

Rotor/Adapter Selection Guide

Centrifuge	Rotor Name	Rotor Type*	Adapter for BD Falcon™ 15 ml Conical Tube	Adapter for BD Falcon™ 50 ml Conical Tube
Beckman Coulter, Inc.				
GS-15/R	S4180	SW	361230	361234
	C0650	FA	—	Dedicated 50 ml rotor
	C1015	FA	Dedicated 15 ml rotor	—
TJ-6/R	TH-4	SW	339276 359487, aerosol, set of 2 358991, aerosol, set of 4	339273 359488, aerosol, set of 2 358992, aerosol, set of 4
GS-6/R/K/KR	GH-3.8	SW	359151, set of 4 359472, set of 2 359487, aerosol, set of 2 358991, aerosol, set of 4	359154, set of 4 359475, set of 2 359488, aerosol, set of 2 358992, aerosol, set of 4
Avanti 30	C0650	FA	—	Dedicated 50 ml rotor
	C1015	FA	Dedicated 15 ml rotor	—
Allegra X-12	SX4750	SW	359472, standard bucket, set of 2 359151, standard bucket, set of 4	393266, standard bucket, set of 2 393267, standard bucket, set of 4
Allegra X-15/R	SX4750A	SW	359487, Aerosolve®, set of 2 358991, Aerosolve, set of 4	359488, Aerosolve, set of 2 358992, Aerosolve, set of 4
		SW	359472, standard bucket, set of 2 359151, standard bucket, set of 4	393266, standard bucket, set of 2 393267, standard bucket, set of 4
		SW	359487, Aerosolve, set of 2 358991, Aerosolve, set of 4	359488, Aerosolve, set of 2 358992, Aerosolve, set of 4
	FX6100	FA	392270	392268
Allegra 6/R Spinchron DLX	GH-3.8	SW	359472, standard bucket, set of 2 359151, standard bucket, set of 4	393266, standard bucket, set of 2 393267, standard bucket, set of 4
		SW	359487, Aerosolve®, set of 2 358991, Aerosolve, set of 4	359488, Aerosolve, set of 2 358992, Aerosolve, set of 4
		SW	359472, standard bucket, set of 2 359151, standard bucket, set of 4	393266, standard bucket, set of 2 393267, standard bucket, set of 4
	GH-3.8A	SW	359487, Aerosolve, set of 2 358991, Aerosolve, set of 4	359488, Aerosolve, set of 2 358992, Aerosolve, set of 4
Allegra 21/R	F0685	FA	392270	392268
Spinchron 15/R	C0650	FA	—	Dedicated 50 ml rotor
	C1015	FA	Dedicated 15 ml rotor	—
	S4180	SW	361230	361234
Allegra X-22/R	SX4250	SW	392257	392258
	F0685	FA	392270	392268
	C0650	FA	—	Dedicated 50 ml rotor
	C1015	FA	Dedicated 15 ml rotor	—
Allegra 25R	TS 5.1-500	SW	368464	368461
	TA 10-250	FA	356964	356966
Allegra 64R	F0485	FA	392270	392268
	F0685	FA	392270	392268
	C0650	FA	—	Dedicated 50 ml rotor
	C1015	FA	Dedicated 15 ml rotor	—
J6 Large Cap.	JS-3.0, JS-4.0, JS-4.2, JS-5.2	SW	359151, set of 4 359472, set of 2	345386 —
	JS-7.5	SW	356964, 4-place adapter	356966, 1-place adapter 362213 (set of buckets), 3-place
J2 Series or Avanti J-25	JS-4.3	SW	359151, set of 4 359472, set of 2 359487, aerosol, set of 2 358991, aerosol, set of 4	359154, set of 4 359475, set of 2 359488, aerosol, set of 2 358992, aerosol, set of 4
		FA	356960, 4-place adapter	356965, 1-place adapter
		FA	—	Dedicated 50 ml rotor
		FA	356964, 4-place adapter	356966, 1-place adapter
	JA-12	FA	—	Dedicated 50 ml rotor
	JA-14	FA	356964, 4-place adapter	356966, 1-place adapter
	JA-18	FA	356962, 1 tube per adapter	356963, 1 tube per adapter
	F15B-8x50C**	FA	—	—
	F13B-14x50CY**	FA	—	—

* SW: Swinging bucket; FA: Fixed angle

** Manufactured by Piramoon Technology Inc.

Rotor/Adapter Selection Guide

Centrifuge	Rotor Name	Rotor Type*	Adapter for BD Falcon™ 15 ml Conical Tube	Adapter for BD Falcon™ 50 ml Conical Tube
Thermo Electron Corporation				
IEC	PR-7000M	FA	—	825, 8x50 ml
	3 Liter GP8F	SW	52165, 48x15 ml	52165, 20x50 ml
	3 Liter GP8F	SW	52285, 48x15 ml	52285, 20x50 ml
	3 Liter GP8	FA	—	825S, 8x50 ml
	1 Liter Multi	FA	—	8850, 8x50 ml
	1 Liter CL3R	SW	52437, TC 20x15 ml	52437, TC 8x50 ml
	1 Liter CL3	SW	—	52151, 4x50 ml***
	1 Liter CL3	FA	—	804S, 4x50 ml***
	1 Liter CL3	FA	58092, 12x15 ml, Centricon	—
	400ML CL2	SW	—	236, 4x50 ml
	400ML CL2	SW	—	52152, 8x50 ml
	400ML CL2	FA	—	58012, 6x50 ml
	400ML CL2	SW	—	52151, 4x50 ml
	400ML CL2	SW	52212, 6x15 ml	—
	400ML CL2	SW	52367, 8x15 ml	52367, 4x50 ml
Sorvall, Inc.				
TC-6	H-400	SW	78033	78035
GLC-2B/3 or	M	FA	03667	—
RC-3	SP/X	FA	00363	None required
	HL-4 w/omnicarrier	SW	00565	00648
	HL-4 w/carrier 00624	SW	—	00630
	HL-4 w/carrier 00634	SW	00363	—
RT-6000B/D or T-6000B/D	A-384	FA	None required	—
	A-500	FA	00363	None required
	A/S-400	FA	11148	None required
	H-1000B	SW	00884 or 11018 with 11152	00438 or 11148 with 11152
RT-7	RTH-250	SW	00884 or 11018 with 11152	00438 or 11148 with 11152
	RTH-750	SW	00447	00436-3 places or 00445-5 places
	SL-50T	FA	00402	—
RC-3B/C or	LA/S-400	FA	11148	None required
RC-3B/C Plus	H-2000B	SW	00884 or 11018 with 11152	00438 or 11148 with 11152
	HG-4L, H-4000, H-6000A	SW	00892	00436-3 places or 00445-5 places
Super T-21	SL-50T, SL-50RT	FA	00402	—
	SL-250T	FA	00456	03072
	STH-750	SW	00447	00436-3 places or 00445-5 places
RC-5B/C or	GSA, SLA-1000, SLA-1500	FA	00456	03072
RC-5BC/Plus or	SLA-600TC	FA	74232	None required
RC-24 or	GS-3, SLA-3000	FA	00456 and 00614	00614 and 03072
RC-26	SH-3000	SW	00447	00436-3 places or 00445-5 places
	HS-4 w/carrier 00479	SW	00456	03072
	HS-4 w/carrier 00481	SW	None required	None required
	HS-4 w/carrier 00480	SW	—	00363
	F15S-8x50C**	FA	—	—
	F13S-14x50CY**	FA	—	—
RC-28S	SL-250T	FA	00456	03072
	F-16/250	FA	00456	03072

3
Tubes

* SW: Swinging bucket; FA: Fixed angle

** Manufactured by Piramoon Technology Inc.

*** Requires 50968 Quick Release Shaft Adapter