

Product Number:	104073
Lot Number:	503867
Expiration:	7-Aug-2024
Platform(s):	HD-1 / HD-X / SR-X



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

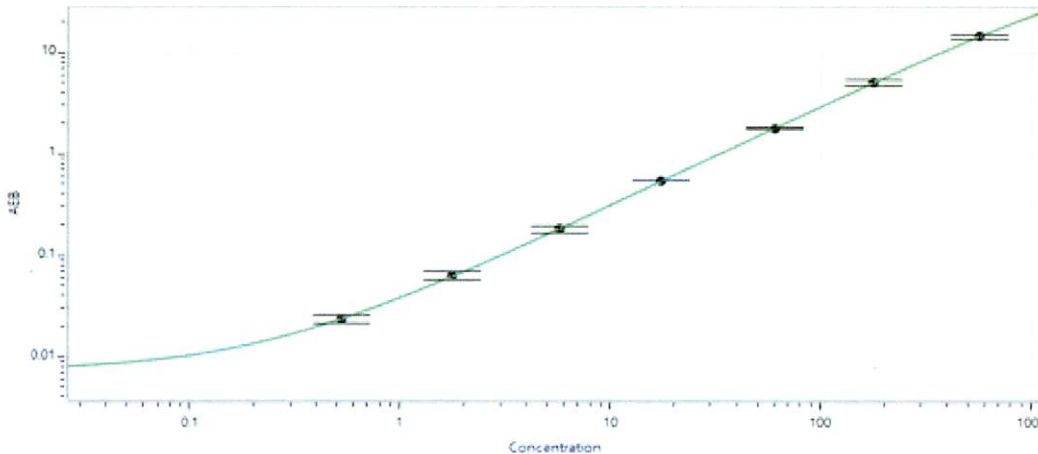
Component	Part Number	Lot Number
Bead Reagent	104067	321910
Detector Reagent	104068	322002
SBG Reagent	104069	322001
Sample Diluent	104071	322606
Calibrators	102255	323302
Control 1	102256	323303
Control 2	102257	323303
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	323303	14.0	10.8 - 17.1
Control 2	323303	612	514 - 771
Panel 1	224402	10.2	7.86 - 11.8
Panel 2	224402	7.02	5.60 - 8.59
Panel 3	224402	79.5	56.1 - 84.1
Panel 4	301701	614	462 - 719

* 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	0.525
C	1.75
D	5.72
E	17.3
F	59.8
G	176
H	564

Review/Approval Sonia Khillan QC Sonia Khillan 07SEP2023
 Name Title Signature/Date