

Product Number:	103186
Lot Number:	502268
Expiration:	12-Feb-2021
Platform(s):	HD-1/HD-x, SRx



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

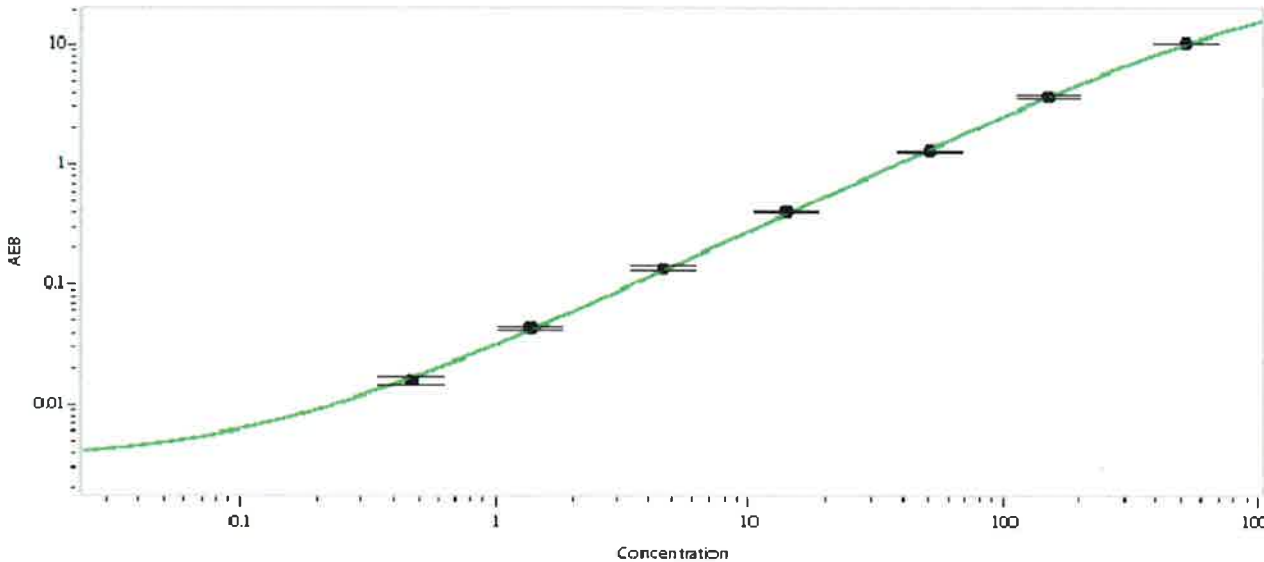
Component	Part Number	Lot Number
Bead Reagent	102246	006304
Detector Reagent	102248	006206
SBG Reagent	102250	006205
Sample Diluent	102252	006310
Calibrators	102255	007001
Controls	102256 - 102257	007002
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	007002	3.95	2.69-4.51
Control 2	007002	153	122-183
Panel 1	922601	18.1	14.1-21.1
Panel 2	922602	6.30	4.94-7.41
Panel 3	922603	37.4	28.4-42.6
Panel 4	922604	160	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	0.474
C	1.36
D	4.58
E	14.1
F	51.4
G	152
H	526

Review/Approval: Patrizia Stadler Assoc. Director, QC *P. Stadler* 17 MAR 2020
Name Title Signature/Date