

Description	Unit	Part Number	Price (\$)
RRHT-1100 Modification Kit For Diode Array Detector	1 kit	5188-5324	2400.00
Kit includes:			
1 ea. 80µL low volume mixer		1 ea. Semi-Micro 5.33µL flow cell	
2 ea. 0.17mm x 400mm Stainless steel capillary		1 pk. 1/16" SS Fittings w ferrules (10)	
1 ea. 0.17mm x 200mm stainless steel capillary		1 ea. 4.6 x 50mm 1.8µm Eclipse XDB-C18 Column	
1 ea. 0.17mm x 105mm stainless steel capillary		1 ea. 4.6 x 50mm, 1.8µm Extend-C18 Column	
		1 ea. 4.6 x 50mm, 1.8µm StableBond C18 Column	

Description	Unit	Part Number	Price (\$)
RRHT-1100 Modification Kit For Diode Array-Mass Spec Detection	1 kit	5188-5328	2400.00
Kit includes:			
1 ea. 80mL low volume mixer		1 ea. Micro 1.7µL flow cell	
1 ea. 0.12mm x 280mm stainless steel capillary		1 pk. 1/16" SS Fittings w ferrules (10)	
1 ea. 0.12mm Needle Seat Std. Autosampler		1 ea. 2.1 x 50mm 1.8µm Eclipse XDB-C18 Column	
1 ea. 0.12mm Needle Seat Well Plate Autosampler		1 ea. 2.1 x 50mm, 1.8µm Extend-C18 Column	
1 ea. 0.12mm x 70mm stainless steel capillary		1 ea. ZDV stainless steel union	
1 ea. 0.12mm x 105mm stainless steel capillary			
1 ea. 0.12mm x 150mm stainless steel capillary			