



## VTi 50.1 Vertical Rotor

### Straight Up Purification

Purpose-built for Optima Series ultracentrifuges, the VTi 50.1 rotor helps you produce more in less time with rapid density gradient separations. Its short pathlength enables faster run times at high resolution, while the durable rotor structure enables higher density applications without de-rating speed.

#### Faster Density Gradient Formation

With vertical rotors like the VTi 50.1, run times are considerably shorter, which allows you to produce more sellable product faster.



#### Common Consumables and Adapters

With the same consumables and adapters as other Beckman Coulter Life Sciences rotors, it fits seamlessly into your rotor library and gives you added workflow versatility.



#### 50% More Volume

With a 468 mL volume capacity (12 tubes at 39 mL each), it's the highest-volume vertical rotor on the market. That means you can purify more samples in less time with each run, ultimately increasing your throughput.

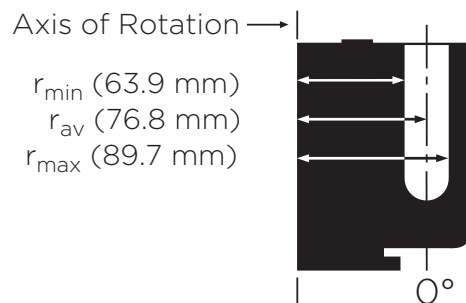


#### Robust Construction

Made of titanium and manufactured in an ISO 9001 facility with a 5-year warranty, this rotor gives you the peace of mind you deserve.



Specification	Value	
Maximum Speed	50,000 RPM	
Density rating at maximum speed	1.7 g/mL	
Relative centrifugal field (RCF) at maximum speed	at $r_{max}$ (89.7 mm)	251,000 x $g$
	at $r_{av}$ (76.8 mm)	215,000 x $g$
	at $r_{min}$ (63.9 mm)	179,000 x $g$
$k$ factor of rotor's cell hole at maximum speed ( $r_{max}$ and $r_{min}$ )	34	
Number of bottle cavities	12	
Nominal tube dimensions (largest tube)	25 x 89 mm	
Nominal bottle capacity	39 mL	
Nominal rotor capacity	468 mL	
Weight of fully-loaded rotor	12.6 kg (27.8 lbs)	
Approximate acceleration time to maximum speed (fully loaded)	13 minutes	
Approximate deceleration time from maximum speed (fully loaded)	14 minutes	
Rotor body material	Titanium	



### Ordering Information

VTi 50.1 Rotor Package	C78759
VTi 50.1 Rotor Package + Software USB*	C82834

\* The USB contains a software update for your instrument to identify the VTi 50.1. Coordinate with your local service team to install the software update.

**VTi 50.1 is compatible with all Beckman Ultracentrifuges classified R and S. Order info below:**

Model	Part Numbers*			
	Instrument		Biosafe Package** <b>BIOE</b>	
	IVD <sup>1</sup>	Non-IVD <sup>2</sup>	IVD <sup>1</sup>	Non-IVD <sup>2</sup>
Optima XPN-100	A99846	A94469	B10053	B10048
Optima XPN-90	A99842	A94468	B10052	B10047
Optima XPN-80	A99839	A95765	B10051	B10046
Optima XE- 100	A99836	A94516	B10050	B10045
Optima XE-90	A99833	A94471	B10049	B10044

\* Availability varies by geography. Contact your local sales representative to determine the correct part number for ordering.

\*\* BioSafe Package includes centrifuge and HEPA filter kit.

<sup>1</sup> In Vitro Diagnostics Use.







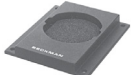




<sup>2</sup> For research use only. Not for use in diagnostic procedures.

## COMPATIBLE LABWARE

Tube			Required Accessory			VTI 50.1
Dimensions/ Nominal Volume	Description	Part Number	Description	Part Number	Tube Rack	Max. Speed/ RCF/ k factor
25 X 89 mm 39 mL	Certified Free & Sterile Ultra-Clear Quick-Seal	C14299 Carton of 48 (6 packs of 8)	Aluminum spacer	342417	348124	50,000 RPM 251,000 X g 34
	Certified Free Ultra-Clear Quick-Seal	C14283 (pkg/50)				
	Standard Ultra-Clear Quick-Seal	344326 (pkg/50)				
25 X 89 mm 39 mL	Certified Free & Sterile Polypropylene Quick-Seal	C14304 Carton of 48 (6 packs of 8)	Aluminum spacer	342417	348124	50,000 RPM 251,000 X g 34
	Certified Free Polypropylene Quick-Seal	C14288 (pkg/50)				
	Standard Polypropylene Quick-Seal	342414 (pkg/50)				
25 X 89 mm 36.2 mL	Polypropylene OptiSeal	362183* (pkg/56)	Gold-anodized aluminum spacer	362204	360542	50,000 RPM 251,000 X g 34
25 X 64 mm 27 mL	Ultra-Clear Quick-Seal	344323 (pkg/50)	Aluminum spacer	342417	348124	50,000 RPM 251,000 X g 34
			Floating spacer	343448		
25 X 64 mm 27 mL	Quick-Seal Polypropylene	343665 (pkg/50)	Aluminum spacer	342417	348124	50,000 RPM 251,000 X g 34
			Floating spacer	343448		
25 X 35 mm 15 mL	Ultra-Clear Quick-Seal	344324 (pkg/50)	Aluminum spacer	342417	348124	50,000 RPM 251,000 X g 34
			Floating spacer	343448 (2 required)		
25 X 35 mm 15 mL	Quick-Seal Polypropylene	343664 (pkg/50)	Aluminum spacer	342417	348124	50,000 RPM 251,000 X g 34
			Floating spacer	343448 (2 required)		

\* Includes disposable plastic plugs

## ACCESSORIES

	Accessory	PN
<b>Included with Purchase</b>		
	12 Rotor Plugs (Red)	355587
	12 Gaskets for Rotor Plug	340825
	2 oz Spinkote Lubricant	306812
	2 Overspeed Discs	330336
	Torque wrench	369791
	Hex plug adapter	355588
	Rotor Vise	332688
<b>Quick-Seal Tube Topper + Accessories</b>		
	Quick-Seal Cordless Tube Topper Kit, 60 Hz (US, Canada, & Japan)	358312
	Quick-Seal Cordless Tube Topper Kit, 50 Hz (Europe)	358313
	Quick-Seal Cordless Tube Topper Kit, 50 Hz (Great Britain)	358314
	Quick-Seal Cordless Tube Topper Kit, 50 Hz (Australia)	358315
	Quick-Seal Tube Rack for 25 mm diameter tubes	348124
	Hemostat (to help remove quickseal tubes)	927208
<b>OptiSeal Tube Accessories</b>		
	OptiSeal Tube Rack for 25 mm diameter tubes	360542
	Tube Extraction Tool	361668