

Product Number:	103186
Lot Number:	502433
Expiration:	27-May-2021
Platform(s):	HD-1/ HD-X



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

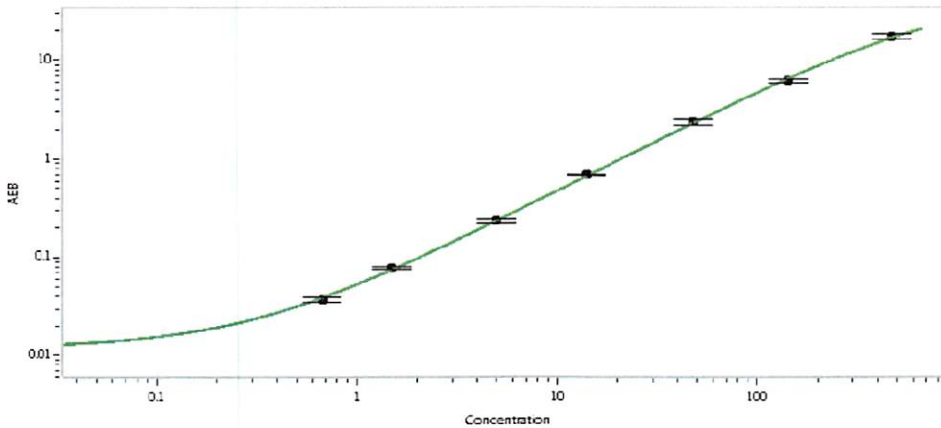
Component	Part Number	Lot Number
Bead Reagent	102246	017009
Detector Reagent	102248	017004
SBG Reagent	102250	017007
Sample Diluent	102252	017008
Calibrators	102255	018301
Control 1	102256	018401
Control 2	102257	018401
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	018401	4.32	3.30-4.95
Control 2	018401	148	117-176
Panel 1	922601	18.8	14.1 - 21.1
Panel 2	922602	6.76	4.94 - 7.41
Panel 3	922603	40.9	28.4 - 42.6
Panel 4	922604	167	129 - 204

* 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	0.673
C	1.48
D	4.94
E	13.9
F	47.3
G	142
H	472

Review/Approval	Shivani Goel	QC Manager	Shivani Goel 13/01/2020
	Name	Title	Signature/Date