



The tools needed to purify viral vectors with density gradient ultracentrifugation



INSTRUMENT

Model	Part Numbers*			
	Instrument		Biosafe Package** BIOE	
	IVD ¹	Non-IVD ²	IVD ¹	Non-IVD ²
Optima XPN-100	A99846	A94469	B10053	B10048
Optima XPN-90	A99842	A94468	B10052	B10047
Optima XPN-80	A99839	A95765	B10051	B10046

* 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.

¹ In Vitro Diagnostics Use.

² For research use only. Not for use in diagnostic procedures.



XPNs can be programmed to have changes in speeds during the sample centrifuge run to optimize purity + time, which can help your density gradient achieve equilibrium faster

ROTORS

Greatest efficiency for density-based separations

Vertical	
Small scale	Large scale
VTi 90	VTi 50.1



VTi 90
362751



VTi 50.1
C78759

Versatility for density- and sized-based separations

Vertical		
Small scale	Large scale	
Type 100 Ti	Type 70 Ti	Type 50.2 Ti



Type 100 Ti
363013



Type 70 Ti
337922



Type 50.2 Ti
337901

- Small scale rotors are ideal for designing experiments with minimal sample
- Large scale rotors are ideal for production and manufacturing
- While swinging bucket rotors can purify viral vectors, the longer run time, lower volume per run, and comparable resolution make it a less attractive rotor for purifying viral vectors at scale

COMMONLY USED TUBES

SMALL SCALE			
	mL	Compatible Rotors	Quick-Seal
			Non-sterile
Polypropylene	2	Type 100 Ti + VTi 90	345829
	3.5	Type 100 Ti + VTi 90	349621
	5.1	Type 100 Ti	362248
	6	Type 100 Ti	344619
Ultra-Clear	5.1	VTi 90	344075

LARGE SCALE					
	mL	Compatible Rotors	OptiSeal	Quick-Seal	
				Non-sterile	Sterile
Polypropylene	15	VTi 50.1, Type 50.2 Ti, Type 70 Ti		343664	
	27	VTi 50.1, Type 50.2 Ti, Type 70 Ti		343665	
	32.4	Type 50.2 Ti, Type 70 Ti	361625		
	36.2	VTi 50.1	362183		
	39	VTi 50.1, Type 50.2 Ti, Type 70 Ti		342414	C14304
Ultra-Clear	15	VTi 50.1, Type 50.2 Ti, Type 70 Ti		344324	
	27	VTi 50.1, Type 50.2 Ti, Type 70 Ti		344323	
	39	VTi 50.1, Type 50.2 Ti, Type 70 Ti		344326	C14299

For a complete list of tubes available for each rotor type, please see the Rotor IFUs or Ultracentrifuge Catalog



© 2022 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the Stylized Logo, and Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries. All other trademarks are the property of their respective owners.

For Beckman Coulter's worldwide office locations and phone numbers, please visit Contact Us at beckman.com
22.03.4533.CENT