

Product Number:	103186
Lot Number:	502186
Expiration:	27-Dec-2020
Platform(s):	HD-1/HD-x SRx



Data below represents results generated on the Simoa™ HD-1 Analyzer.

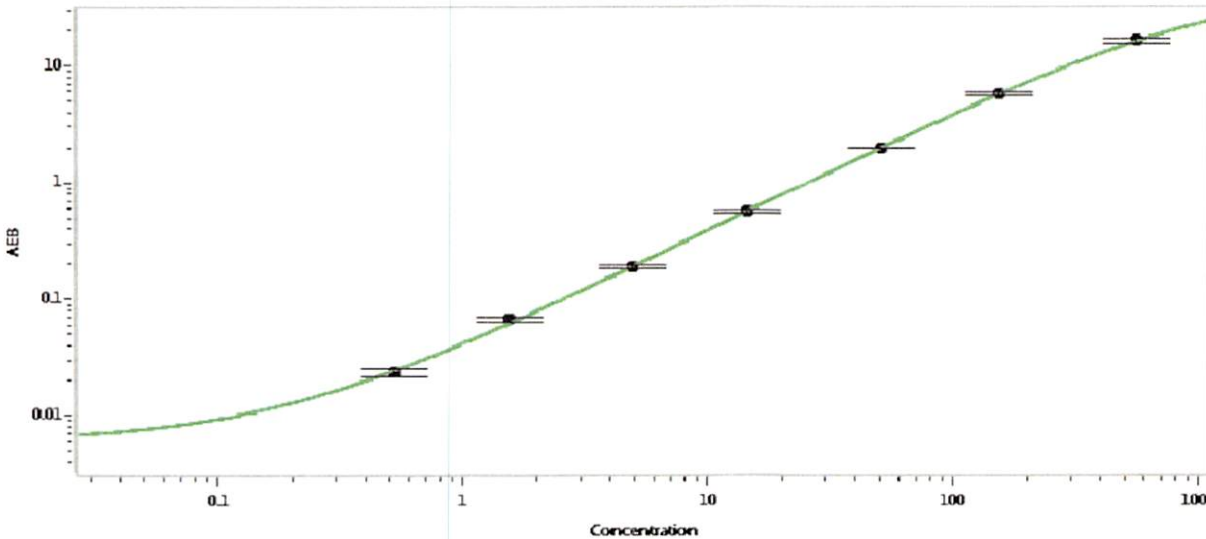
Component	Part Number	Lot Number
Bead Reagent	102246	001005
Detector Reagent	102248	000703
SBG Reagent	102250	000702
Sample Diluent	102252	000705
Calibrators	102255	000801
Controls	102256 - 102257	000901
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	000901	4.41	3.08-5.03
Control 2	000901	165	127-190
Panel 1	922601	19.8	14.1-21.1
Panel 2	922602	7.19	4.94-7.41
Panel 3	922603	42.0	28.4-42.6
Panel 4	922604	190	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	0.525
C	1.54
D	4.91
E	14.5
F	51.1
G	154
H	561

Review/Approval: Patrizia Stadler Assoc. Director, QC *P. Stadler* 14 Jan 2020
Name Title Signature/Date