

Sample tubes that can be used with the QIAAsymphony SP (software version 2.1) for user-validated protocols

 Recommended: These tubes are for use in combination with this protocol.
 User defined: User is responsible for defining and/or validating as the user deems appropriate.
 Not recommended: These tubes are not for use in combination with this protocol.
x /  Minimum sample volume (µl) required per sample per protocol (denoted by x) / Clot detection possible.
x /  Minimum sample volume (µl) required per sample per protocol (denoted by x) / Clot detection not possible.

Material	Supplier	Example cat. no.	Name in QIAAsymphony touchscreen	Insert	User-validated protocols										
					Cellfree protocols		Complex protocols								
					Cellfree500_V2	Cellfree200_FIX_V2	Complex200_V4	Complex 400_V2	Complex 800_V4	Complex200_FIX_V2	Complex400_FIX_V2	Complex800_FIX_V2	Complex200_OBL_V2	Complex400_OBL_V2	Complex800_OBL_V2
17 x 100 mm, 14 ml polystyrene round-bottom test tube	BD	352051	BD#352051 FalconPP 17x100	No insert needed	750 / ■		600 / □	800 / □	1200 / □						
Micro tube 2 ml with cap; Micro tube 2 ml, PP, SKIRTED	Sarstedt	72694	SAR#72.694 T2.0 ScrewSkirt	Insert 3B	600 / □	220 / □	300 / □	500 / □	950 / □	220 / □	420 / □	820 / □	630 / □ [†]	1040 / □ [†]	
Micro tube 2 ml with cap; Micro tube 2 ml, PP, NON-SKIRTED	Sarstedt	72.693	SAR#72.693 T2.0 Screw	Insert 3B/03	600 / □	220 / □	300 / □	500 / □	950 / □	220 / □	420 / □	820 / □	630 / □ [†]	1040 / □ [†]	
16 x 100 mm x 10 ml BD Vacutainer whole blood tube, K2-EDTA	BD	367525	BD#367525 VacutainerPP 16x100	No insert needed	□ [†]										
PAXgene Blood RNA Tube	BD	762165	BD#762165 PAXgeneRNA 16x100	No insert needed											
13 x 75 mm x 4 ml BD Vacutainer whole blood tube, K2-EDTA*	BD	367839	BD#367839 VacutainerK2 13x75	Insert 01/02/1A											
13 x 100 mm x 6 ml BD Vacutainer whole blood tube, K2-EDTA*	BD	367864	BD#367864 VacutainerK2 13x100	Insert 01/02/1A	□ [†]										
2.7 ml BD Citrate Tube 9NC	BD	363095	BD#363095 Vacutainer9NC 13x75	Insert 01/02/1A											
BD Vacutainer® PPT™ Plasma Preparation Tube for Molecular Diagnostic Test Methods	BD	362788	BD#362788 PPT 13x100	Insert 01/02/1A	□ [†]										
BD Microtainer™ Tubes with Microgard™ Closure	BD	365963; 365974; 365965; 365992	BD#365963 MicrotainerTube	Insert 3B											
Copan UTM 16 x 100 Tube 10 ml with 3 ml of UTM-RT medium*	Copan	330c; 302c; 303c; 328c; 335c; 346c; 307c; 305c; 338c	COP#330c UTM 16x100	No insert needed			600 / □	800 / □	1200 / □						
Copan E_Swab Tube	Copan	Part of E Swab Kits 480CE; 481CE; 482CE; 483CE; 484CE	COP#480CE E-SwabTube	Insert 01/1A			350 / □	550 / □							
Copan UTM 12 x 80 Tube with 1ml UTM Medium (with beads), 1 Regular FLOCKED Swab, Molded Breakpoint	Copan	359C; 361C; 360C	COP#359C UTM 12x80	Insert 01/1A			450 / □	650 / □							
Oragene DNA Self Collection Kit	DNA genotek	OG-500	DGT#OG-500 Oragene DANN	No insert needed			500 / □	700 / □	1100 / □						
GreinerBioOne454382Vacuette9NC13x75	Greiner Bio-One	454382	GR#454382 Vacuette9NC 13x75	Insert 01/02/1A											
Vacuette K3EDTA, 9 ml, 16 x 100 mm	Greiner Bio-One	455036	GR#455036 VacuettePP 16x100	No insert needed	■ [†]										

Table continued on next page

Table continued from previous page

Material	Supplier	Example cat. no.	Name in QIAAsymphony touchscreen	Insert	User-validated protocols										
					Cellfree protocols		Complex protocols								
					Cellfree500_V2	Cellfree200_FIX_V2	Complex200_V4	Complex 400_V2	Complex 800_V4	Complex200_FIX_V2	Complex 400_FIX_V2	Complex 800_FIX_V2	Complex200_OBL_V2	Complex400_OBL_V2	Complex800_OBL_V2
Vacurette K2-EDTA, 3 ml, 13 x 75 mm	Greiner Bio-One	454020	GR#454020 VacuretteK2 13x75	Insert 01/02/1A											
Vacurette K2-EDTA, 6 ml, 13 x 100 mm	Greiner Bio-One	456043	GR#456043 VacuretteK2 13x100	Insert 01/02/1A	□ [†]										
Nunc Cryo Tube 3.6 ml 12.5 x 72 mm	Nunc	379189	NU#379189 3.6Cryo 12.5x72	Insert 02/1A			450 / □	650 / □	1150 / □						
Nunc Cryo Tube 12.5 x 92 (4.5 ml)	Nunc	363452	NU#363452 4.5mlCryo 12.5x92	Insert 02/1A	700 / □		400 / □	600 / □	1000 / □					1040 / □ [‡]	1855 / □ [‡]
QIAGEN Collection Tube (2 ml)	QIAGEN	19201	QIA#19201 Collection	Insert 3B/03	□										
digene Hybrid Capture Specimen Transport Medium 16 x 73*	QIAGEN	5001-0461; 5001-0795; Part of 5123-1220 & 5122-1220	QIA#5001-0461 DigeneHC 16x73	No insert needed			300 / □	500 / □							
Roche STM (Specimen Transport Medium) 13 x 83 Tube	Roche	753920	RO#753920 STM 13x83	Insert 01/1A			450 / □	650 / □							
S-Monovette 4 ml, 15 x 75 mm, K2-EDTA	Sarstedt	3.1068.001	SAR#31068001 Monov 15x75	No insert needed											
S-Monovette 7.5 ml, 15 x 92 mm, Lithium-Heparin	Sarstedt	1.1608.001	SAR#11608001 Monov 15x92	No insert needed	■ [†]										
S-Monovette 9 ml, 16 x 92 mm, K2-EDTA	Sarstedt	2.1066.001	SAR#21066001 Monov16x92	No insert needed	■ [†]										
S-Monovette 2.6 ml, 13 x 65 mm, K2-EDTA	Sarstedt	4.1901.001	SAR#41901001 MonovK2 13x65	Insert 01/1A											
S-Monovette 4.9 ml, 13 x 90 mm, K2-EDTA	Sarstedt	4.1931.001	SAR#41931001 MonovK2 13x90	Insert 01/1A											
S-Monovette 2.7 ml, 11 x 66 mm, K2-EDTA	Sarstedt	5.1167.001	SAR#51167001 MonovK3 11x66	Insert 02/1A											
Sarstedt Tube round bottom13x82 7 ml reagent and centrifuge tube	Sarstedt	60550100	SAR#60550100 13x82 Round	Insert 01/1A	□		□	□	□						
Reagent and centrifuge tube 16.8 x 82	Sarstedt	55524	SAR#55524 Tube 16.8x82	No insert needed	800 / □		500 / □	700 / □	1100 / □						
Reagent and centrifuge tube 16.8 x 67	Sarstedt	55533	SAR#55533 Tube 16.8x67	No insert needed	800 / □		500 / □	700 / □	1100 / □						
Urine Monovette 10 ml 102 x 15 mm*	Sarstedt	10.252 or 10.253	SAR#10.252 UrineMonov	No insert needed			350 / □	550 / □	950 / □						
Micro tube 1,5 ml with cap	Sarstedt	72607	SAR#72.607 T1.5 Screw	Insert 3B	□		□	□	□						
Sarstedt polypropylene tube, 10 ml, round bottom	Sarstedt	62551201	SAR#62.551.201 T10.0 ScrewSkirt	No insert needed	■		450 / □	650 / □	1150 / □						
Reagent and centrifuge tube 3,5ml 66x11,5mm	Sarstedt	60549001	SAR#60.549.001 T3.5 ScrewSkirt	Insert 02/2A	700 / □		450 / □	650 / □	1150 / □						
Reagent and centrifuge tube 10ml 97x16PS+Cap	Sarstedt	609921829	SAR#60.9921.829 T10.0 ScrewSkirt V	No insert needed	800 / □		500 / □	700 / □	1100 / □						
1.5 ml Plain Skirted Tube, Natural Standard Screw Cap	Starlab	E1415-2240	SL#E1415-2240 T1.5 ScrewSkirt	Insert 3B	□		□	□	□						
Venosafe K2-EDTA, 9 ml	Terumo	VF-109SDK	TER#VF-109SDK Venosafe 16x100	No insert needed	■ [†]										
Venoject K3EDAT, 3 ml	Terumo	VT-053STK	TER#VT-053STK VenojectK2 13x75	Insert 02/1A											
Venosafe K2-EDTA 5.5 ml	Terumo	VF-076SDK	TER#VF-076SDK VenosafeK2 13x100	Insert 02/1A/2A											

* This tube type may fit loosely in the tube insert(s) 01 or 02, which can result in movement of the tubes during sample input. We recommend the new insert 1A or 2A.

[†] This tube must be filled with the nominal volume for blood, as recommended by the manufacturer. Centrifugation must be carried out in a swing-rotor, according to the manufacturer's instructions. In some cases, the liquid level above the layer of gel or above the interface may not be sufficient for sample transfer, although it may appear so by eye. In this case, the plasma must be manually transferred to a secondary tube that can then be placed onto the QIAAsymphony SP.

[‡] For the OBL (Off-board lysis) protocols the required volume is defined by the lysate volume generated during manual lysis.



For up-to-date licensing information and product-specific disclaimers, see the respective QIAGEN kit handbook or user manual. QIAGEN kit handbooks and user manuals are available at www.qiagen.com or can be requested from QIAGEN Technical Services or your local distributor.

© 2010 QIAGEN, all rights reserved.