

Product Number:	104073
Lot Number:	503808
Expiration:	7-Feb-2024
Platform(s):	HD-1 / HD-X / SR-X



Data below represents results generated on the Simoa HD-X® Analyzer.

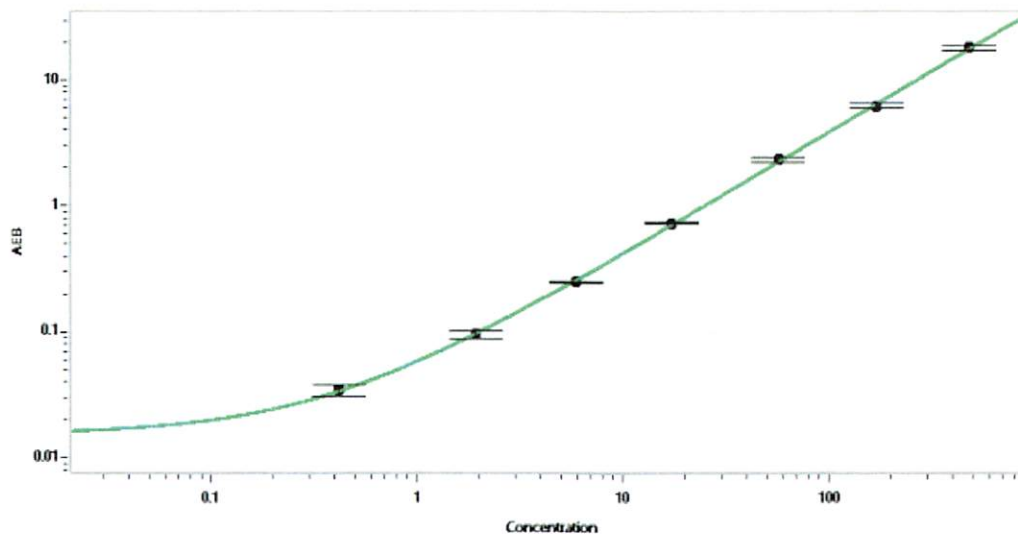
Component	Part Number	Lot Number
Bead Reagent	104067	303809
Detector Reagent	104068	303905
SBG Reagent	104069	304010
Sample Diluent	104071	304005
Calibrators	102255	303214 / 307909
Control 1	102256	303215
Control 2	102257	303215
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	303215	5.43	4.42 - 6.63
Control 2	303215	247	190 - 286
Panel 1	224402	9.56	8.84 - 13.3
Panel 2	224402	6.49	6.38 - 9.57
Panel 3	224402	71.4	63.0 - 94.5
Panel 4	212511	262	195 - 356

* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.416
C	1.94
D	5.90
E	17.4
F	56.9
G	170
H	482

Review/Approval Matt Koulouris QC Supervisor [Signature] 29 Mar 2023
 Name Title Signature/Date