



6. Spare Parts

Sorvall:

ST 8 - series

Thermo Scientific:

SL 8 - series,
Megafuge 8 – series
(formerly known as Heraeus Megafuge 8 – series)

Fisher Scientific:

AccuSpin 24C

VWR:

MegaStar600

ThermoFisher
SCIENTIFIC



ThermoFisher
SCIENTIFIC
The world leader in serving science



Overview

- [Front Panel](#) → [Required Tools](#)
- [Front View](#) → [Required Tools](#)
- [Back View](#) → [Required Tools](#)
- [Top View](#) → [Required Tools](#)
- [Drive](#) → [Required Tools](#)



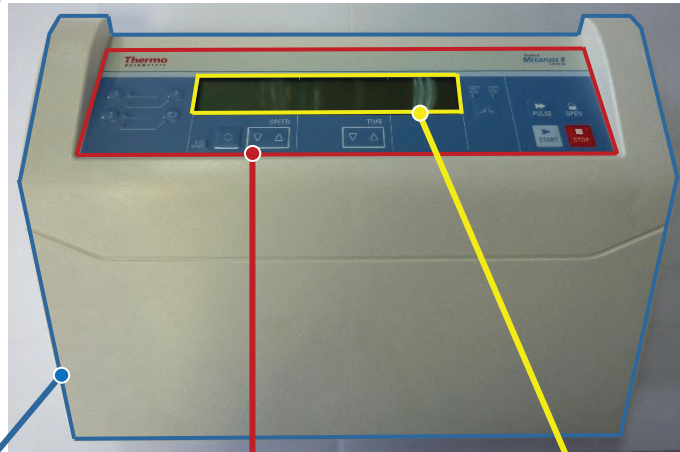
6.1

19.07.2022

Megafuge 8 – series, ST 8 -, SL 8 -, accuSpin 24C, MegaStar600

ThermoFisher
S C I E N T I F I C

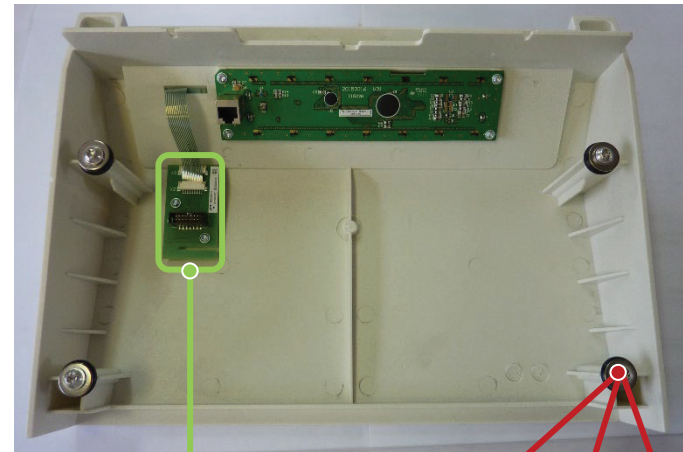
Front Panel



20059014
Front Panel

Keypad Overlay

20150316
Display



20150285
Keypad Adaptor

20290706
O-Ring

20480320
Screw

20480317
Washer

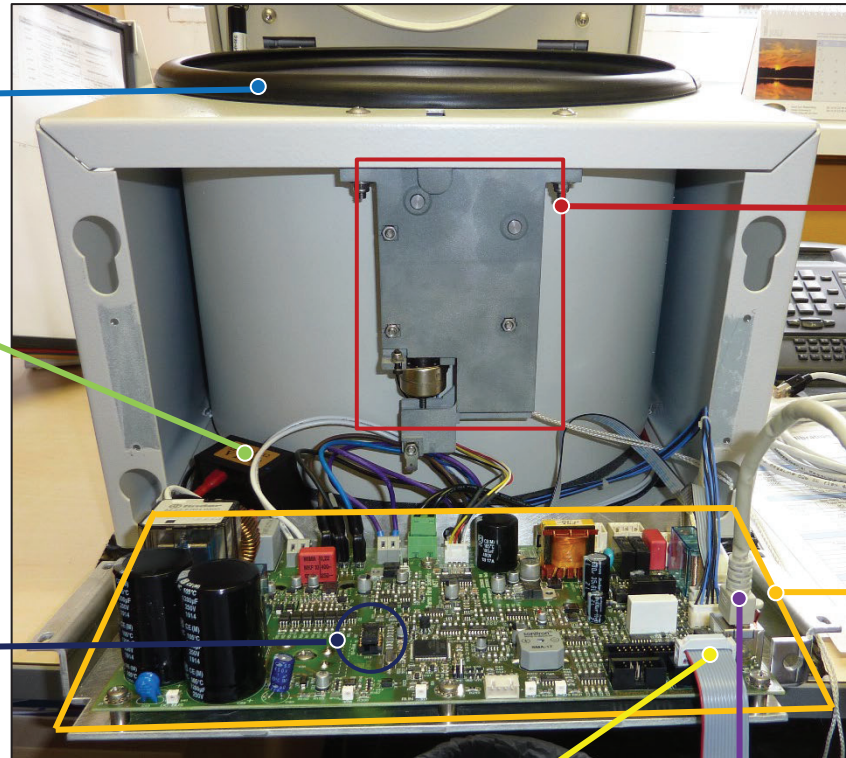
Version	Part number
Thermo Megafuge 8	20059701
Thermo SL 8	20059705
Sorvall ST 8	20059703
Fisher	20059707
VWR	20059801

6.2

Front Panel: Required Tools for Disassembly

Part Nr.:	Description	Required Tools	Additional Recommended Tools
20059014	Front Panel	4mm Allen	
20150316	Display	4mm Allen, Torx 15	
20059701	Keypad Overlay Thermo Megafuge 8	4mm Allen	
20059705	Keypad Overlay Thermo SL 8		
20059703	Keypad Overlay Sorvall ST 8		
20059707	Keypad Overlay Fisher	4mm Allen	
20150285	Keypad Adaptor	4mm Allen, Torx 15	
20290706	O-Ring	4mm Allen, Torx 30	
20480320	Screw		
20480317	Washer		

Front View



20058389
Gasket

200150324
Choke

70903020
Lid Latch

20150319
Main PCB

20059094
Keypad Cable

20059093
Display Cable

NV-RAM

Version	Part number
230V	70904751
120V	70904753
100V	70904755
120V Fisher 230V Fisher	70904783 70904789
230V VWR 120V VWR	70904784 70900172

6.4

Front View: Required Tools for Disassembly

Part Nr.:	Description	Required Tools	Additional Recommended Tools
20058389	Gasket		
70903020	Lid Latch	4mm Allen	
20059164	Choke		
20150319	Main PCB	4mm Allen	small screwdriver
20059093	Display Cable	4mm Allen	
20059094	Keypad Cable	4mm Allen	
70904751	NV-RAM 230V	4mm Allen	Small screwdriver
70904753	NV-RAM 120V		Small screwdriver
70904755	NV-RAM 100V		Small screwdriver
70904783	NV-RAM 120V Fisher		Small screwdriver
70904789	NV-RAM 230V Fisher		Small screwdriver
70904784	NV-RAM 230V VWR		Small screwdriver
70900172	NV-RAM 120V VWR		Small screwdriver

6.5

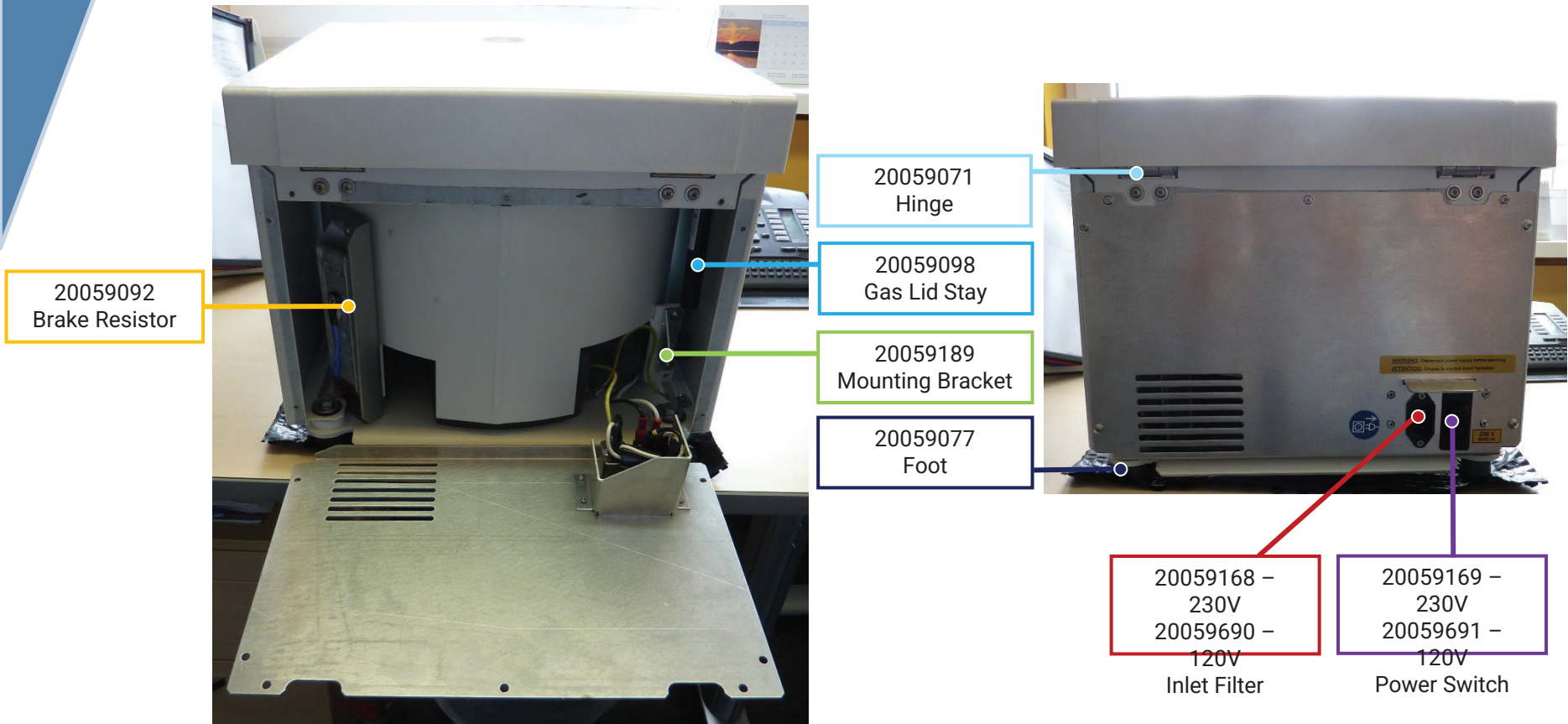
19.07.2022

Megafuge 8 – series, ST 8 -, SL 8 -, accuSpin 24C, MegaStar600

[Overview](#)

ThermoFisher
S C I E N T I F I C

Back View



6.6

19.07.2022

MegaFuge 8 – series, ST 8 -, SL 8 -, accuSpin 24C, MegaStar600

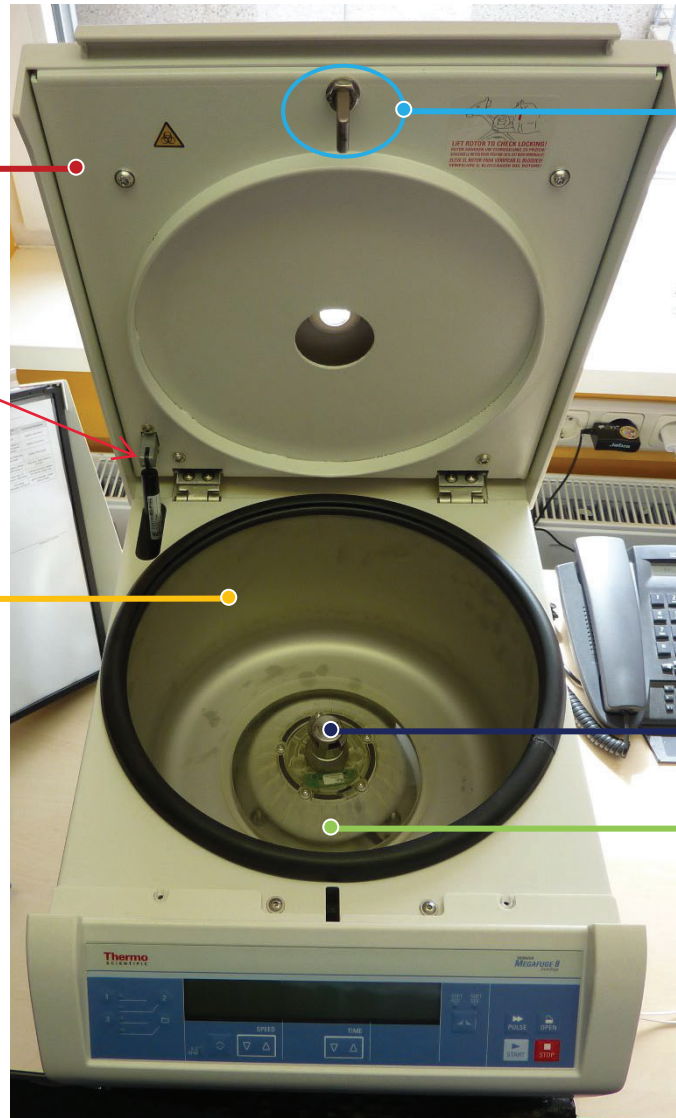
[Overview](#)

ThermoFisher
SCIENTIFIC

Back View: Required Tools for Disassembly

Part Nr.	Description	Required Tools	Additional Recommended Tools
20059071	Hinge	4mm Allen, Torx 30	striker pin, hammer, pipe wrench
20059169	Power Switch 230V	4mm Allen	small screwdriver
20059691	Power Switch 120V	4mm Allen	small screwdriver
20059077	Foot	4mm Allen, 13mm jaw wrench	striker pin, hammer, pipe wrench, small screwdriver, small side cutter
20059168	Inlet Filter 230V	4mm Allen, 8mm socket wrench	
20059690	Inlet Filter 120V	4mm Allen, 8mm socket wrench	
20059098	Gas Lid Stay		striker pin, hammer, pipe wrench
20059189	Mounting Bracket	4mm Allen, 32mm jaw wrench, Torx 30, 10mm socket wrench	striker pin, hammer, pipe wrench
20059092	Brake Resistor	4mm Allen	

Top View



70903021
Lid

20053064
Lid Hook

Gas Spring Holder = 20056650

20059004
Bowl

50140042
AutoLock

70903034
Sensor Cap

6.8

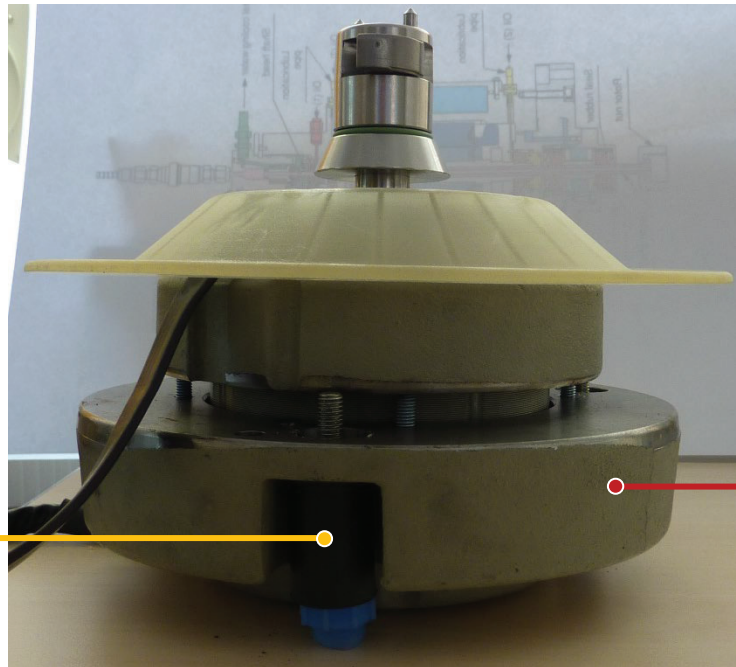
19.07.2022

Megafuge 8 – series, ST 8 -, SL 8 -, accuSpin 24C, MegaStar600

Top View: Required Tools for Disassembly

Part Nr.:	Description	Required Tools	Additional Recommended Tools
70903021	Lid	32mm jaw wrench, Torx 30, 4mm Allen	
20053064	Lid Hook	17mm jaw wrench, pipe wrench	
70903034	Sensor Cap	32mm jaw wrench, Torx 30	
20059004	Bowl	Philips 2	
50140042	AutoLock	AutoLock Tool Kit	

Drive



70902001
3x sAV-mount
incl. 6x sand-
paper discs

20059030
Motor

6.10

19.07.2022

Megafuge 8 – series, ST 8 -, SL 8 -, accuSpin 24C, MegaStar600

[Overview](#)

ThermoFisher
SCIENTIFIC

Drive: Required Tools for Disassembly

Part Nr.:	Description	Required Tools	Additional Recommended Tools
20059030	Motor	AutoLock Tool Kit, 3mm Allen, Torx 20, 13mm socket wrench	small side cutter
	Service Kit AV-Mount	13mm socket & jaw wrench, Torx 20	AutoLock Tool Kit