

Product Number:	104111
Lot Number:	503769
Expiration:	9-Jan-2024
Platform(s):	HD-1 / HD-X & SR-X



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

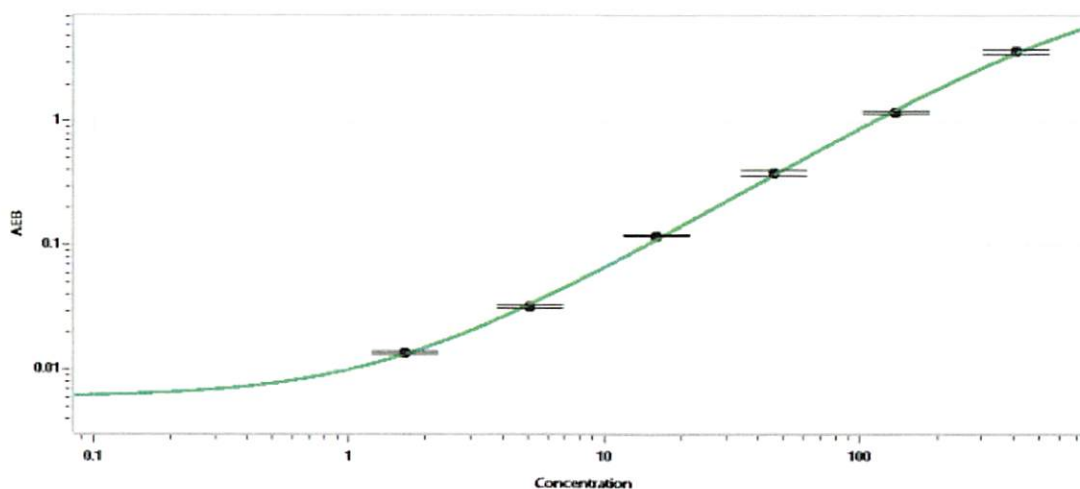
Component	Part Number	Lot Number
Bead Reagent	104098	300908
Detector Reagent	104099	309606
SBG Reagent	104100	301202
Sample Diluent	104101	301203
Reference Calibrator	104106	301806
Control 1	104107	301807
Control 2	104108	303102
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	301807	40.4	31.5 - 47.2
Control 2	303102	785	646 - 968
Panel 1	301702	16.8	12.7 - 20.9
Panel 2	301702	16.2	12.3 - 20.9
Panel 3	214602	647	535 - 802
Panel 4	301702	86.4	57.9 - 125

* 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	1.66
C	5.13
D	16.0
E	45.6
F	138
G	408

Review/Approval

Sonia Khillan

Name

QC

Title

Sanjiv Kella 22 MAY 2023

Signature/Date