

Product Number:	103186
Lot Number:	502183
Expiration:	10-Dec-2020
Platform(s):	HD-1/HD-x



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

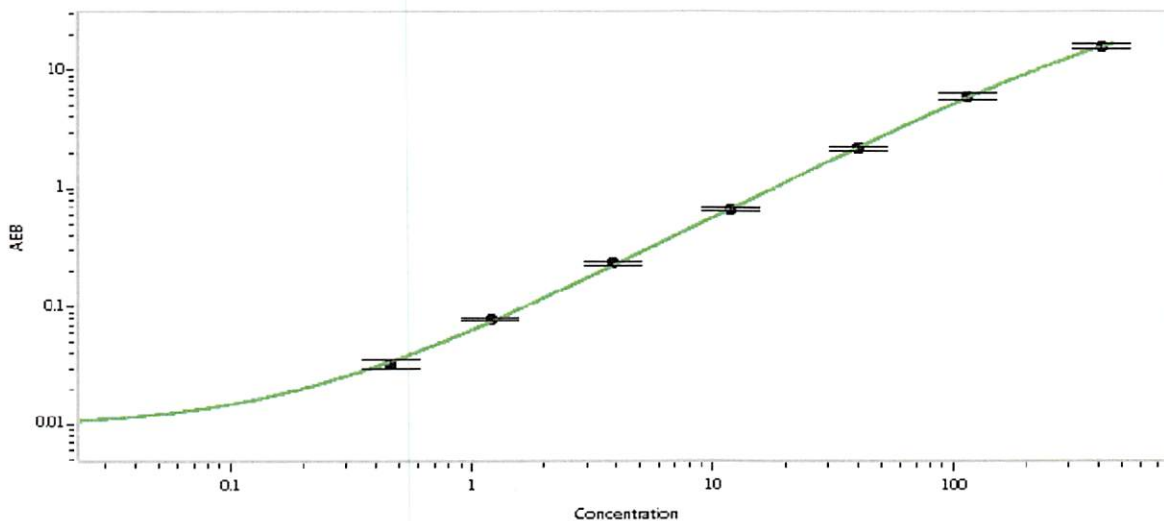
Component	Part Number	Lot Number
Bead Reagent	102246	934507
Detector Reagent	102248	934405
SBG Reagent	102250	934404
Sample Diluent	102252	934403
Calibrators	102255	935001
Controls	102256 - 102257	935701
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	935701	3.88	3.21-4.86
Control 2	935701	159	134-201
Panel 1	922601	16.2	14.1-21.1
Panel 2	922602	6.10	4.94-7.41
Panel 3	922603	33.9	28.4-42.6
Panel 4	922604	153	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	0.461
C	1.20
D	3.87
E	11.8
F	40.2
G	115
H	415

Review/Approval: Patrizia Stadler Assoc. Director, QC *p. Stadler* 13 Jan 2020  
 Name Title Signature/Date