	
Calcium Chloride	M	U	S	S	S	S	S	S	S	S	S	S	S	S	M	S	-	S	S	S	S	S	M	S	S	S	S	
Calcium Hypochlorite	M	-	U	-	S	M	M	S	-	M	-	S	-	S	M	S	-	S	S	S	M	S	M	U	S	-	S	
Kerosene	S	S	S	-	S	S	S	U	S	M	U	S	U	M	M	S	-	M	M	M	S	S	U	S	S	U	S	
Sodium Chloride (10%)	S	-	S	S	S	S	S	-	-	-	-	S	S	S	S	S	-	S	S	S	S	-	S	S	M	-	S	
Sodium Chloride (sat'd)	U	-	S	U	S	S	S	-	-	-	-	S	S	S	S	S	-	S	S	-	S	-	S	S	M	-	S	
Carbon Tetrachloride	U	U	M	S	S	U	M	U	S	U	U	S	U	M	U	S	S	M	M	S	M	M	M	M	U	S	S	
Aqua Regia	U	-	U	U	-	-	U	-	-	-	-	-	U	U	U	U	U	U	U	-	-	-	-	-	S	-	M	
Solution 555 (20%)	S	S	S	-	-	-	S	-	S	S	S	S	S	S	S	-	-	S	S	S	-	S	S	S	S	S	S	
Magnesium Chloride	M	S	S	-	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	M	S	S	S	
Mercaptoacetic Acid	U	S	U	-	S	M	S	-	S	M	S	U	U	U	U	-	S	U	U	S	M	S	U	S	S	S	S	
Methyl Alcohol	S	S	S	U	S	S	M	S	S	S	S	S	U	S	U	M	S	S	S	S	S	S	S	M	S	M	U	
Methylene Chloride	U	U	U	U	M	S	S	U	S	U	U	S	U	U	U	U	U	U	M	U	U	U	S	S	M	U	U	

A Chemical Compatibility Chart

CHEMICAL	MATERIAL	ALUMINUM	ANODIC COATING for ALUMINUM	BUNA N	CELLULOSE ACETATE BUTYRATE	POLYURETHANE ROTOR PAINT	COMPOSITE Carbon Fiber/Epoxy	DELRIN®	ETHYLENE PROPYLENE	GLASS	NEOPRENE	NORYL®	NYLON	PET®, POLYCLEAR®, CLEARCRIMP®, CCCCLEARCRIMP®	POLYALLUMER	POLYCARBONATE	POLYESTER, GLASS THERMOSET	POLYETHERIMIDE	POLYETHYLENE	POLYPROPYLENE	POLYSULFONE	POLYVINYL CHLORIDE	RULON A®, TEFLON®	SILICONE RUBBER	STAINLESS STEEL	TITANIUM	TYGON®	VITON®
Methyl Ethyl Ketone		S	S	U	U	S	S	M	S	S	U	U	S	U	S	U	U	U	S	S	U	U	S	S	S	S	U	U
Metrizamide®		M	S	S	-	S	S	S	-	S	S	S	S	-	S	S	-	-	S	S	S	S	S	S	M	S	S	S
Lactic Acid (100%)		-	-	S	-	-	-	-	-	M	S	U	-	S	S	S	M	S	S	-	M	S	M	S	S	-	S	
Lactic Acid (20%)		-	-	S	S	-	-	-	-	M	S	M	-	S	S	S	S	S	S	S	M	S	M	S	S	-	S	
N-Butyl Alcohol		S	-	S	U	-	-	S	-	S	M	-	U	S	M	S	S	S	S	M	M	S	M	-	S	-	S	
N-Butyl Phthalate		S	S	U	-	S	S	S	-	S	U	U	S	U	U	M	-	U	U	S	U	S	M	M	S	U	S	
N, N-Dimethylformamide		S	S	S	U	S	M	S	-	S	S	U	S	U	S	U	U	-	S	S	U	U	S	M	S	S	U	
Sodium Borate		M	S	S	S	S	S	S	S	S	S	S	U	S	S	S	S	-	S	S	S	S	S	M	S	S	S	
Sodium Bromide		U	S	S	-	S	S	S	-	S	S	S	S	S	S	S	S	-	S	S	S	S	S	M	S	S	S	
Sodium Carbonate (2%)		M	U	S	S	S	S	S	S	S	S	S	S	S	U	S	S	S	S	S	S	S	S	S	S	S	S	
Sodium Dodecyl Sulfate		S	S	S	-	S	S	S	-	S	S	S	S	S	S	-	S	S	S	S	S	S	S	S	S	S	S	
Sodium Hypochlorite (5%)		U	U	M	S	S	M	U	S	S	M	S	S	S	M	S	S	S	S	M	S	S	M	U	S	M	S	
Sodium Iodide		M	S	S	-	S	S	S	-	S	S	S	S	S	S	S	-	-	S	S	S	S	S	M	S	S	S	
Sodium Nitrate		S	S	S	-	S	S	S	S	S	S	S	S	S	S	S	S	-	S	S	S	S	S	U	S	S	S	
Sodium Sulfate		U	S	S	-	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	M	S	S	S	
Sodium Sulfide		S	-	S	S	-	-	S	-	-	-	S	S	S	U	U	-	-	S	-	-	-	S	S	M	-	S	
Sodium Sulfite		S	S	S	-	S	S	S	M	S	S	S	S	S	S	M	-	S	S	S	S	S	S	S	S	S	S	
Nickel Salts		U	S	S	S	S	S	-	S	S	S	-	-	S	S	S	S	-	S	S	S	S	M	S	S	S	S	
Oils (Petroleum)		S	S	S	-	-	-	S	U	S	S	S	S	U	U	M	S	M	U	U	S	S	U	S	S	S	S	
Oils (Other)		S	-	S	-	-	-	S	M	S	S	S	S	U	S	S	S	S	U	S	S	S	S	-	S	M	S	
Oleic Acid		S	-	U	S	S	S	U	U	S	U	S	S	M	S	S	S	S	S	S	S	S	M	U	S	M	M	
Oxalic Acid		U	U	M	S	S	S	U	S	S	S	S	U	S	U	S	S	S	S	S	S	S	S	U	M	S	S	
Perchloric Acid (10%)		U	-	U	-	S	U	U	-	S	M	M	-	-	M	U	M	S	M	M	-	M	S	U	-	S	-	
Perchloric Acid (70%)		U	U	U	-	-	U	U	-	S	U	M	U	U	M	U	U	U	M	M	U	M	S	U	U	S	U	
Phenol (5%)		U	S	U	-	S	M	M	-	S	U	M	U	U	S	U	M	S	M	S	U	U	S	U	M	M	S	
Phenol (50%)		U	S	U	-	S	U	M	-	S	U	M	U	U	U	U	U	S	U	M	U	U	S	U	U	M	S	
Phosphoric Acid (10%)		U	U	M	S	S	S	U	S	S	S	S	U	-	S	S	S	S	S	S	S	S	U	M	U	S	S	
Phosphoric Acid (conc.)		U	U	M	M	-	-	U	S	-	M	S	U	U	M	M	S	S	M	S	M	S	U	M	U	-	S	
Physiologic Media (Serum, Urine)		M	S	S	S	-	-	S	-	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	
Picric Acid		S	S	U	-	S	M	S	S	S	M	S	U	S	S	S	U	S	S	S	S	U	S	U	M	S	M	

CHEMICAL	MATERIAL																										
	ALUMINUM	ANODIC COATING for ALUMINUM	BUNA N	CELLULOSE ACETATE BUTYRATE	POLYURETHANE ROTOR PAINT	COMPOSITE Carbon Fiber/Epoxy	DELRIN®	ETHYLENE PROPYLENE	GLASS	NEOPRENE	NORYL®	NYLON	PET®, POLYCLEAR®, CLEARCRIMP®, CIRCLECRIMP®	POLYALLUMER	POLYCARBONATE	POLYESTER, GLASS THERMOSET	POLYETHERIMIDE	POLYETHYLENE	POLYPROPYLENE	POLYSULFONE	POLYVINYL CHLORIDE	RULON A®, TEFLON®	SILICONE RUBBER	STAINLESS STEEL	TITANIUM	TYGON®	VITON®
Pyridine (50%)	U	S	U	U	S	U	U	-	U	S	S	U	U	M	U	U	-	U	S	M	U	S	S	U	U	U	U
Rubidium Bromide	M	S	S	-	S	S	S	-	S	S	S	S	S	S	S	-	-	S	S	S	S	S	S	M	S	S	S
Rubidium Chloride	M	S	S	-	S	S	S	-	S	S	S	S	S	S	S	-	-	S	S	S	S	S	S	M	S	S	S
Sucrose	M	S	S	-	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S
Sucrose, Alkaline	M	S	S	-	S	S	S	-	S	S	S	S	S	S	U	S	S	S	S	S	S	S	S	M	S	S	S
Sulfosalicylic Acid	U	U	S	S	S	S	S	-	S	S	S	U	S	S	S	-	S	S	S	-	S	S	S	U	S	S	S
Nitric Acid (10%)	U	S	U	S	S	U	U	-	S	U	S	U	-	S	S	S	S	S	S	S	S	S	M	S	S	S	S
Nitric Acid (50%)	U	S	U	M	S	U	U	-	S	U	S	U	U	M	M	U	M	M	M	S	S	S	U	S	S	M	S
Nitric Acid (95%)	U	-	U	U	-	U	U	-	-	U	U	U	U	M	U	U	U	U	M	U	U	S	U	S	S	-	S
Hydrochloric Acid (10%)	U	U	M	S	S	S	U	-	S	S	S	U	U	S	U	S	S	S	S	S	S	S	S	U	M	S	S
Hydrochloric Acid (50%)	U	U	U	U	S	U	U	-	S	M	S	U	U	M	U	U	S	S	S	S	M	S	M	U	U	M	M
Sulfuric Acid (10%)	M	U	U	S	S	U	U	-	S	S	M	U	S	S	S	S	S	S	S	S	S	S	U	U	U	S	S
Sulfuric Acid (50%)	M	U	U	U	S	U	U	-	S	S	M	U	U	S	U	U	M	S	S	S	S	S	U	U	U	M	S
Sulfuric Acid (conc.)	M	U	U	U	-	U	U	M	-	-	M	U	U	S	U	U	U	M	S	U	M	S	U	U	U	-	S
Stearic Acid	S	-	S	-	-	-	S	M	S	S	S	S	-	S	S	S	S	S	S	S	S	S	M	M	S	S	S
Tetrahydrofuran	S	S	U	U	S	U	U	M	S	U	U	S	U	U	U	-	M	U	U	U	U	S	U	S	S	U	U
Toluene	S	S	U	U	S	S	M	U	S	U	U	S	U	U	U	S	U	M	U	U	U	S	U	S	U	U	M
Trichloroacetic Acid	U	U	U	-	S	S	U	M	S	U	S	U	U	S	M	-	M	S	S	U	U	S	U	U	U	M	U
Trichloroethane	S	-	U	-	-	-	M	U	-	U	-	S	U	U	U	U	U	U	U	U	U	S	U	-	S	-	S
Trichloroethylene	-	-	U	U	-	-	-	U	-	U	-	S	U	U	U	U	U	U	U	U	U	S	U	-	U	-	S
Trisodium Phosphate	-	-	-	S	-	-	M	-	-	-	-	-	-	S	-	-	S	S	S	-	-	S	-	-	S	-	S
Tris Buffer (neutral pH)	U	S	S	S	S	S	S	-	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S
Triton X-100®	S	S	S	-	S	S	S	-	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S
Urea	S	-	U	S	S	S	S	-	-	-	-	S	S	S	M	S	S	S	S	-	S	S	S	M	S	-	S
Hydrogen Peroxide (10%)	U	U	M	S	S	U	U	-	S	S	S	U	S	S	S	M	U	S	S	S	S	S	S	M	S	U	S
Hydrogen Peroxide (3%)	S	M	S	S	S	-	S	-	S	S	S	S	S	S	S	S	M	S	S	S	S	S	S	S	S	S	S
Xylene	S	S	U	S	S	S	M	U	S	U	U	U	U	U	U	M	U	M	U	U	U	S	U	M	S	U	S
Zinc Chloride	U	U	S	S	S	S	U	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S
Zinc Sulfate	U	S	S	-	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S	S
Citric Acid (10%)	M	S	S	M	S	S	M	S	S	S	S	S	S	S	S	S	M	S	S	S	S	S	S	S	S	S	S

A Chemical Compatibility Chart

*Polyethyleneterephthalate

Key

S Satisfactory

M = Moderate attack, may be satisfactory for use in centrifuge depending on length of exposure, speed involved, etc.; suggest testing under actual conditions of use.

U Unsatisfactory, not recommended.

-- Performance unknown; suggest testing, using sample to avoid loss of valuable material.

Chemical resistance data is included only as a guide to product use. Because no organized chemical resistance data exists for materials under the stress of centrifugation, when in doubt we recommend pretesting sample lots.

Contact Information

United States	866-9-THERMO
	+1 866 984 3766

Canada	+1 866 984 3766
--------	-----------------

Austria	+43 1 801 400
---------	---------------

Belgium	+32 2 482 30 30
---------	-----------------

Germany	08001 536 376
	+49 6184 90 6940

France	+33 2 2803 2180
	+33 2 2803 2000

Italy	+39 02 02 95059 341
-------	---------------------

Netherlands	+31 76 571 4440
-------------	-----------------

Nordic / Baltic Countries	+35 89 329 100
---------------------------	----------------

Russia	+7 (812) 703 42 15
--------	--------------------

B Contact Information

Spain +34 932 23 09 18

Portugal +34 932 23 09 18

Switzerland +41 44 454 12 12

UK / Ireland +44 870 609 9203

China +86 21 6865 4588

+86 10 8419 3588

India +91 22 6716 2200

Japan +81 45 453 9220

Other Asian Countries +852 2885 4613

Latin America +1 866 984 3766

Other Countries +49 6184 90 6940

+33 2 2803 2180
