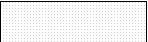


Bravo Vacuum Filtration Station Labware Compatibility Matrix

This matrix lists labware types that Agilent Technologies has tested with the Bravo Vacuum Filtration Station assembled in Configuration A or Configuration B. For each combination of collection plate and filter plate, the matrix specifies the requirements for the collar gasket type (blue or black) and the number of spacer inserts to include in the base, if any, to adjust the spacing of the collection plate. This matrix shows the mechanical hardware compatibility only.

Collection Plates (Part Number)	Filter Plates (Part Number)																
	Agilent				Millipore				MN		Qiagen		Waters	Pall			
	Captiva (A5969045)	Bond Elut Plexa (A4968030)	Bond Elut Square (A3960150)	Versaplate (75401025)	Solvinert 96 PP (MSRLN04)	Solvinert 96 PS (MSHVN45)	Solvinert Deep Well (MDRPNP4)	Multiscreen HTS 384 (MZHV)	Chromafil (738662.M)	Nucleospin (740625.1)	Qiaprep (120012)	Turbofilter (120025)	uElution (186001828BA)	Acroprep 1mL (8247)	Acroprep 350uL (8036)	Acroprep 384 Short Tip	Acroprep 384 Long
Captiva (Agilent A8690D1000)	Collar: Tall Insert: None Gasket: Black Configuration: B	Collar: Tall Insert: None Gasket: Black Configuration: B	Collar: Tall Insert: 1 Tall, 1 Short Gasket: Single Black Configuration: B (Configuration A not recommended)	Collar: Tall Insert: 2 Short Gasket: Black Configuration: B	Collar: Tall Insert: None Gasket: Blue Configuration: A (Configuration B not recommended)	Collar: Tall Insert: None Gasket: Blue Configuration: A (Configuration B not recommended)			Collar: Tall Insert: None Gasket: Black Configuration: B	Collar: Tall Insert: None Gasket: Black Configuration: B			Collar: Tall Insert: None Gasket: Black Configuration: B		Collar: Tall Insert: None Gasket: Blue Configuration: A (Configuration B not recommended)		
Eppendorf 96/500 (0030 507.103)	Collar: Tall Insert: 1 Tall, 1 Short Gasket: Black Configuration: B	Collar: Tall Insert: 1 Tall, 1 Short Gasket: Black Configuration: B	Collar: Short Insert: None Gasket: Single Black Configuration: B (Configuration A not recommended)		Collar: Tall Insert: 1 Tall, 1 Short Gasket: Blue Configuration: A (Configuration B not recommended)	Collar: Tall Insert: 1 Tall, 1 Short Gasket: Blue Configuration: A (Configuration B not recommended)	Collar: Tall Insert: 1 Tall, 1 Short Gasket: Blue Configuration: A (Configuration B not recommended)		Collar: Tall Insert: 1 Tall, 1 Short Gasket: Black Configuration: B	Collar: Tall Insert: 1 Tall, 1 Short Gasket: Black Configuration: B	Collar: Tall Insert: 1 Tall 1 Short Gasket: Black Configuration: B	Collar: Tall Insert: 1 Tall 1 Short Gasket: Black Configuration: B	Collar: Tall Insert: 1 Tall, 1 Short Gasket: Black Configuration: B		Collar: Tall Insert: 1 Tall, 1 Short Gasket: Blue Configuration: A (Configuration B not recommended)		
Eppendorf 96/1000 (0030 507.200)	Collar: Tall Insert: None Gasket: Black Configuration: B	Collar: Tall Insert: None Gasket: Black Configuration: B	Collar: Tall Insert: 6 Short Gasket: Single Black Configuration: B	Collar: Tall Insert: 2 Short Gasket: Black Configuration: B					Collar: Tall Insert: None Gasket: Black Configuration: B	Collar: Tall Insert: None Gasket: Black Configuration: B			Collar: Tall Insert: None Gasket: Black Configuration: B				
Nunc (260252)	Collar: Tall Insert: 1 Tall Gasket: Black Configuration: B	Collar: Tall Insert: 1 Tall Gasket: Black Configuration: B		Collar: Tall Insert: 1 Tall, 1 Short Gasket: Black Configuration: B	Collar: Tall Insert: 5 Short Gasket: Blue Configuration: A (Configuration B not recommended)	Collar: Tall Insert: 5 Short Gasket: Blue Configuration: A (Configuration B not recommended)	Collar: Tall Insert: 4 Short Gasket: Blue Configuration: A (Configuration B not recommended)		Collar: Tall Insert: 1 Tall Gasket: Black Configuration: B	Collar: Tall Insert: 1 Tall Gasket: Black Configuration: B			Collar: Tall Insert: 1 Tall Gasket: Black Configuration: B		Collar: Tall Insert: 5 Short Gasket: Blue Configuration: A (Configuration B not recommended)		
Costar (Corning 3959)	Collar: Tall Insert: None Gasket: Black Configuration: B	Collar: Tall Insert: None Gasket: Black Configuration: B		Collar: Tall Insert: 3 Short Gasket: Black Configuration: B	Collar: Tall Insert: None Gasket: Blue Configuration: A (Configuration B not recommended)	Collar: Tall Insert: None Gasket: Blue Configuration: A (Configuration B not recommended)			Collar: Tall Insert: None Gasket: Black Configuration: B	Collar: Tall Insert: None Gasket: Black Configuration: B	Collar: Tall Insert: None Gasket: Black Configuration: B	Collar: Tall Insert: None Gasket: Black Configuration: B	Collar: Tall Insert: None Gasket: Black Configuration: B		Collar: Tall Insert: None Gasket: Blue Configuration: A (Configuration B not recommended)		
Greiner U-bottom (650101)	Collar: Short Insert: None Gasket: Black Configuration: B (Limited to 100 µL)	Collar: Short Insert: None Gasket: Black Configuration: B (Limited to 100 µL)	Collar: Short Insert: 1 Tall Gasket: Single Black Configuration: B (Limited to 100 µL)	Collar: Short Insert: 2 Short Gasket: Black Configuration: B	Collar: Short Insert: None Gasket: Blue Configuration: A (Configuration B not recommended)	Collar: Short Insert: None Gasket: Blue Configuration: A (Configuration B not recommended)	Collar: Short Insert: 4 Short Gasket: Blue Configuration: A (Configuration B not recommended)		Collar: Short Insert: None Gasket: Black Configuration: B (Limited to 100 µL)	Collar: Short Insert: None Gasket: Black Configuration: B (Limited to 100 µL)			Collar: Short Insert: None Gasket: Black Configuration: B (Limited to 100 µL)	Collar: Tall Insert: 4 Short Gasket: Blue Configuration: A	Collar: Short Insert: None Gasket: Blue Configuration: A (Configuration B not recommended)		
Greiner 384 (781101)								Collar: Short Insert: None Gasket: Blue Configuration: A (Configuration B not recommended)								Collar: Short Insert: None Gasket: Blue Configuration: A (Configuration B not recommended)	Collar: Short Insert: None Gasket: Blue Configuration: A (Configuration B not recommended)



Not supported



Configuration A cannot be auto-assembled for this labware combination. The configuration must be pre-assembled at the Bravo deck location. In addition, the Bravo device profile must specify the **Always move to safe Z before each process** option.

Notes: The blue gasket in the collar may be replaced by two black gaskets to accommodate filter plates that cannot be seated in the blue gasket, such as the Agilent Captiva and Bond Elut plates. In this case, one black gasket adheres to the top of the collar and a second black gasket adheres to the underside of the collar.

The Waters HLB Extraction plate (part number 186000309) and Pall Acroprep 2 mL plate (part number 8130) are not compatible with Configurations A and B of the Bravo Vacuum Filtration Station.