

Thermo Scientific

TZ-32 Zonal Rotor

for the Thermo Scientific WX+ Centrifuge

Instruction Manual

50146926-a • 01 / 2015

Table of Contents

1. Preface	4
1.1. Intended Use	4
1.2. Items Supplied.	5
1.2.1. Rotor Assembly	5
1.2.2. Seal Attachment Assembly	7
1.3. Precautions	9
1.4. Symbols used in the manual	9
2. Technical Specifications	10
3. Construction	11
3.1. High-Speed Rotation	11
3.2. Loading and Unloading	11
4. Assembly	12
4.1. Rotor Assembly	12
4.1.3. Rotor body and cover	12
4.1.4. Septa	12
4.1.5. Shaft	13
4.1.6. Stopper	13
4.1.7. Rulon seal	13
4.1.8. Cap assembly	13
4.2. RPZ-S Seal Attachment Assembly	14
4.2.9. Seal assembly	14
4.2.10. Guard plate assembly	16
4.2.11. Shield plate	16
5. Operation	17
5.1. General Description	20
5.2. Preparation	21
5.2.1. Prepare the Centrifuge	21
5.2.2. Connect a tube to the seal assembly	25
5.2.3. Confirm test operation	26
5.3. Operation	27
5.3.1. Prepare the Rotor	27

5.3.2.	Preparation for spinning the rotor	30
5.3.3.	Spinning.	31
5.3.4.	Injection of gradient solution and sample.	31
5.3.5.	High speed spinning.	35
5.3.6.	Deceleration and extraction.	36
5.4.	Operational Checkpoints.	37
5.4.1.	Setting of sample temperature and rotor temperature	37
5.4.2.	Precautions on loading and unloading	37
5.5.	Capacity in Rotor.	39
5.6.	Separation Characteristic	39
5.7.	Rotor Life	39
6.	Maintenance	40
6.1.	Cleaning of Rotor and Septa.	40
6.2.	Sterilization of Rotor	40
6.3.	Inspection of Rotors	41
6.4.	Handling of Rulon Seal and Renewal.	42
6.5.	Disassembling and Cleaning of Cap Assembly.	43
6.6.	Disassembling and Cleaning of Seal Assembly	43
6.7.	Cleaning of Ball Bearing	44
6.8.	Consumable Parts of the Rotor	45
6.9.	Service Parts List for the Cap Assembly	45
6.10.	Service Parts List for the Seal Assembly.	46
7.	Troubleshooting	47
7.1.	Check the Rulon Seal	48
7.2.	Other Check Points	49

Table of Figures

Rotor Components	8
Seal attachment components (Type RPZ-S. Part cord. 901306)	10
Construction – High Speed Rotation	14
Construction –Loading and Unloading.	14
Rotor body and cover.	15
Septa	15
Shaft, Stopper and Rulon Seal	16
Cap assembly	16
Seal assembly	17
Construction of the seal assembly	18
Guard plate assembly	19
Shield plate	19
Overspeed decal View from the bottom.	20
Mounting the supports.	24
Centering the guard plate	25
Loosening the lock nut.	25
Adjusting the base.	26
Height adjustment	26
Tightening the three setscrews.	27
Tube connection	28
Setting the seal packing in detail	30
Setting the seal packing.	31
Assembling the Rulon seal and O-ring on the shaft.	31
Shaft into septa.	32
Assembled septa into rotor	32
Assemble the rotor cover	33
Mount the guard plate	33
Mount the seal assembly	34
Mount the shield plate	35
Feed a gradient solution by pump	35
Injection of sample and overlay	36
How to avoid bubbles while injecting a density gradient solution and sample	36
Preparation of the pump	37
Removal of air bubbles	37
Place the cap	38
Checking the cap	38
Deceleration and extraction	39
Relation between volume and radius	42
Removal of the metal ring	43
Rotor inspection points	44
Polishing the Rulon seal.	45
Removal of the Rulon seal	45
Parts of the cap assembly	46
Construction of the seal assembly	47
Service parts list for the cap assembly	48
Service parts list for the seal assembly.	49

Table of Tables

Rotor Accessories	8
Seal Attachment Accessories	10
Technical Specifications	12
Employed materials	20
Chemical Resistance of Noryl	21
Separation Procedure	22
Operation modes	33
Increase of Volume	39
Separation characteristic	41
Rotor Life	41
Sterilization methods	43
Troubleshooting	49
Defective phenomena and corrective action sealed sections	50

1. Preface

Zonal centrifugation has drawn attention in the biological, medical, and agricultural fields. Swinging bucket rotors have mainly been used for separating cellular substances, virus, and other fine particles. The advantages of the zonal rotor over the swinging bucket rotors exist in reduction of convective disturbance due to “wall effect”, large capacity, and capability of dynamic continuous loading, unloading and analytical operation of samples while the rotor is spinning. Because of the high performance and reliability, the TZ-32 Zonal rotor can be used in many fields of laboratory work and commercial production.

1.1. Intended Use

The rotor is used in combination with the according centrifuge as a laboratory product designed to separate components by generation of Relative Centrifugal Force (RCF).

If these rotors are used in a manner not specified by the manufacturer, the protection provided by the equipment may be impaired.

These rotors should be operated by trained specialists only.

1.2. Items Supplied

1.2.1. Rotor Assembly

Rotor Components







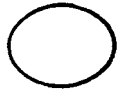







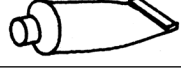
Figure 1
Rotor
Components



CAUTION The TZ-32 zonal rotor assembly does not include the RPZ-S seal attachments assembly. It has to be ordered separately. See "Figure 2" on page 7.

Rotor Accessories

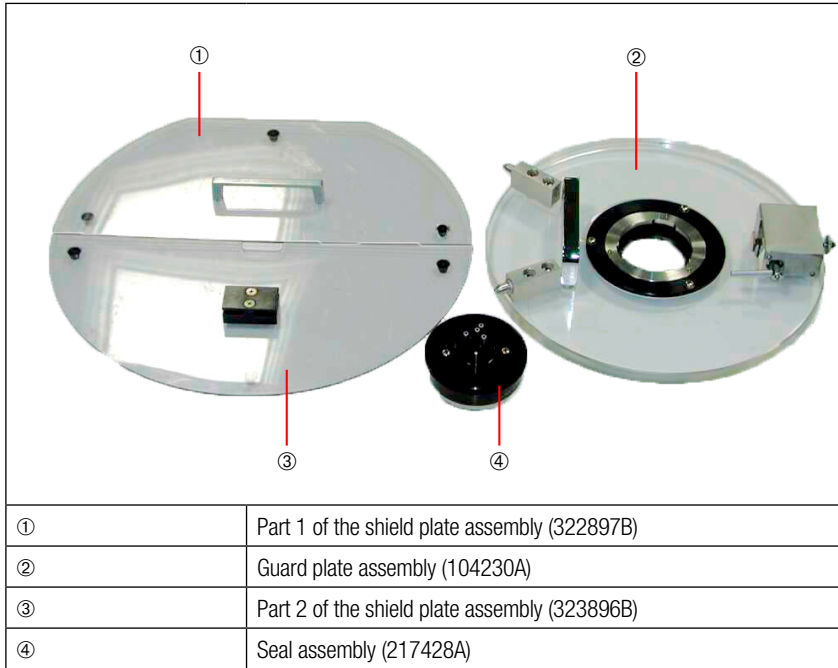
Table 1
Rotor Accessories

Item	Article No.	Note	Quantity	Picture
Rulon seal	439345		3	
Rotor stand	216715		1	
Seal packing	451153	TZ-32	1	
O-ring	8062222	1AS22 (JIS-B-2401) For rotor	2	
O-ring	8062014	1AP14 (JIS-B-2401) For Rulon seal	2	
O-ring	8062003	1AP3 (JIS-B-2401) For Rulon seal	2	
O-ring	8062018	1AP18 (JIS-B-2401) For shaft	6	
O-ring	8062218	1AS18 (JIS-B-2401) For shaft	2	
Stopper	455897		2	
Instruction manual	50146926		1	
Over speed decal	34644414		2	
Screwdriver	464912		1	
Tool box	320717		1	
Circlip remover	84850103		1	
Lubricant for screws	84810201		1	

1.2.2. Seal Attachment Assembly









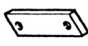
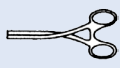



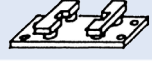

Seal Attachment Components

Figure 2
Seal attachment
components
(Type RPZ-S. Part
cord. 901306)



Seal Attachment Accessories

Table 2
Seal Attachment
Accessories

Item	Article No.	Note	Quantity	Picture
Hexagon socket head bolt	70017308	M6 x 30	6	
D6 spring washer	80112078	D6	6	
D6 flat washer	80202088	D6	6	
Tee connector	484286		4	
Rubber tube clamp	466544		6	
Tygon tube	662905	R-3603 1/8" x 1/4" x 5 m	1	
Ball bearing	60000178		1	
Support	S408766		3	
Spacer	S408765		3	
Extractor	479965	For Rulon seal	1	
Seal rubber	S202661		1	
Snap ring remover	84850109		1	
Hexagon bar wrench	60000122	For M4 setscrew	1	
Holder base	343920A		1	
Handle	323663A		1	

1.3. Precautions

- Do not spin a rotor beyond the allowable speed.
 - » The maximum speed marked on a rotor refers to an allowable one when the maximum density of the sample or density gradient solution is less than 1.2.
 - » When a sample or density gradient solution of more than 1.2 in maximum density is used, the maximum speed must be reduced according to the instruction manual for the rotor.
 - » Upon reaching the end of the service life of a rotor specified in its instruction manual, the allowable speed should be reduced by 10%.
- Set a rotor gently and securely onto the drive shaft hub.
 - » Practice setting repeatedly to acquire the knack.
- Confirm that the rotor is free from corrosion and scores.
 - » If white corrosion is recognized on a tube hole bottom of an angle rotor, or a hole bottom of a bucket for swing rotor, immediately stop using it and contact the nearest agent of Thermo Fisher Scientific
 - » When operation is over, store the rotor after cleaning and drying.
- Make it a practice to keep a logbook for rotors.
 - » Keeping a logbook for rotors is indispensable for controlling the life or when the warranty clause is involved. Keep the logbook up to date and store it properly.
- Do not rotate this rotor at a speed of more than 5000 rpm unless all components (rotor assembly, cap assembly) are assembled.

Do not rotate this rotor at a speed of more than 5000 rpm unless the rotor is filled with liquid.

If the above cautions are not observed, septa could be damaged.
- Always ground the ultracentrifuge before use.

1.4. Symbols used in the manual



This symbol refers to general hazards.

CAUTION means that material damage could occur.

WARNING means that injuries or material damage or contamination could occur.



This symbol refers to biological hazards.

Obey the safety information contained in the instruction manual to keep you and your environment safe.

2. Technical Specifications

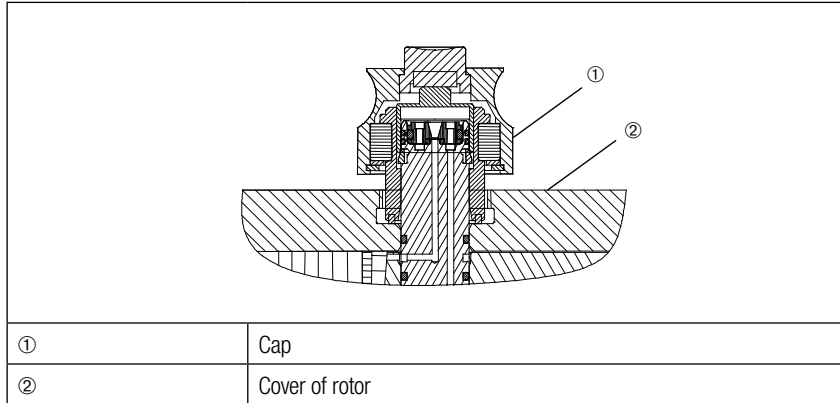
Table 3
Technical
Specifications

Maximum Speed	32000 rpm
Maximum Centrifugal Force	102000 x g
K-factor	363
Capacity	1690 ml
Maximum Density of Gradient Solution	1.2 g/ml
Maximum Radius	8.89 cm
Rotor Material	Titanium Alloy
Weight	15.4 kg

3. Construction

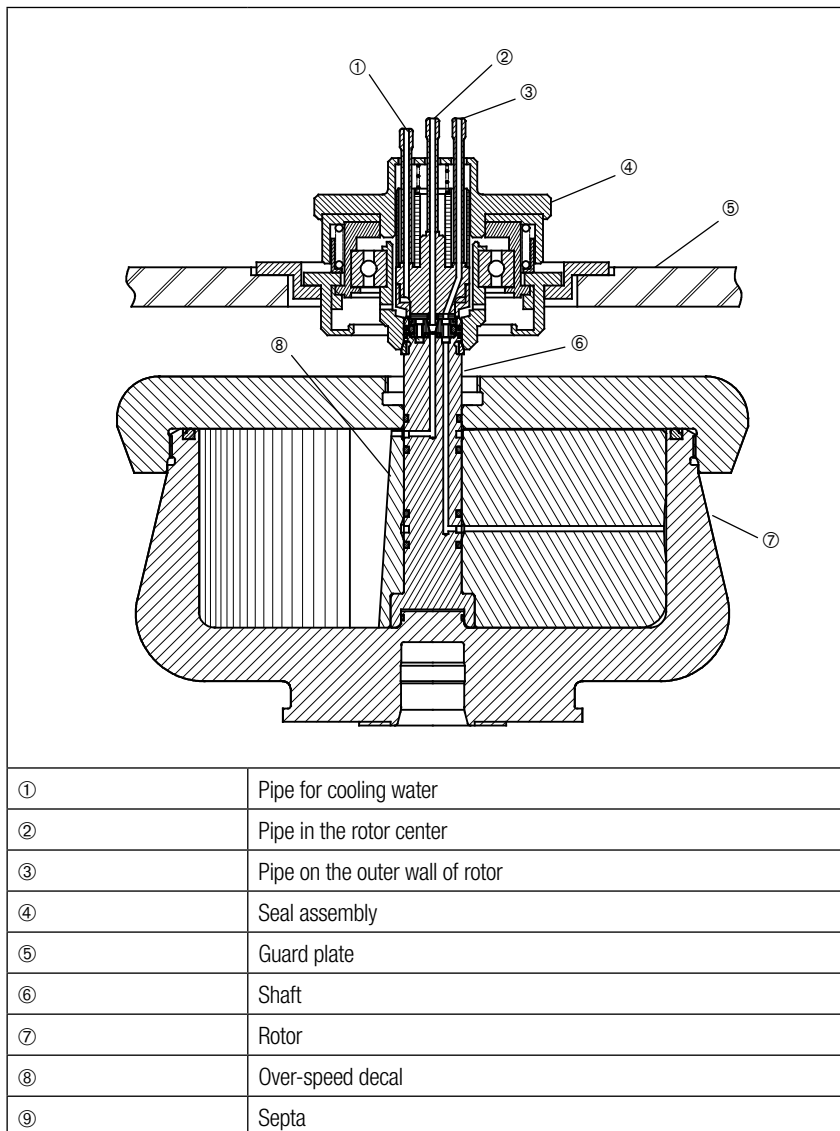
3.1. High-Speed Rotation

Figure 3
Construction –
High Speed
Rotation



3.2. Loading and Unloading

Figure 4
Construction –
Loading and
Unloading



4. Assembly

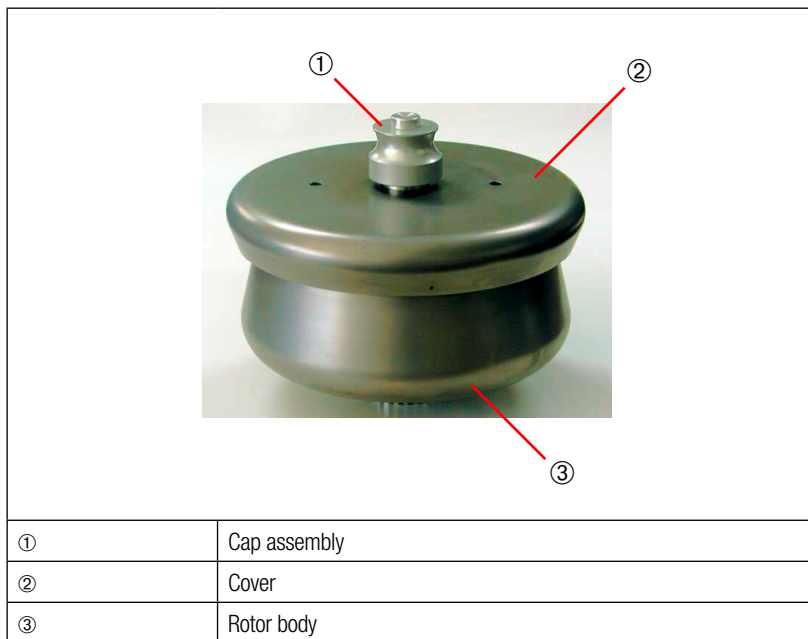
4.1. Rotor Assembly

The rotor is composed of a rotor body, cover, septa, shaft, cap assembly and Rulon™ seal attached to the shaft. The TZ-32 rotor has a capacity of 1690 ml.

4.1.3. Rotor body and cover

The rotor body and cover of the TZ-32 rotor use, as material, titanium alloy of an excellent corrosion resistance. The seal packing is fluorine rubber. The metal ring is titanium alloy.

Figure 5
Rotor body and cover



4.1.4. Septa

The septa has four vanes that divide the bowl into four sector-shaped compartments, each vane containing a channel to carry fluid to the outer edge of the rotor bowl. The solution injected through the seal assembly is led to the outer edge of the rotor from the four holes located outside the Rulon seal through the shaft. The septa has slanting surfaces, the top of each being provided with a hole communicating from the center of the seal assembly with the hole in the center of the Rulon seal. Otherwise, the septa is made of a plastic (Noryl) for which the chemical resistance, cautions in spinning, and the maintenance will be described later.

Figure 6
Septa



4. 1. 5. Shaft

A shaft is inserted in the central part of septa, and form a sample passage way. And, the tops of the shaft become the installation of the cap and the connection part of seal assembly. The material of the shaft uses titanium alloy.

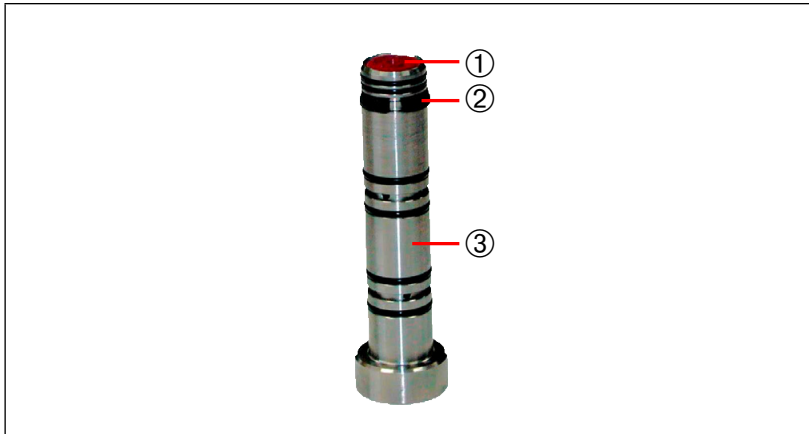
4. 1. 6. Stopper

A stopper is the coming off stop part when a cap is installed, and material is Noryl.

4. 1. 7. Rulon seal

The Rulon seal mouted on the top of shaft. This part is important component in loading and unloading, and material is fluorine plastic.

Figure 7
Shaft, Stopper and
Rulon Seal



4. 1. 8. Cap assembly

The cap is mounted onto the shaft with the seal assembly removed after injecting a sample into the rotor. It seals the sample in the rotor when evacuating the rotor chamber and spinning the rotor at a high speed.

The cap can be easily and securely removed by pushing the extruded portion (push button) atop the cap.

Figure 8
Cap assembly



4.2. RPZ-S Seal Attachment Assembly

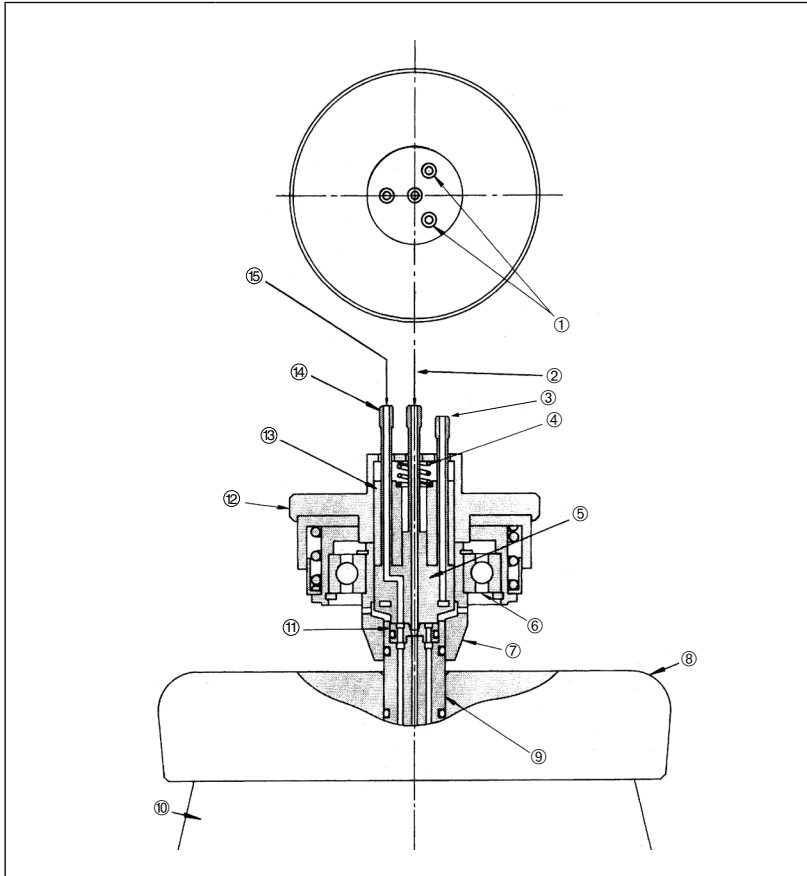
4.2.9. Seal assembly

The seal assembly is mounted onto the rotor shaft as shown in [Figure 10](#). The rotor, shaft, and adapter rotate, but the other components do not rotate. The outer side of the holder cover is supported by the guard plate which will be described later. The solution injected through the pipe is introduced through the body and Rulon seal into the rotor being spinned. One of the two higher pipes projecting from the body, which is located on the outer wall side of the rotor, is led to the outer groove of the Rulon seal and the outer wall of the rotor from the four septa vanes through the four holes of the Rulon seal. The pipe in the rotor center is led to the four holes on top of the septa through the center of the shaft. The two lower pipes are used for circulating cooling water into the interior of the body.

Figure 9
Seal assembly



Figure 10
Construction of the
seal assembly

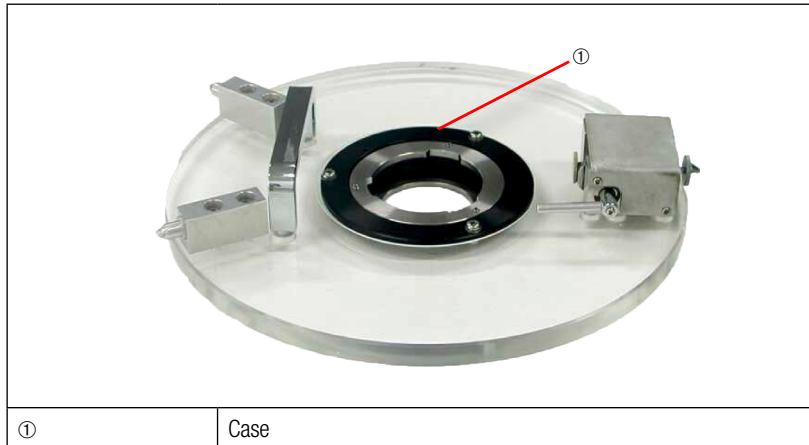


①	Body cooling water pipes
②	Pipe in the rotor center
③	Cooling water pipe
④	Coil spring (B)
⑤	Body
⑥	Seal Assembly
⑦	Ball bearing
⑧	Adaptor
⑨	Cover
⑩	Shaft
⑪	Rotor body
⑪	Rulon seal
⑫	Holder cover
⑬	Guide body
⑭	Pipe
⑮	Pipe on the outer wall side of rotor

4. 2. 10. Guard plate assembly

The center case supports the seal assembly and maintains the contact between the body and the Rulon seal for injecting solution into and extracting it, from the spinning rotor. It is mounted by using the three supports leading from the rotor chamber of the centrifuge. It is made of a transparent acrylic resin plate for the spinning rotor to be visible. It also prevents tools, etc. from dropping into the rotor chamber while it is spinning.

Figure 11
Guard plate assembly



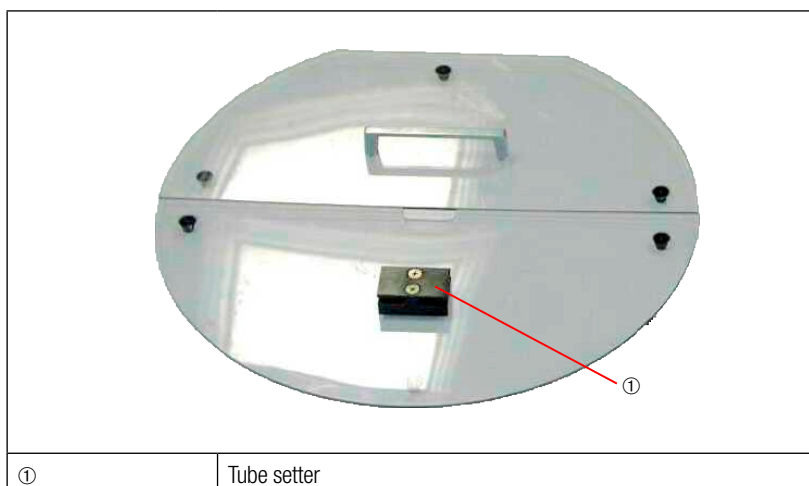
4. 2. 11. Shield plate

It is mounted on top of the rotor chamber to suppress frosting in the chamber when injecting gradient solution or samples into or extracting them from the rotor which is being cooled.


Since the shield plate is also made of a transparent acrylic resin plate, the rotor chamber is clearly visible through this plate.


A tube setter is provided for retaining the tubes projecting from the seal assembly. Since the shield plate is divided into two semicircular portions, it is easily mountable or removable.

Figure 12
Shield plate



5. Operation

	CAUTION
<p>Never spin the rotor at a speed over 5,000 rpm unless it is filled with the liquid. If this is neglected, the septa will be damaged.</p>	

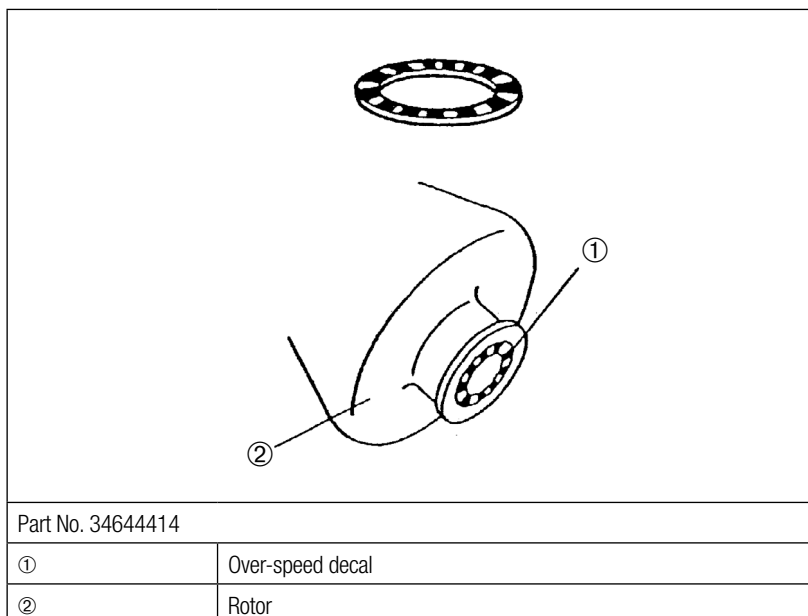
	CAUTION
<p>The over-speed decal, if corroded or discolored must be replaced immediately. The over speed decal is important to control the rotor revolution speed. The over-speed decal must match the maximum speed of the rotor. When replacing be very careful not to damage the over-speed decal and rotor body.</p>	

Replacing over-speed decal

To replace the over-speed decal follow this procedure:

1. Preparation
 - » Prepare a new over-speed decal and knife.
 - » Wash and then dry the rotor well.
2. Hold the rotor upside down.
3. Pry the edge of the over-speed decal with the knife and remove the decal.
Be very careful not to damage the rotor.
4. Clean the decal hole on the rotor by using alcohol.
5. Remove the backing paper from the new over-speed decal.
Place the over-speed decal so that it fits into the groove of the decal hole on the rotor.
Make sure the decal does not move.

Figure 13
Overspeed decal
View from the
bottom



Maximum Speed

The maximum speed of the rotor inscribed on the rotor cover surface is determined using the maximum density of gradient solution below 1.2 g/ml.

If the maximum density of gradient solution exceeds 1.2 g/ml, decrease the maximum speed of the rotor according to the following equation:

$$\text{Allowable maximum speed} = \text{Maximum speed} \times \sqrt{\frac{1.2}{\text{Maximum density of gradient solution (g/ml)}}}$$

Employed Materials

The following materials are used on the passage of the sample:

Table 4
Employed materials

Part name	Material
Tygon tube	PVC (specially processed)
Tee	Polypropylene
Metallic section of seal attachment assembly etc.	Stainless steel
Stationary seal surface (seal assembly side)	Tungsten carbide
Rulon seal	Tetrafluorethylene resin (Rulon)
O-Ring	NBR
Rotor, cover, shaft, metal ring	Titanium alloy
Septa	Noryl (polyphenylene oxide resin)
Seal packing	Fluorine rubber

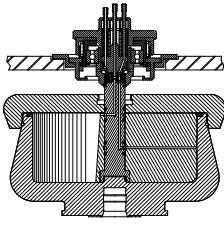
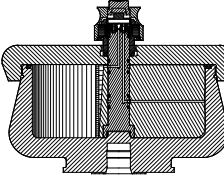
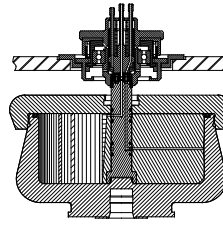
Table 5
Chemical
Resistance of Noryl

CAUTION		
Do not use the following substances with Noryl		
Benzene	Fluorine	Petroleum ether
Carbon tetrachloride	Freon	Polypropylene glycol
Chlorine	Furfural	Pyridine
Chloroform	Heptane	Pyridine sulfate
Dibutyl phthalate	Hexanol	Toluene (toluol)
Ethylene glycol	Kerosene	
The following substances may be used with Noryl at 26 °C or below		
Acetone	Copper sulfate	Mineral oil
Acetic acid	Cyclohexane	Phenol
Arsenic acid	Ethyl acetate	Potassium bichromate
Alcohol	Ethyl chloride	Potassium carbonate
Alkalies	Ferric chloride	Potassium chlorate
Ammonium chloride	Formic acid	Potassium chloride
Ammonium hydroxide	Hydrobromic acid	Potassium hydroxide
Ammonium phosphate	Hydrochloric acid	Potassium nitrate
Ammonium sulfate	Lactic acid	Potassium sulfate
Amyl acetate	Nitric acid	Sodium bicarbonate
Amyl alcohol	Ortho-phosphoric acid	Sodium bisulfate
Borax	Picric acid	Sodium carbonate
Bromine	Phosphoric acid	Sodium chlorate
Butyl alcohol	Stearic acid	Sodium cyanide
Butyl cellosolve	Sulfuric acid	Sodium hydroxide
Butyl mercaptan	Ferrous sulfate	Sodium nitrate
Benzoic acid	Formaldehyde	Sodium perborate
Boric acid	Glycerin, glycerol	Sodium peroxide
Butyric acid	Hydrogen peroxide	Sodium phosphate
Calcium chloride	Iodine	Sodium pyrophosphate
Calcium hydroxide	Isopropanol	Sodium silicate
Carbon bisulfide	Lead acetate	Sodium sulfate
Cesium chloride	Lithium chloride	Sodium sulfite
Chloroacetic acid	Magnesium chloride	Sodium thiosulfate
Chlorosulfonic acid	Magnesium hydroxide	Stannous chloride
Citric acid	Methanol	Sucrose
Copper chloride	Methyl cyclopentane	Tetranitromethane
Copper nitrate		

5.1. General Description

The zonal rotor is employed in the following manner:

Table 6
Separation
Procedure

Item \ Procedure	Loading	Centrifugal Separation	Unloading
Solution flow			
Speed	3,000 rpm	Specified speed	3,000 rpm
Procedure	Inject density gradient, sample, overlay solution	Separation	Inject high density extrusion and recover separated solution

Mount the seal assembly on the rotor being turned at a speed of 3000 rpm, and load a density gradient solution, a sample and overlay solution under the atmospheric pressure with the door of the centrifuge open. Then, remove the seal assembly and place the cap assembly. After closing the door of the centrifuge, start the rotor and raise the rotor speed up to a predetermined speed.

After the centrifugation reduce the rotor speed to 3000 rpm remove the cap assembly, mount the seal assembly again and unload the solution from the rotor by feeding heavy solution into the edge of the rotor. The gradient solution containing sedimented particles is unloaded by feeding a higher-density solution at the edge of the rotor. Also, the absorbancy of the solution thus unloaded is fractionated by a fraction collector while it is being measured by a spectrophotometer. In a common use, prepare the density gradient in the order of high to low density layers from the center to the edge of the rotor. Feed as "cushion" to the gradient solution of higher density a solution of a density equivalent to or higher than the maximum density of the solution. This prevents the heaviest particles from reaching the edge of the rotor when a sample is centrifuged.

After loading the "cushion" stop the pump. Load a sample to be centrifuged beginning with the center of the rotor after load the gradient solution and also load "overlay" or a lower - density solution, so that the sample is layered cylindrically in the rotor. (The sample is loaded at the edge of the rotor as the case may be.) Then, a portion of the cushion previously loaded is discharged in an amount equivalent to the amount of the sample and overlay thus loaded.

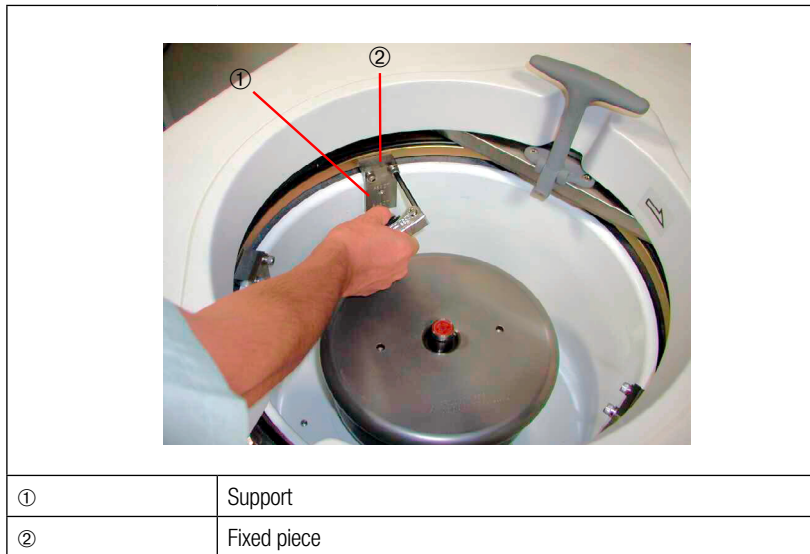
5.2. Preparation

5.2.1. Prepare the Centrifuge

Mount the supports

- a. Mount the three supports onto the fixed pieces in the rotor chamber of the centrifuge as shown in Fig.5-1.
- b. Center and adjust height of guard plate.

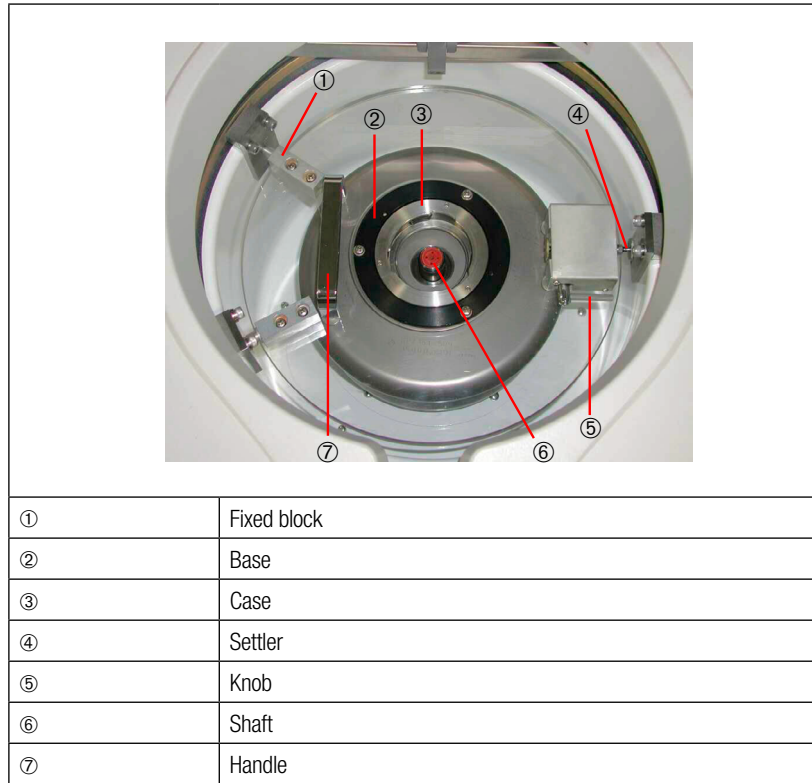
Figure 14
Mounting the
supports



Mounting the guard plate

Put the rotor on the spinning shaft and mount the guard plate. (In TZ-32, mount it to the upper hole in the support.) The guard plate can be fixed by inserting first the end of the fixed block into the left hole on the support, erecting the right knob and then turning it to horizontal.

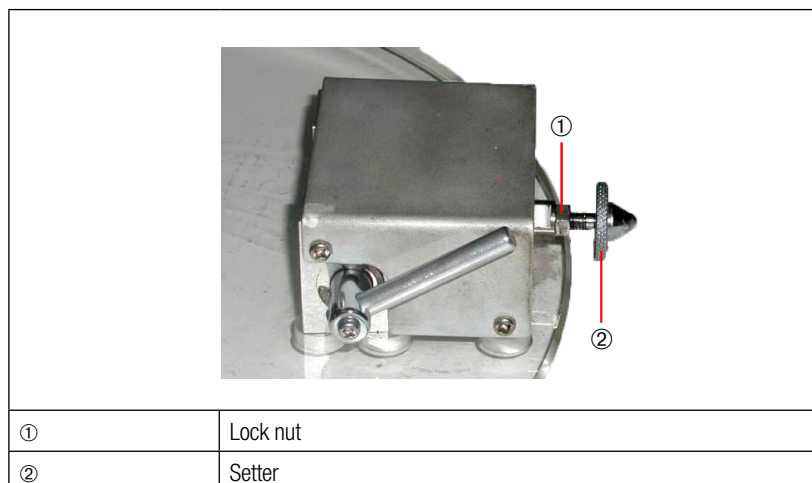
Figure 15
Centering the guard plate



Adjusting the base

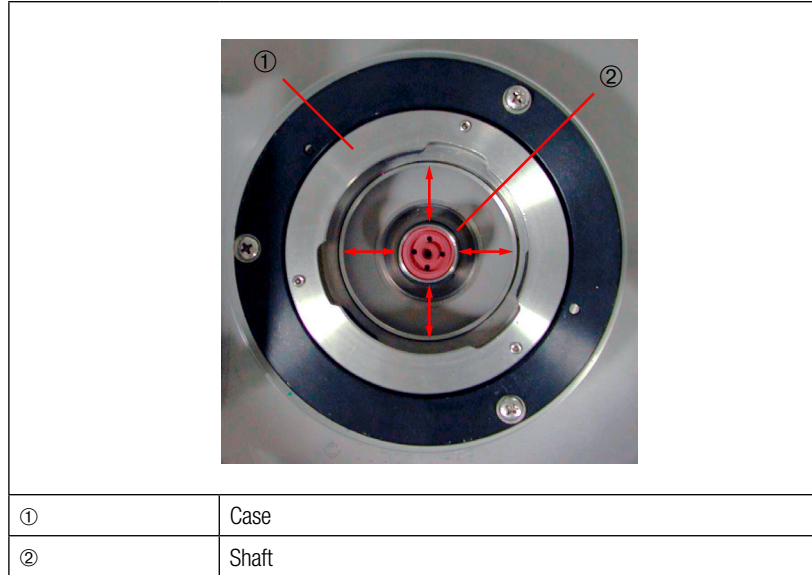
- a. Loosen the four upper screws in the fixed block to make it movable. Then, loosen the lock nut of the setter. Move the fixed block and setter back and forth (by turning the setter) until the acrylic plate is aligned with the rotor center. Tighten the screws to fix them.

Figure 16
Loosening the lock nut



- b. Loosen the four screws in the base, and adjust until the center of the rotor shaft coincides with that of the base. Tighten the screws and fix the base. Adjust the case, while visual checking, until it is concentric with at the center of the shaft.

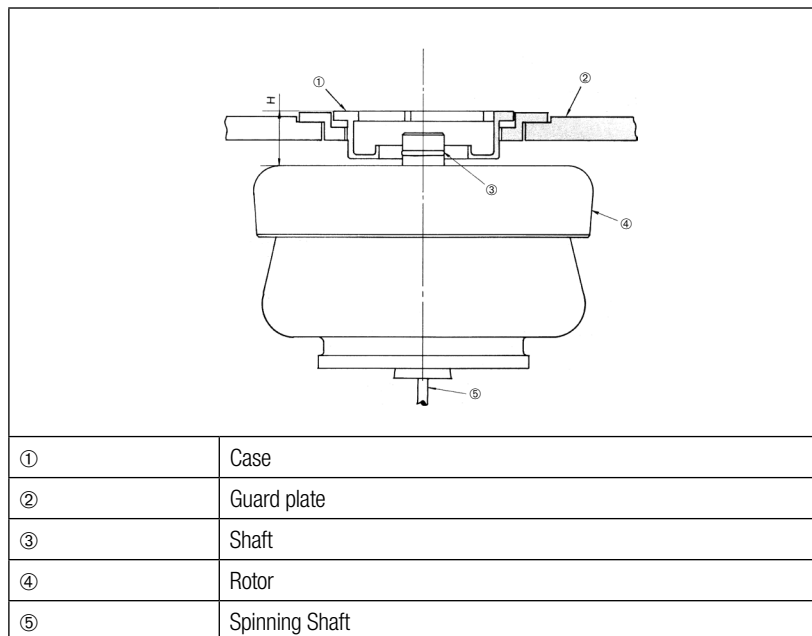
Figure 17
Adjusting the base



Height adjustment of case

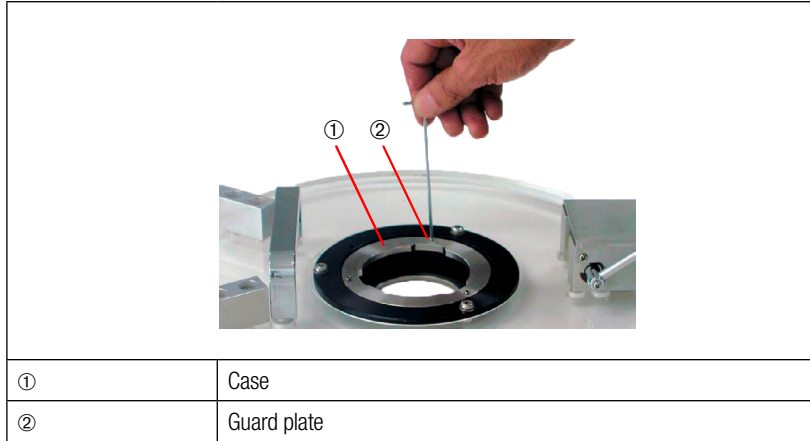
Adjust the height of the rotor and case as shown in [Figure 18](#).

Figure 18
Height adjustment



The height of the case is adjusted as follows; Loosen the three setscrews by the hex. bar wrench accessory to this rotor, and the case can be turned. Turning clockwise lowers it, while turning counterclockwise lifts it. After adjustment, tighten the three setscrews to fix the case.

Figure 19
Tightening the three
setscrews

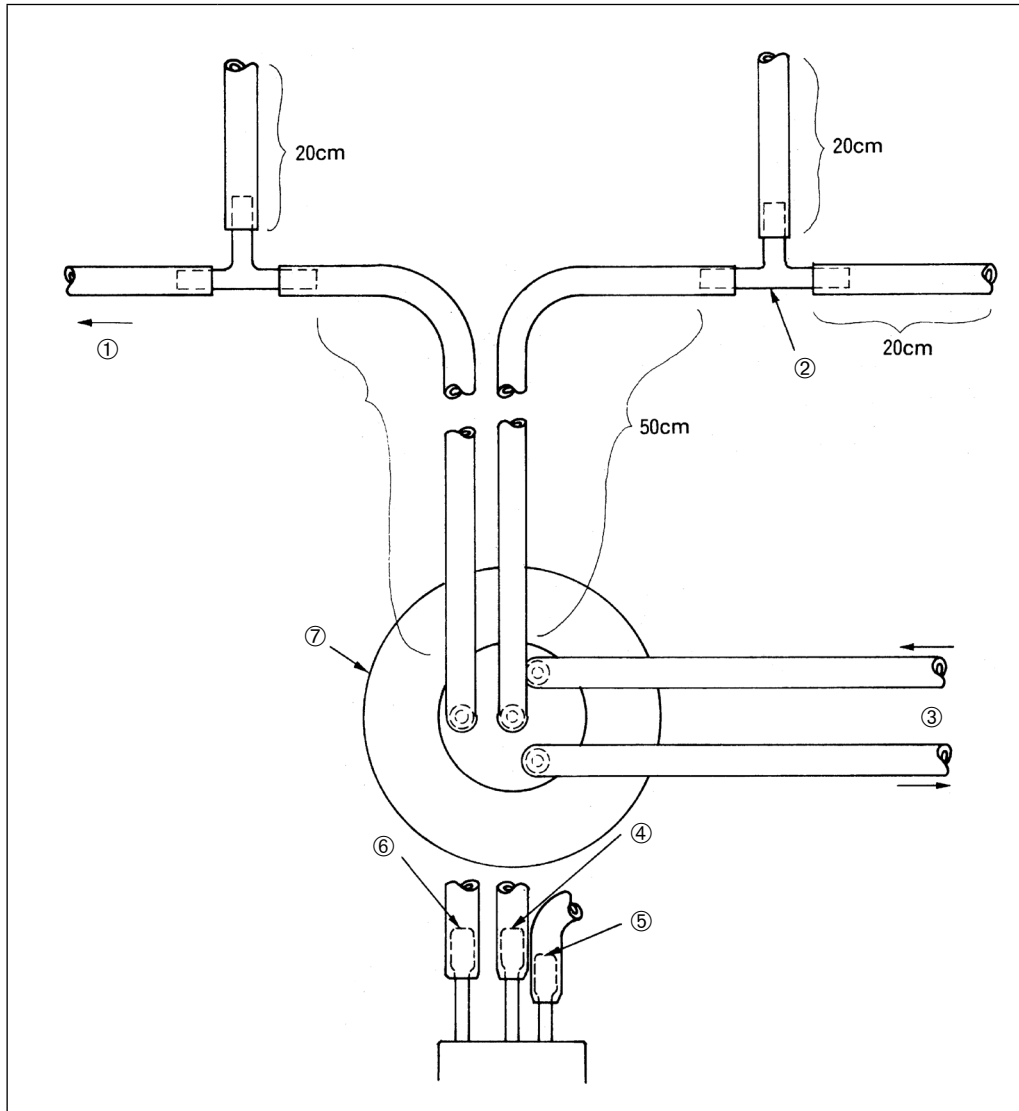


5.2.2. Connect a tube to the seal assembly

Tube connection

For connecting the tubes, see Figure 20. The two higher ones of the four pipes led from the seal assembly are used for injection into or extraction from the rotor. The outer one is led to the outer wall of the rotor and the center one is led to the center of the rotor. The two remaining pipes are used for cooling gradient solution and samples when necessary. Use pipes with an inner diameter of 3.2 millimeter (1/8").

Figure 20
Tube connection



①	Pump
②	Tee
③	Cooling water
④	Pipe on the rotor center
⑤	Cooling pipes (lower at the top end than the other two pipes)
⑥	Pipe on the outer wall of the rotor
⑦	Seal assembly

Cooling water

Cooling water is needed for cooling the contact body, which touches the Rulon seal. Use city water as coolant in case centrifuge is to be done at normal temperature. However use cooled water when centrifuge is to be done at a temperature below 5 °C. The minimum flow rate required is 450 ml/min.

5. 2. 3. Confirm test operation

Check the seal assembly for leakage before injecting a gradient solution into the rotor. This check should be done by injecting distilled water into the rotor by the pump or the injector. The operation and handling are the same as in injecting gradient solution.

Check the seal assembly in the following procedure.

- a. Set the rotor and guard plate, and regulate the rotor speed to 3000 rpm.
- b. Mount the seal assembly with the tubes connected.
- c. Inject about 100 ml of distilled water by the pump or injector through the pipe leading to the outer wall of the rotor.

Leakage from the seal assembly, if any, can be detected from much splashed injected solution sticking to the rotor chamber of the centrifuge. If any leakage is detected, check as follows before performing the preliminary operation again.

Refer to item 7 for the detailed check.

- » Check that the Rulon seal is mounted onto the shaft normally.
(Check that the O-ring is mounted or if there is any scratch on the top surface.)
- » Check that the seal assembly body operates smoothly.
(Check the contact between the Rulon seal and contact body of seal assembly, and also the spring which depresses the guide body.)
- » Check if the outlet pipe on the rotor center blocks normal flow of solution.

For repair of the Rulon seal surface, see „6. 4. Handling of Rulon Seal and Renewal“ on page 44.

5.3. Operation

5.3.1. Prepare the Rotor

- Put the seal packing and metal ring into the rotor. Put the O-ring into the extrusion on the center bottom of the rotor. Apply a thin coat of vacuum grease to the seal packing and O-ring.

CAUTION Set the seal packing and metal ring correct. If they are set incorrect, the seal packing comes out of the seal assembly, thus possibly causing the leakage of solution and damaging the rotor.

CAUTION Since O-ring and seal packing which are used in the rotor or the connecting portion of the seal section will come into contact with sample, always rinse them before use. When these parts are replaced, some residue generated when rubber was formed may remain on the surfaces of new parts. Rinse them away before use.

Figure 21
Setting the seal packing in detail

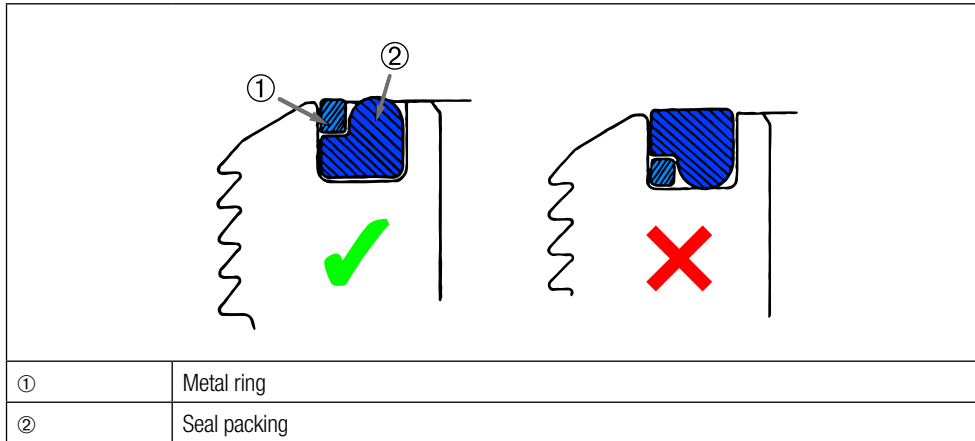
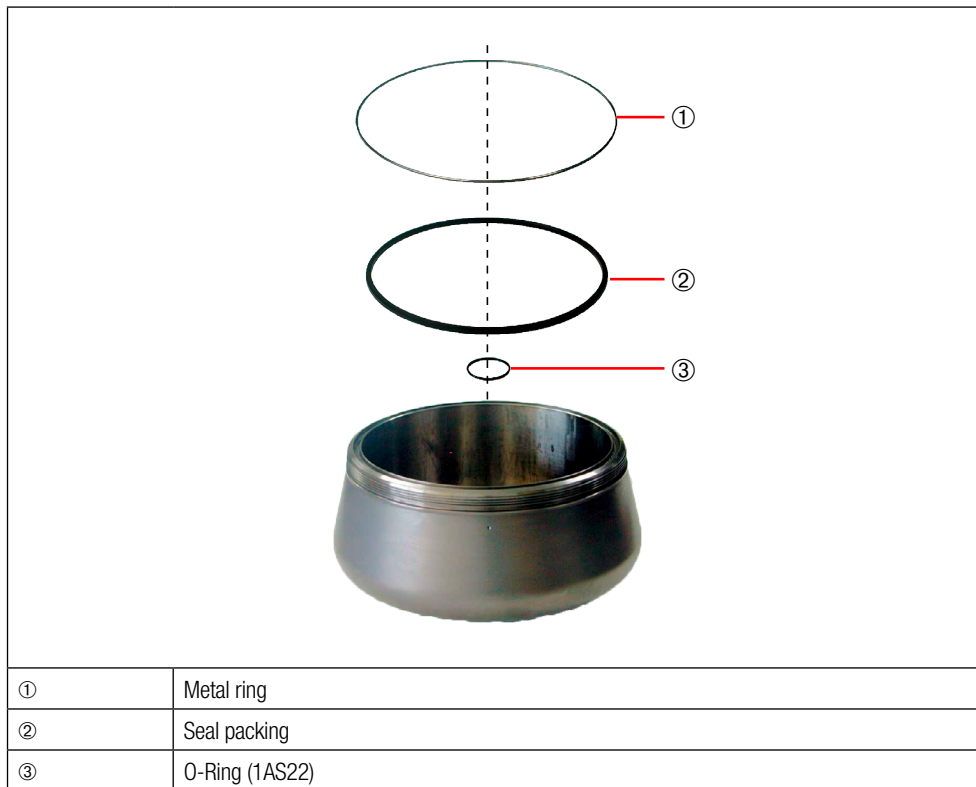


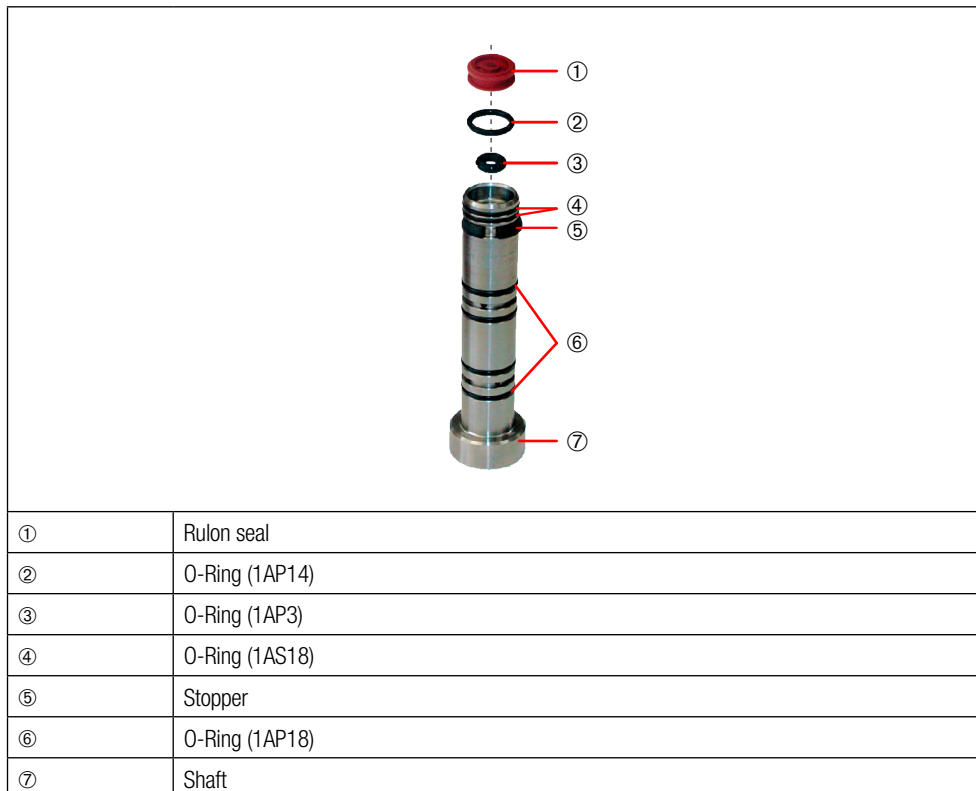
Figure 22
Setting the seal packing



- Put the Rulon seal and O-Ring into the shaft.

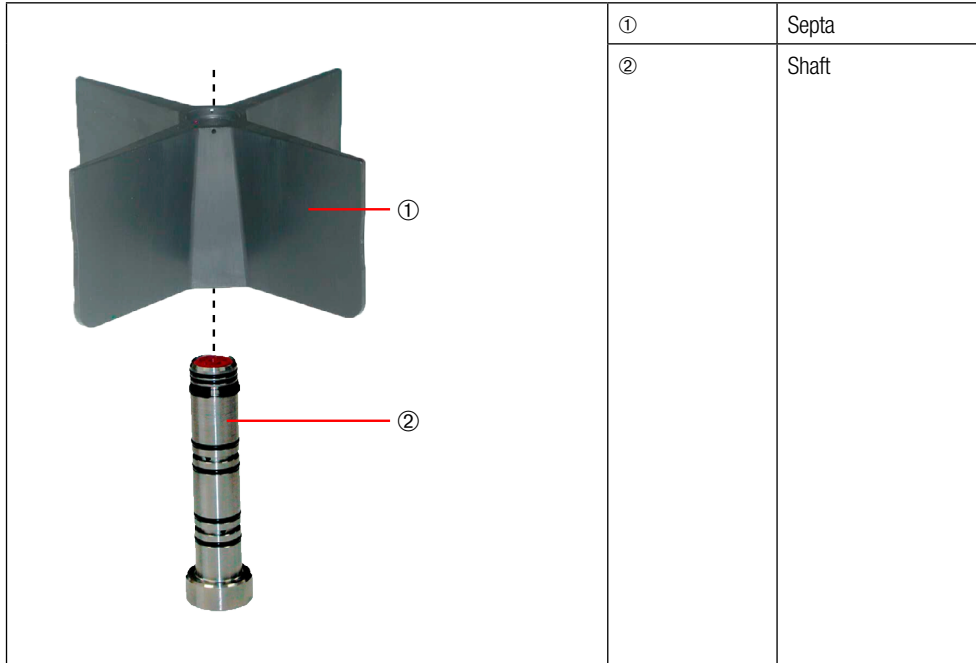
CAUTION Be careful not to scratch the Rulon seal surface by a fingernail. When inserting the seal, carefully push it with the soft portion of finger applied. Apply a thin coat of vacuum grease to the O-ring.

Figure 23
Assembling the Rulon seal and O-ring on the shaft



- Put the shaft into the septa.

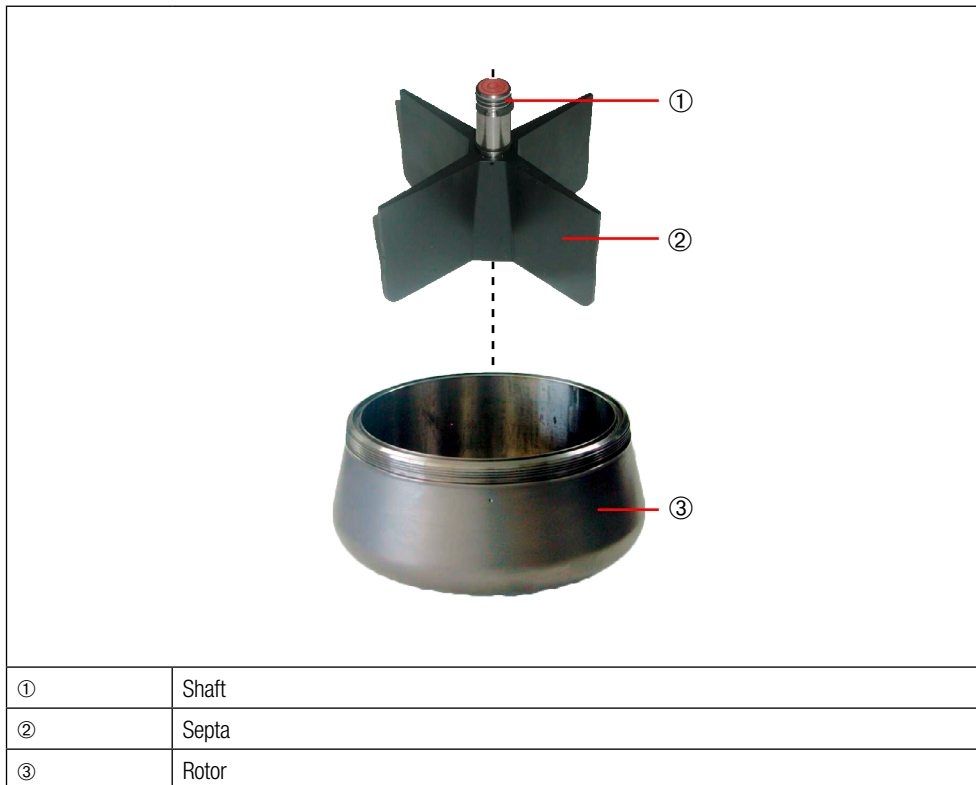
Figure 24
Shaft into septa



- Put the assembled septa into the rotor.

Fit the hole in the bottom end of the shaft onto the extrusion on the rotor bottom where the O-ring (1AS22) is inserted. Thus, the upper surfaces of both rotor and septa will be flush with each other.

Figure 25
Assembled septa into rotor

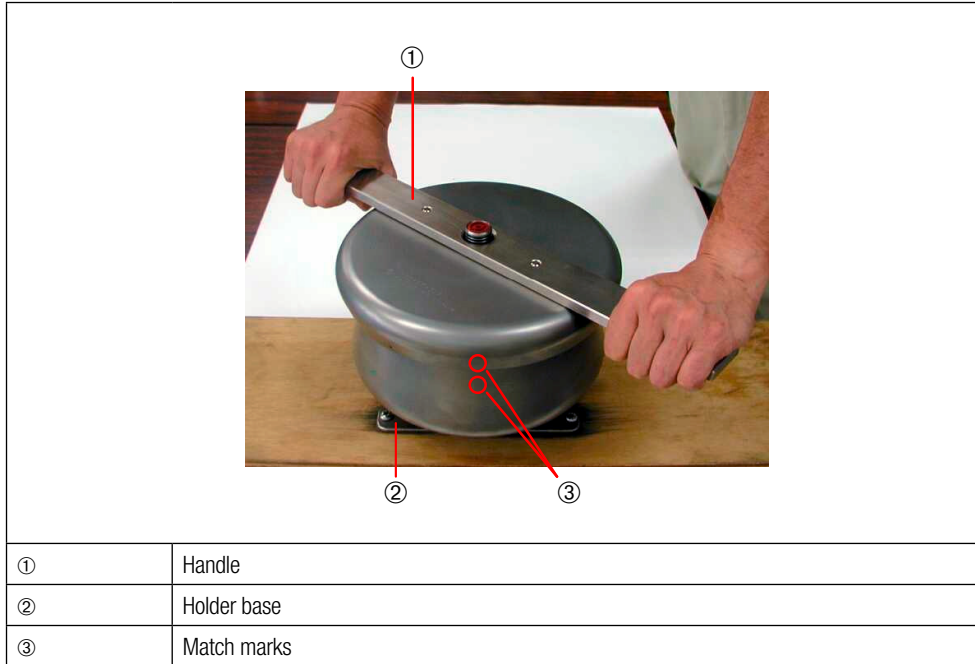


- Put the cover onto the rotor containing the septa.

Put the rotor on the fixed holder base and screw in the cover by hand. Mount the handle as shown in the figure to fasten the cover.

CAUTION Fasten the cover until the match mark on the cover meets that on the rotor body. As the rotor is used over and over again, the match mark on the cover will go over that on, the rotor body, which, however, will result in no trouble.

Figure 26
Assemble the rotor cover



5. 3. 2. Preparation for spinning the rotor

- Mount the rotor on the spinning shaft in the rotor chamber of the centrifuge.
- Mount the guard plate.

Secure the fixed piece on the left support first and fix it by turning over the right-hand support toward the right support.

For setting the supports use the marked upper support hole. Neither mark A nor mark B is employed in the zonal rotor.

Figure 27
Mount the guard plate



5.3.3. Spinning

CAUTION Never raise the rotor speed to more than 5000 rpm with the rotor not filled with the solution to full extent (less than 90%). High speed spinning of the rotor not full of the solution will cause the septa to be damaged. The procedure of spinning the zonal rotor depends on the type of ultracentrifuge used with the rotor. For spinning the rotor carefully read and follow the instructions given below and in the instruction manual of the centrifuge.

Table 7
Operation modes

Speed set to 3000 rpm by acceleration	High speed spinning	Speed set to 3000 rpm by deceleration
1. Set to ZONAL operation mode from panel. 2. Set operating factors for high speed spinning. 3. Press the START switch.	1. Close the door. 2. Press the VACUUM switch. 3. Press the START switch.	1. When the timer counts down to "0", the rotor decelerates to 3000 rpm. 2. Press the VACUUM switch. NOTE When it is desired to stop the rotor, set to the normal operation mode on the panel.

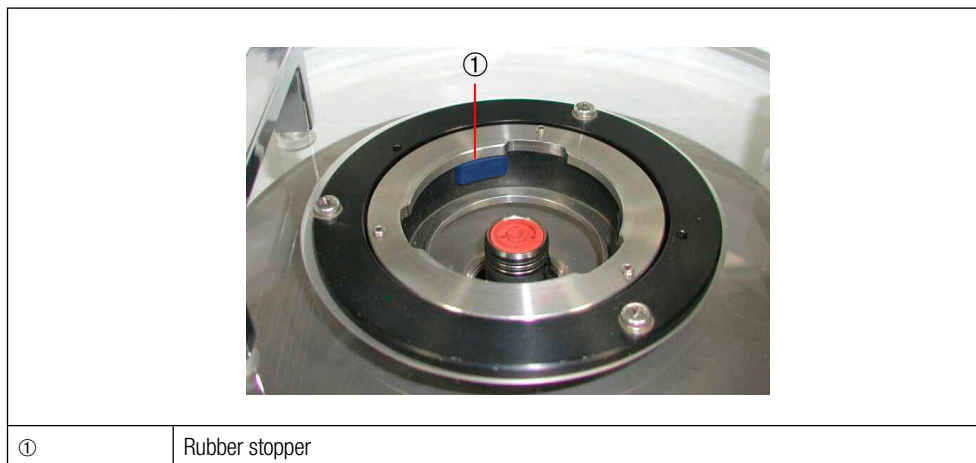
For details on operating from the panel, refer to the instruction manual of the centrifuge.

5.3.4. Injection of gradient solution and sample

1. Mount the seal assembly.

Depress the three extrusions of the seal assembly by fitting them into the three relief grooves. Then, the guide at the end of the seal assembly enters the rotor shaft. Turn the seal assembly clockwise and release it when it touches the rubber stopper.

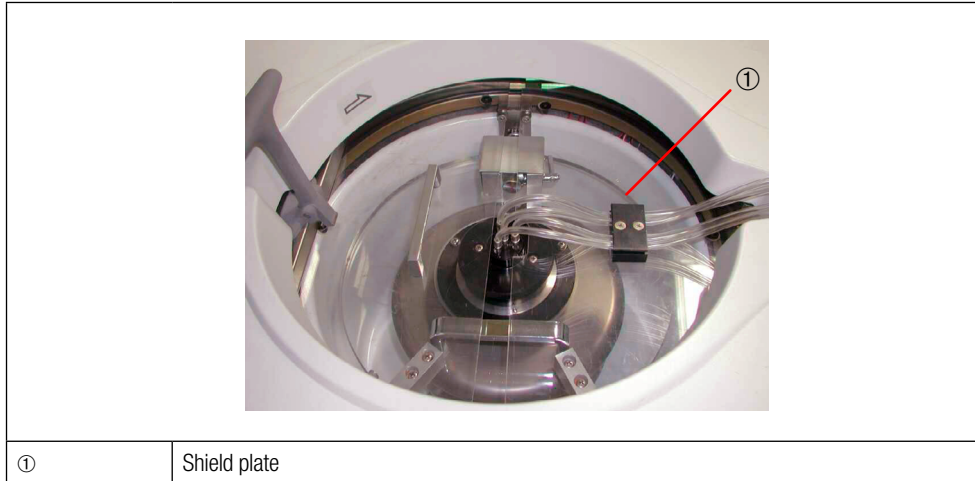
Figure 28
Mount the seal assembly



2. Mount the shield plate.

Put the tube setter side of the shield plate to right or left end of the rotor chamber in the centrifuge. Insert the tubes from the shield plate into the tube setter through the notch of the shield plate. Mount another shield plate, and position it so that there is no clearance at the center.

Figure 29
Mount the shield plate



3. Feed a gradient solution by pump.

a. Pipe on the outer wall of the rotor

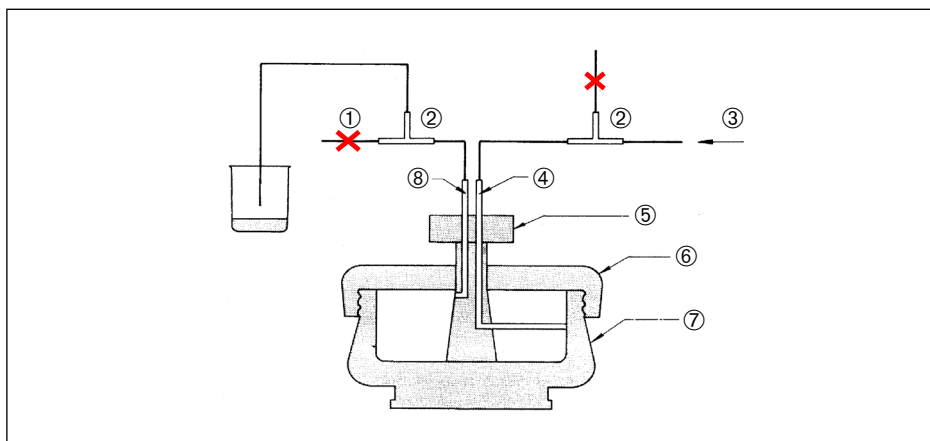
Connect one end to the pump and the other one to the clamp.

b. Pipe on the rotor center

Put one end of the pipe into a beaker containing water. Make sure gradient solution is infected. Clamp the other end.

CAUTION at the time of infecting gradient solution. Select such an infection speed of gradient solution that infection pressure of the pump (pipe pressure at infecting side of the seal is below 1 kg/cm²). Special care should be exercised in selecting an infection speed for a high-viscosity liquid such as sucrose (For 60% solution of sucrose the infection speed is limited to approximately 35 ml/min.)

Figure 30
Feed a gradient solution by pump



4. Injection of sample and overlay

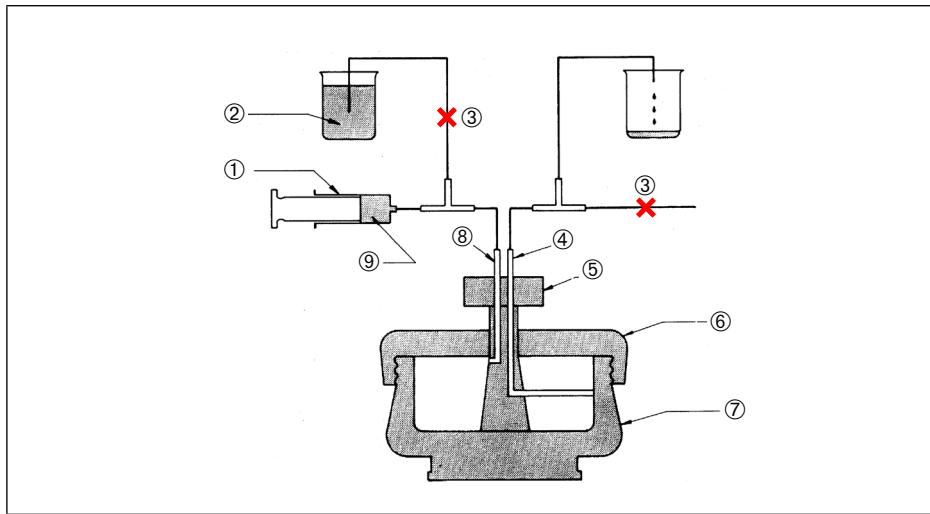
a. Pipe on the outer wall of the rotor

Clamp one end of the pipe at an outer position from the tee on the pump side, and put the other end into a beaker for receiving the "cushion" discharged due to injection of sample and overlay.

b. Pipe on the rotor center

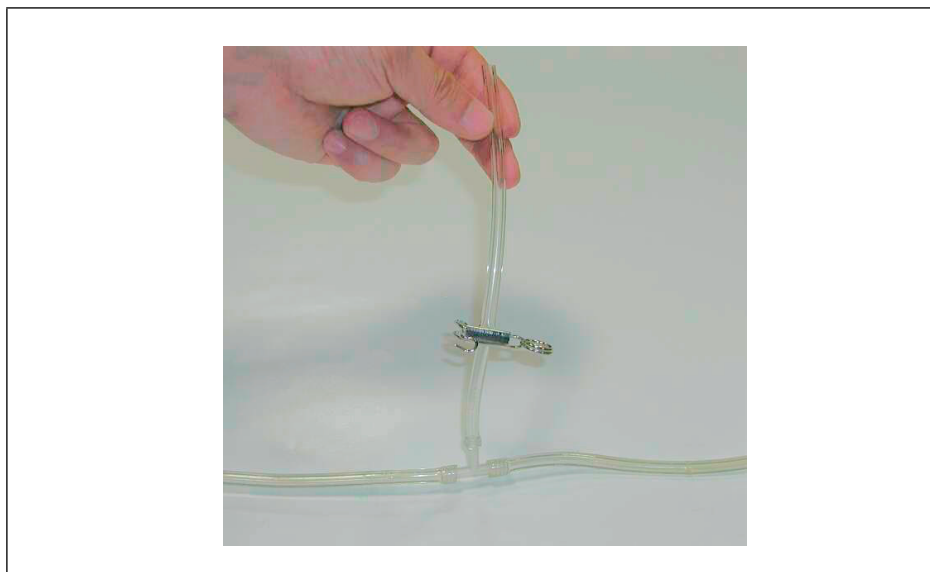
When the lightest gradient solution has filled the tee, clamp both ends of the pipe and stop the pump. Inject a sample through one end by an injector. Then feed an overlay of 75 ml into the rotor.

Figure 31
Injection of sample and overlay



CAUTION Do not feed bubbles while injecting a density gradient solution and sample. Otherwise the gradient solution and sample will be disturbed. If bubbles are about to enter, lead them to another tube from the tee by setting upright the tube containing bubbles before they passthrough.

Figure 32
How to avoid bubbles while injecting a density gradient solution and sample



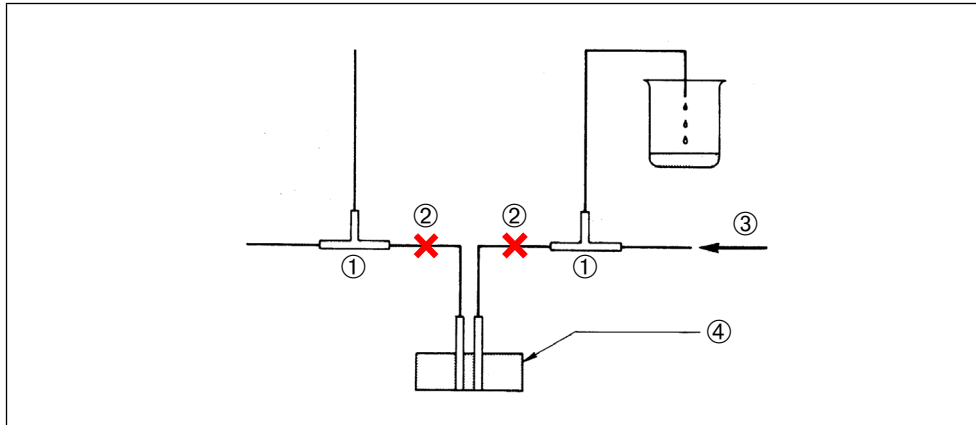
5. Preparation of the pump

Pipe on the outer wall of the rotor and rotor center

Clamp each pipe between the tee and the seal assembly, and remove the seal assembly. Flow a heavy extracting solution through the outerwall of rotor line of the clamped pipe and eliminate bubbles from the pipe on the outer wall of rotorline.

Figure 33

Preparation of the pump



6. Removal of air bubbles

a. Pipe on the outer wall of the rotor

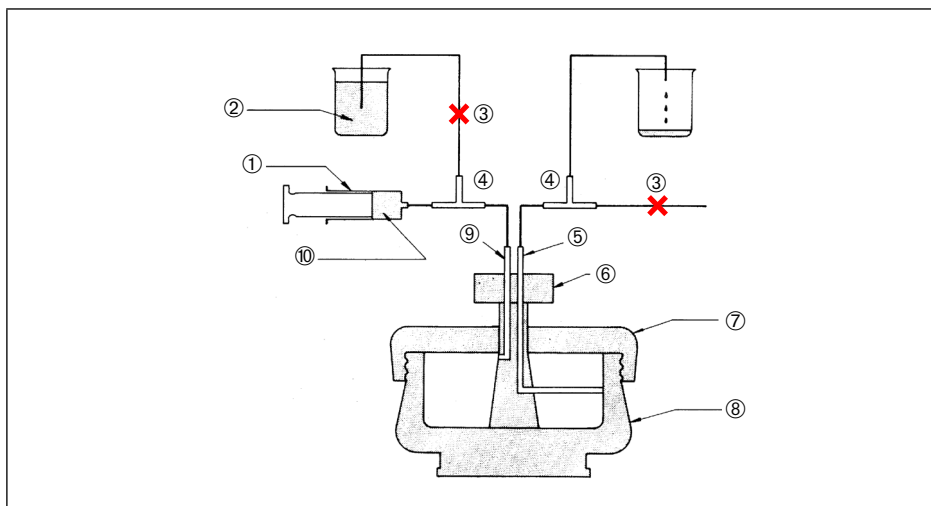
Clamp the pipe on the pump side.

b. Pipe on the rotor center

Prepare about 15 ml of an overlay, disconnect the clamp on the rotor side, and clamp the other end for injection. Bubbles will be removed through the pipe on the rotor center.

Figure 34

Removal of air bubbles



7. Remove the shield plate and remove the seal assembly.

Remove the tubes from the tube setter and then the shield plate. The seal assembly can be pulled ipwards when it touches the stopper while turning it counterclockwise by hand.

5.3.5. High speed spinning

1. Place the cap

Hold the upper fringe as shown in the figure, and push it onto the shaft downwards.

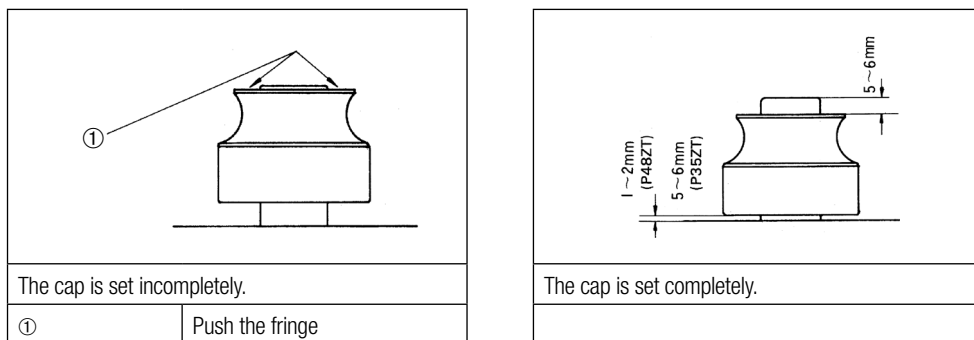
CAUTION Make sure that the cap, once set completely, will not come off the shaft. Even when it is pulled holding the fringe. It is recommended to try setting the cap into the rotor as preliminary practice before spinning the rotor.

Figure 35
Place the cap



2. Referring to the illustrations, check that the cap is set completely.

Figure 36
Checking the cap



3. For the high speed spinning, see „5.3.3. Spinning“ on page 33.

5.3.6. Deceleration and extraction

1. Regulation of the rotor speed to 3000 rpm by deceleration, see „5.3.3. Spinning“ on page 33.

2. Push the R. PUMP button and stop the vacuum pump to let air into the rotor chamber.

3. Remove the cap.

Hold the running cap by both hands. Pull it upwards while pushing the center extrusion.

4. Mount the seal assembly.

See “5.3.4. Injection of gradient solution and sample” on page 33.

5. Unloading the separated solution by pump

- a. Pipe on the outer wall of rotor

Feed by pump with a heavy solution having a density equivalent to or higher than the maximum density of the

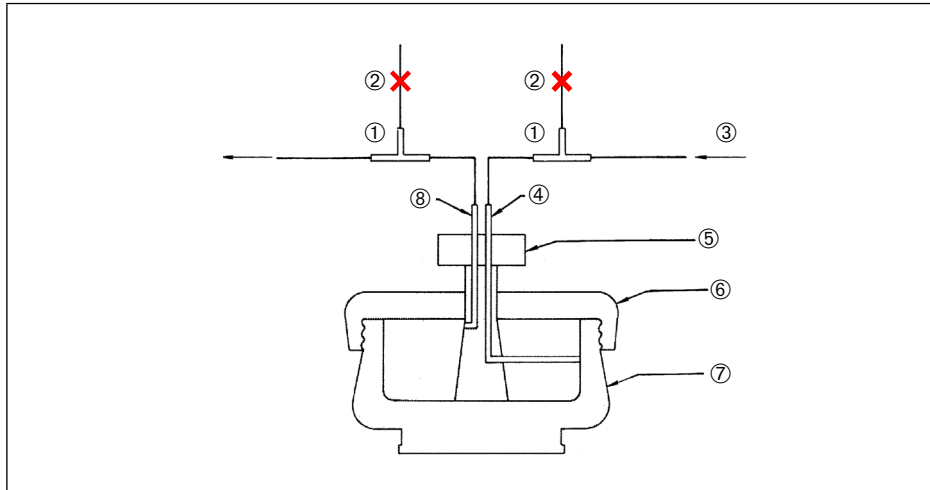
gradient solution.

Clamp other end.

- a. Pipe on the rotor center

Clamp one end. Lead the other end to a fraction collector through a spectrophotometer flow cell.

Figure 37
Deceleration and
extraction



5.4. Operational Checkpoints

5.4.1. Setting of sample temperature and rotor temperature

Make the temperature of injecting sample and density gradient solution and the setting temperature of a rotor when spinning as equal as possible. If not, following problems may occur:

1. When the temperature of injecting solution is lower than the temperature of the spinning rotor:

Because of the temperature rise of the solution injected into the rotor, the volume expands and the pressure inside the rotor increases. The push button of the cap assembly cannot be pushed down and the cap cannot be removed. As the result, it is impossible to unload. Following table shows the quantities of increased volumes when using the TZ-32 rotor and water as sample.

Table 8
Increase of Volume

Temperature Rise	Increase of Volume
5 °C → 15 °C, 10 K	1.5 ml
5 °C → 25 °C, 20 K	5.0 ml

Countermeasure

While operating at 3000 rpm, cool the rotor to the temperature when the solution is injected to the rotor. Then try to remove the cap again. If it is still difficult, stop running the rotor and push down the push button hard, then remove the cap.

2. When the temperature of injecting solution is higher than the temperature of the spinning rotor:

Because of the temperature fall of the solution injected into the rotor the volume decreases and the inside of the rotor becomes a negative pressure status. The cylinder of the cap assembly and the pushbutton are sucked up to the rotor shaft. As the result, the cap cannot be removed. The decreases of volumes are the same values of the increases of volumes shown in the above table.

Countermeasure

Remove the cap while operating at 3000 rpm. If still difficult, rise the temperature of the rotor to the temperature when the solution is injected while operating at 3000 rpm and try to remove the cap again.

Precautions on mounting the seal assembly Wipe the spinning Rulon seal surface with a tissue paper moistened with water lightly just before mounting the seal assembly. This prevents the Rulon seal from wearing abnormally and makes the cap removal easier. Keep the seal assembly as horizontal as possible and center it visually. Lower the seal assembly vertically onto the Rulon seal surface and mount it quickly taking care not to hit the adapter on the Rulon seal surface. Never incline the seal assembly when mounting. If the seal assembly is mounted in inclined condition, the tip of the adapter made of Delrin may hit the Rulon surface and the Rulon may be faulty. The flow resistance of extraction pipe must not be large. The diameter of the extraction pipe must not be small and the length must not be long. The flow cell of the UV monitor must not be a micro flow cell .

5.4.2. Precautions on loading and unloading

1. Do not feed bubbles into the rotor at loading
 - » In the pipe connection, provide bubble eliminating equipment to the position just before the seal
 - » Be sure to clamp the pipe to keep the no-bubble status up to the exit of the contact body of the seal assembly on all occasions and setting is carried out.

- » When the seal assembly is removed due to an injection trouble and the connection is carried out again.
 - a. When the rotor is full

Inject a heavy solution from "OUTER WALL" side to push the bubbles out of the shaft at "CENTER" side.

Open the end of the pipe for eliminating bubbles and inject a heavy solution (normally 20 to 40 ml) with an injector from "OUTER WALL" side until all the bubbles are eliminated from the solution.
 - b. When the rotor is not full

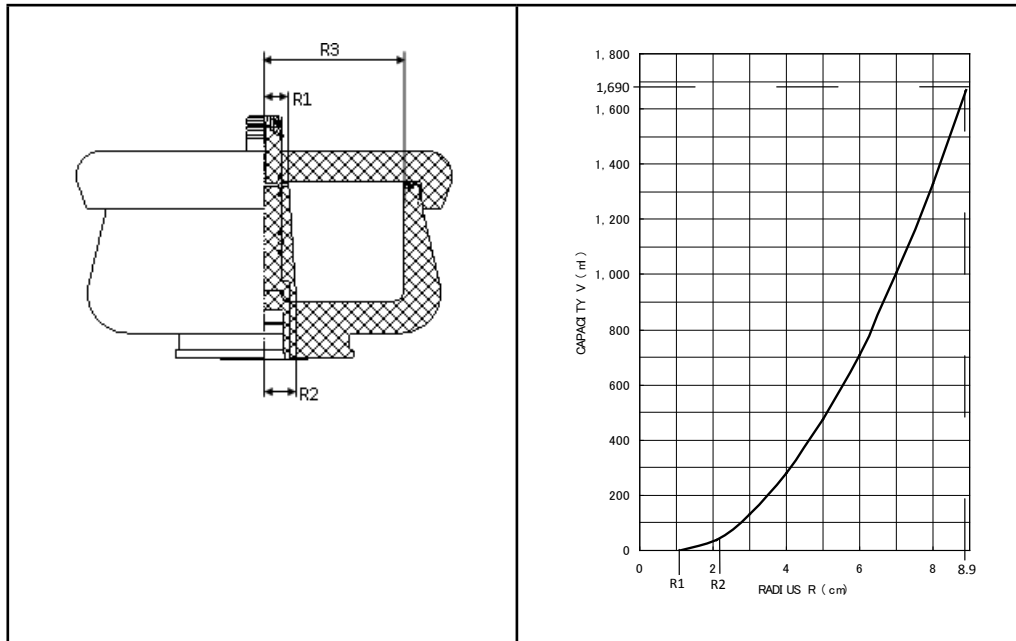
Because there is no way to eliminate bubbles from the injections, continue injection. If not, abnormal injection pressure may be applied to the Rulon seal and cause leakage of solution. In addition when containing bubbles not only the density gradients but also the fractionated patterns may be disturbed in the extraction process.
- 2. Do not feed bubbles into the rotor at unloading

When mounting the seal assembly for unloading after centrifugation, eliminate bubbles at "OUTER WALL" side.

Provide the 'air cushion', that is an air layer in the pipe for eliminating bubbles to absorb the shock of flow against the seal assembly.
- 3. Control the pressure for injecting density gradient solution.
 - » The higher the concentration of solution, the slower the injection speed.
 - » It is permitted to measure the injection pressure with a pressure gauge to keep the injection pressure at 1 kg/cm² or less.
- 4. Precautions on mounting the seal assembly
 - » Wipe the spinning Rulon seal surface with a tissue paper moistened with water lightly just before mounting the seal assembly. This prevents the Rulon seal from wearing abnormally and makes the cap removal easier.
 - » Keep the seal assembly as horizontal as possible and center it visually. Lower the seal assembly vertically onto the Rulon seal surface and mount it quickly taking care not to hit the adapter on the Rulon seal surface. Never incline the seal assembly when mounting. If the seal assembly is mounted in inclined condition, the tip of the adapter made of Delrin may hit the Rulon surface and the Rulon may be faulty.
- 5. The flow resistance of extraction pipe must not be large.
 - » The diameter of the extraction pipe must not be small and the length must not be long.
 - » The flow cell of the UV monitor must not be a micro flow cell

5.5. Capacity in Rotor

Figure 38
Relation between
volume and radius



5.6. Separation Characteristic

Table 9
Separation
characteristic

Speed (rpm)	(rpm) ²	Centrifugal Force (x g)			K Factor
		R = R _{min} (2.05 cm)	R = R _{av} (5.475 cm)	R = R _{max} (8.89 cm)	
10000	1000 x 10 ⁵	2529	6120	9940	3716
15000	2250 x 10 ⁵	5160	13800	22400	1652
20000	4000 x 10 ⁵	9170	24500	39800	929
22500	4840 x 10 ⁵	11100	29600	48100	768
25000	6250 x 10 ⁵	14300	38200	62100	595
27500	7563 x 10 ⁵	17300	46200	75200	491
30000	9000 x 10 ⁵	20600	55000	89500	413
32000	10240 x 10 ⁵	23500	62700	102000	363

5.7. Rotor Life

NOTE Keep the rotor operation log regularly. The rotor operation log is necessary for control of the rotor life.

Table 10
Rotor Life

Operations	Time
1000	2500 h

While using the rotor repeatedly, its strength decreases gradually due to fatigue and creep of material, by which the rotor life is determined. Sum up the number of operations and number of operating hours, and if either value reaches the figures (primary life) shown in Table 7, decrease the maximum speed of the rotor by 10%.

If the operations or operating hours of rotor reach the figures in Table 10 after decreasing the maximum speed by 10% (secondary life), don't use the rotor any more. This specification is applied to the rotor body assembly. The Rulon seal, o-ring, stopper and cap assembly are consumable parts.

6. Maintenance

6.1. Cleaning of Rotor and Septa

After use, disassemble and clean the septa, shaft, metal ring, seal packing, and O-ring.

Cleaning procedure:

Wash it with distilled water at 40 – 50 °C after use, and wipe off water with a soft cloth. After fully drying it, coat it with stopcock grease.

Immediately after a corrosive sample was used, immerse the rotor in neutral detergent (excluding those containing chlorine) diluted with lukewarm water and rinse it sufficiently in distilled water.

Removal of metal ring

Remove it with the accessory screwdriver or the like which has a thin flat end.

Figure 39
Removal of the
metal ring



6.2. Sterilization of Rotor

Sterilize the rotor appropriately following Table 6-1.

Do not autoclave the rotor or sterilize it in boiling water because its material may deteriorate and its strength may decrease.

Table 11
Sterilization
methods

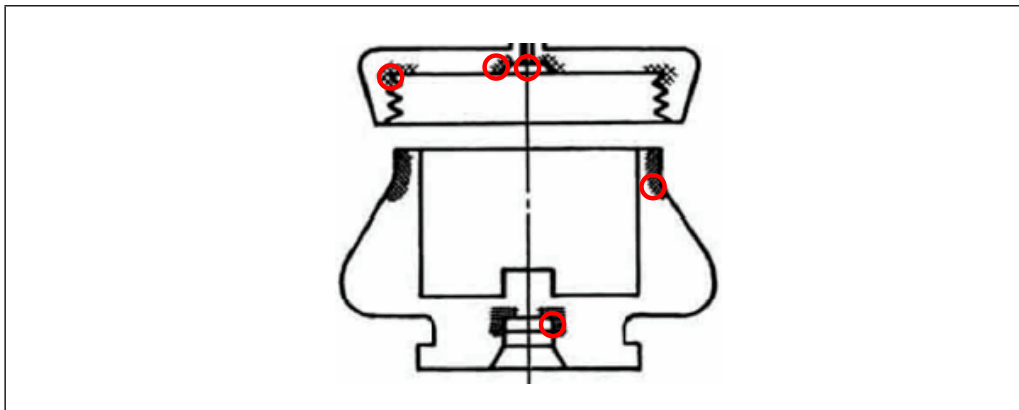
Sterilize Method	Condition	Rotor	
		Usable: ✓	Not usable: ✗
Autoclaving	121 °C (1.0 kg / cm ²) for 20 min		✗
Boiling	15-30 min		✗
Ultraviolet rays	200-300 nm		✓
Gas	Ethylene oxide		✓
	Formaldehyde		✓
Chemical solution	Ethanol (70%)		✓
	Hydrogen peroxide (3%)		✓
	Formalin		✓

6.3. Inspection of Rotors

Check the rotor as follows after every use.

- If the optical adapter/disk is unclean, wipe it with a soft cloth but do not damage it.
- Replacements of packings and O-rings are available.
- If the rotor appears to be corroded, deformation, cracks or discoloration is observed, do not use the rotor and call a Thermo Fisher Scientific authorized sales representative. The rotor may be irreparable depending on the degree of corrosion and wear.

Figure 40
Rotor inspection
points



6.4. Handling of Rulon Seal and Renewal

Rulon seal is a soft plastic with excellent lubrication developed by mixing tetrafluoroethylene resin with a special filler, and is easily scratched by a fingernail. Do not scratch the upper surface, or solution may leak due to poor sealing.

- Renewal of Rulon seal

Slightly scratched or worn Rulon seal can be reconditioned by polishing. The flat portion of the zonal rotor cover can be used for polishing the Rulon seal. Some technique is required for finishing.

Put fine sandpaper (No. 800 is recommendable) on a rotor cover, and polish the contact surface until it becomes even.

Do this carefully while removing polishing agent from the Rulon seal surface and sandpaper. Finish it after making sure no scratch due to the sandpaper remains on the Rulon seal surface.

Figure 41
Polishing the Rulon seal



- Removal of Rulon seal

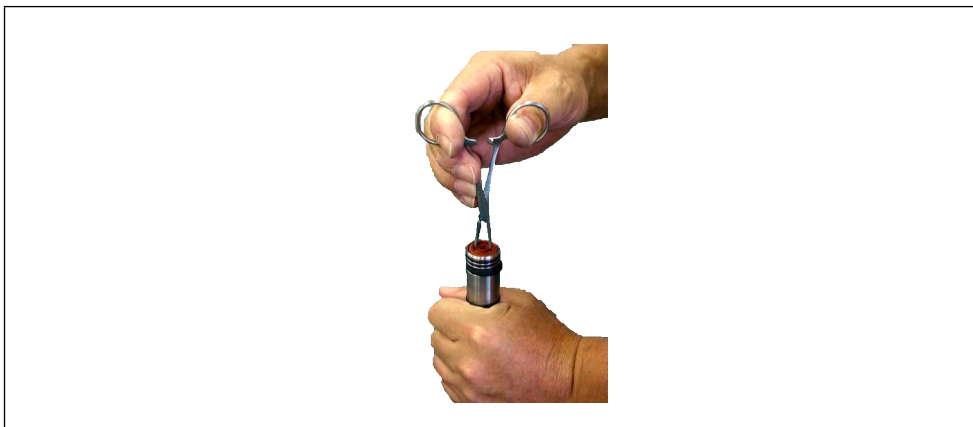
The Rulon seal can be washed and sterilized together with the shaft. Remove the Rulon seal if the Rulon seal is scratched. Be careful not to scratch the Rulon seal surface.

Removing procedure

- Insert the accessory extractor into two holes.
- Turn it clockwise and counterclockwise while lightly depressing it.
- Carefully pull it upwards while turning.

Before resetting it, apply a thin coat of vacuum grease to the O-ring and insert it carefully with soft portions of applied fingers.

Figure 42
Removal of the Rulon seal

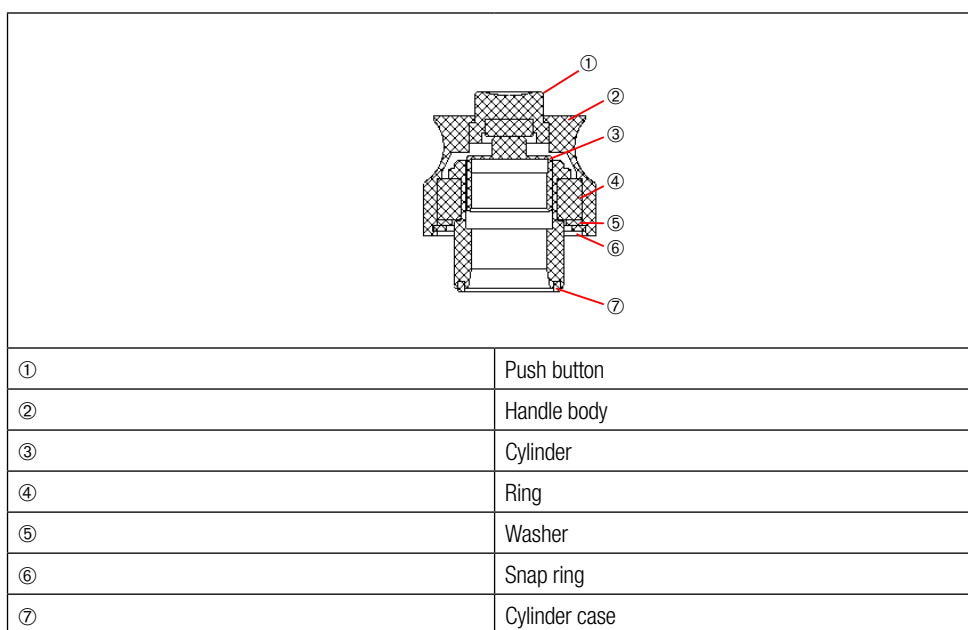


6.5. Disassembling and Cleaning of Cap Assembly

The cap can be disassembled by removing the snap ring using the accessory circlip remover. Be careful not to scratch it since carbon is used as a sliding agent. Wash it in water or clean it by boiling. [Figure 43](#) shows the parts of the cap assembly.

- Check of cap assembly
 - » The cylinder (③) in the cap must move up and down smoothly.
 - » The operation should be almost normal after disassembling and cleaning.
 - » Replace the ring (④) with new one if it has too much play due to wear of ring.
 - » Replace the cap with new one if the corner that catches the stopper is not square.

Figure 43
Parts of the cap assembly



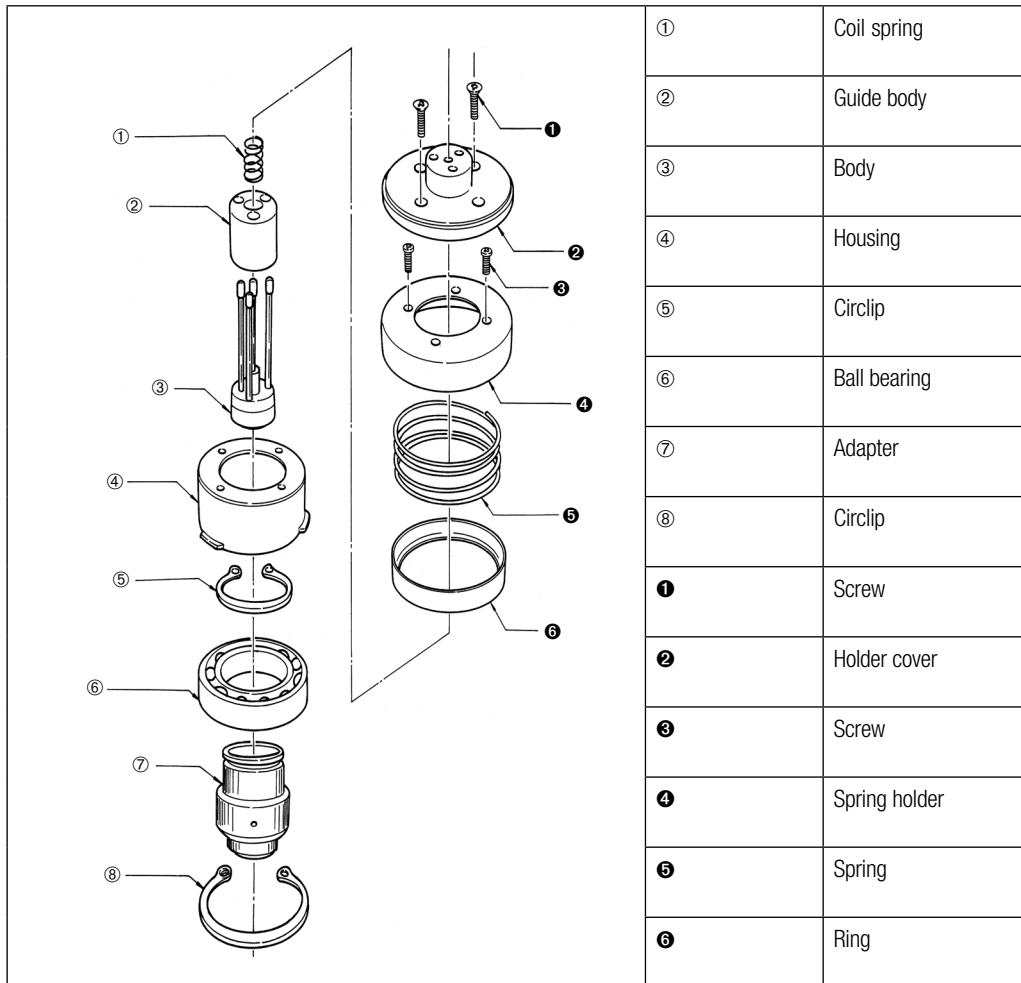
6.6. Disassembling and Cleaning of Seal Assembly

Disassemble the seal assembly by the following procedure. See [“Figure 44”](#) on page 46.

1. Loosen screws (●) and detach the holder cover.
2. Loosen screws (●) and remove the spring holder, spring (B), and ring.
3. Remove the body, guide body, and coil spring (①).
4. Remove the circlip from the housing, and pull out the adapter and ball bearing.

Fully clean the body, housing and ball bearing. Wash the ball bearing in lukewarm water, dry it, and apply the accessory aluminum lubricant to the balls.

Figure 44
Construction of the seal assembly



6.7. Cleaning of Ball Bearing

Don't leave aqueous sucrose or the other in the ball bearing.

Fully wash it in water or lukewarm water use, and dry it sufficiently with an air blower or dryer.

Slightly apply aluminum screw lubricant. After drying, to the balls to the contacts of the balls with between the inner and outer rings.

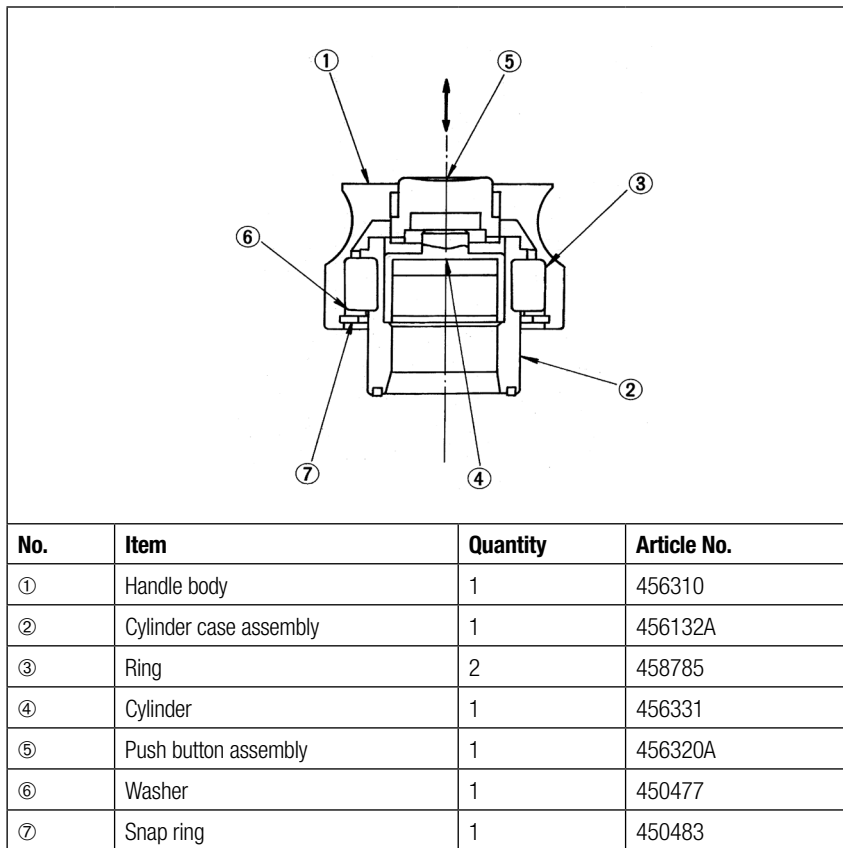
Apply a small amount of lubricant to the inner and outer rings at three positions or so, and turn the bearing so that the lubricant can be applied evenly.

6.8. Consumable Parts of the Rotor

- Rulon Seal
- O-Ring
- Cap assembly
- Seal packing
- Stopper

6.9. Service Parts List for the Cap Assembly

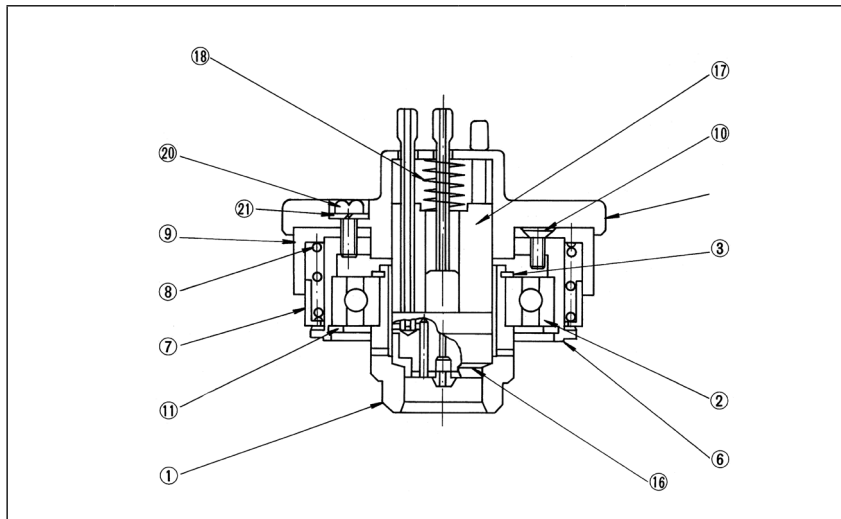
Figure 45
Service parts list for
the cap assembly



6.10. Service Parts List for the Seal Assembly

Figure 46

Service parts list for the seal assembly



No.	Item	Quantity	Article No.
1	Adapter (2)	1	481841
2	BB	1	60000178
3	Circlip	1	481840
6	Housing	1	437842
7	Ring	1	437844
8	Spring	1	437845
9	Spring holder	1	437846
10	M5 x 10 flat head screw	2	80155598
11	Circlip	1	4650213
16	Contact body assembly	1	439825A
17	Guide body	1	437863
18	Coil spring	1	452958
20	M5 x 12 cross-recessed pan head screw	2	80152408
21	M5 spring washer	2	80112058

7. Troubleshooting

Table 12
Troubleshooting

Symptom	Possible Cause	Corrective Action
Pressure is 1.0 kgf/cm ² or higher when density gradient solution is injected.	<ol style="list-style-type: none"> 1. The injection flow rate is too high. 2. The injection passage is blocked. 	<ol style="list-style-type: none"> 1. The appropriate flow rate is approx. 50 ml/min. 2. Disassemble and check. Blocking is likely to occur in the seal assembly, shaft and septa.
Bubbles enter the sample line.	<ol style="list-style-type: none"> 1. The pressure rises. 2. The solution stops flowing. 	<ol style="list-style-type: none"> 1. If the pressure does not rise much, it is possible to continue operation. 2. If the pressure is too high or the sample stops flowing, remove bubbles. See "5. 4. 2. Precautions on loading and unloading" on page 39.
Pressure drops abruptly during loading and unloading.	The sample injection line is bypassed.	See "Table 6" on page 39.
Sample leakage and vibration of seal assembly.	<ol style="list-style-type: none"> 1. The installation condition of the guard plate. 2. The condition of the seal assembly. 	<ol style="list-style-type: none"> 1. Level the guard plate. Rattle the three supports. Center the rotor and guard plate. See "5. 2. Preparation" on page 23. Adjust the height of rotor and case. 2. Check the surface of the contact body for the Rulon seal for scratches and wear. Check the bonded portion of the rear of the contact piece for peeling. Check the adapter inside for scratches and wear. Spins the bail bearing smooth? Clean and grease the bail bearing.
Temperature of the rotor rises.	Poor degree of vacuum.	Precool the rotor body and sample to separate at low temperatures. See "5. 4. 1. Setting of sample temperature and rotor temperature" on page 39.

Table 13
Defective
phenomena and
corrective action
sealed sections

Symptom	Location		Probable Cause	Corrective Action
	Part Name	Portion		
Bypassing and leak of sample and solution.	Seal assembly (Contact body)	WC sealing face (in contact with seal)	1. Poor sealing due to surface roughness or injury.	Replace the contact body assembly.
			2. Poor follow-up in vertical direction of contact body assembly with seal.	Check the spring which depresses the guide body.
Bypassing and leak of sample and solution.	Rulon seal	Contact face	Poor sealing due to surface roughness, injury or step between inner and outer rings.	Replace the Rulon seal.
Bypassing		Insertion recess for O-ring	Poor sealing due to injury in recess portion in contact with O-ring.	Replace the Rulon seal.
		O-ring (1AP3)	Poor sealing due to degradation or injury.	Replace the O-ring.
Outer wall line solution line leaks		O-ring (1AP14)	1. Poor sealing due to degradation or injury.	
			2. Poor sealing due to insufficient grease.	Apply grease.

7.1. Check the Rulon Seal

There are most problem occurrences of Rulon seal in the defect in the operation of zonal rotor.

Explain the point of the check of Rulon seal.

- Check whether the Rulon seal is faulty or not
 - » Scratches or dents on surface
 - » Abnormal wear on surface due to previous sample leakage

Is the height of the surface of the outside circle and the surface of the inside circle the same?

CAUTION Be extremely careful when removing the Rulon seal because the removal of the Rulon seal from the rotor may influence the damage (deformation) to the Rulon seal greatly.

When mounting a new Rulon seal, check it according to the above procedure.

- Check of Rulon seal for stable placement
 - The O-rings must be mounted to the perimeter and the bottom of the Rulon seal correctly.
 - When mounting the Rulon seal to the rotor shaft, press the Rulon seal comparatively hard so that it is mounted deeply without inclination.
 - Apply a thin coat of vacuum grease to both the big and small O-rings.

7.2. Other Check Points

- Check of O-rings of rotor shaft for wear
Remove the O-rings and check them for wear or crack by stretching a little. If worn or cracked, replace the O-rings with new ones.
- Wear or elongation of stopper
Check the lower corners of the stopper for wear to prevent the cap from falling off.
Replace the stopper with new one every six to twelve months.
- Check of other O-rings mounted to the shaft for scratch
Be careful when inserting the shaft into the septa because the O-rings are susceptible to scratches and cuts. Before inserting, apply a thin coat of vacuum grease to the shaft and O-ring for smooth insertion.
- Check of sample passage of septa for plugging
Insufficient maintenance after use may result in a partial plugging.

Thermo Electron LED GmbH
Zweigniederlassung Osterode
Am Kalkberg, 37520 Osterode
Germany

thermoscientific.com/centrifuge

© 2015 Thermo Fisher Scientific Inc. All rights reserved.

Delrin, TEFLON, and Viton are registered trademarks of DuPont. Noryl is a registered trademark of SABIC. POLYCLEAR is a registered trademark of Hongye CO., Ltd. Hypaque is a registered trademark of Amersham Health As. RULON A and Tygon are registered trademarks of Saint-Gobain Performance Plastics. Alconox is a registered trademark of Alconox. Ficoll is a registered trademark of GE Healthcare. Haemo-Sol is a registered trademark of Haemo-Sol. Triton X-100 is a registered trademark of Sigma-Aldrich Co. LLC. Valox is a registered trademark of General Electric Co.

All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.
Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

United States/Canada +1 866 984 3766
Latin America +1 866 984 3766
Austria +43 1 801 40 0
Belgium +32 53 73 42 41
France +33 2 2803 2180
Germany national toll free 0800 1 536 376
Germany international +49 61 84 90 6000
Italy +39 02 95059 552

Netherlands +31 76 579 55 55
Nordic/Baltic States/CIS +358 9 329 10200
Russia +7 812 703 42 15
Spain/Portugal +34 93 223 09 18
Switzerland +41 44 454 12 22
UK/Ireland +44 870 609 9203
India +91 22 6716 2200

China +800 810 5118 or
+400 650 5118
Japan +81 3 5826 1616
Other Asian Countries +852 2885 4613
Australia +61 39757 4300
New Zealand +64 9 980 6700
Countries not listed +49 6184 90 6000 or
+33 2 2803 2180

Thermo
SCIENTIFIC
Part of Thermo Fisher Scientific