

Product Number:	103400
Lot Number:	503285
Expiration:	27-Jun-2022
Platform(s):	SR-X

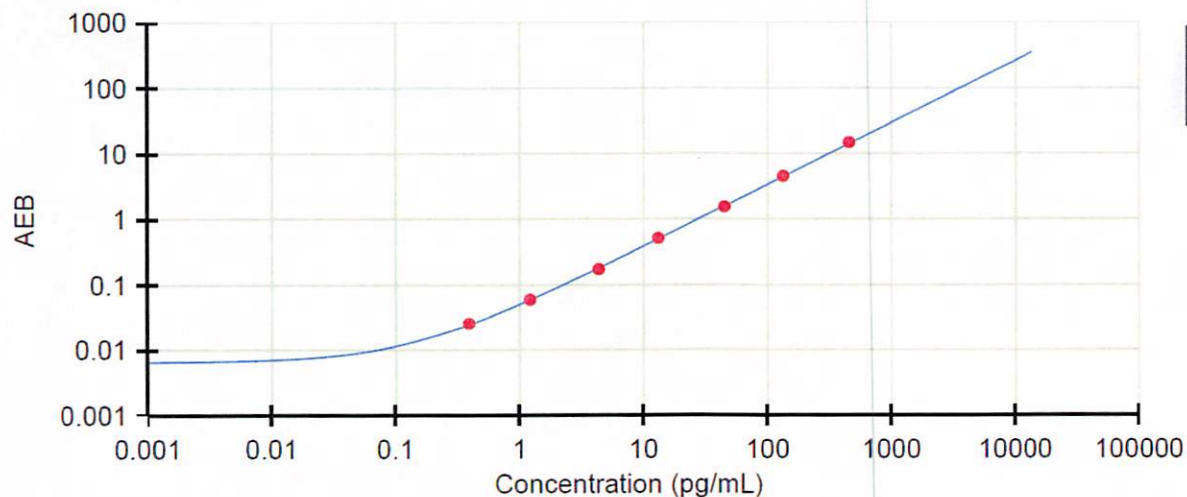


Data below represents results generated on the Quanterix SR-X™ Ultra-Sensitive

Component	Part Number	Lot Number
Bead Reagent	102246	117801
Detector Reagent	102248	117803
SBG Reagent	102250	117808
Sample Diluent	103399	135010
Calibrators	102255	134704
Control 1	102256	134705
Control 2	102257	134705
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

Release Materials				Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Control 1	134705	3.74	3.20-4.80	
Control 2	134705	290	224-336	
Panel 1	134208	6.13	3.95-9.43	
Panel 2	134208	2.95	1.86-7.15	
Panel 3	134208	60.5	49.9-74.8	
Panel 4	135801	368	274-411	



Review/Approval Shivam Goel QC Shivam Goel 13 Jun 2022