

Product Number:	103186
Lot Number:	502296
Expiration:	27-Jan-2021
Platform(s):	HD-1/HD-x, SRx



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

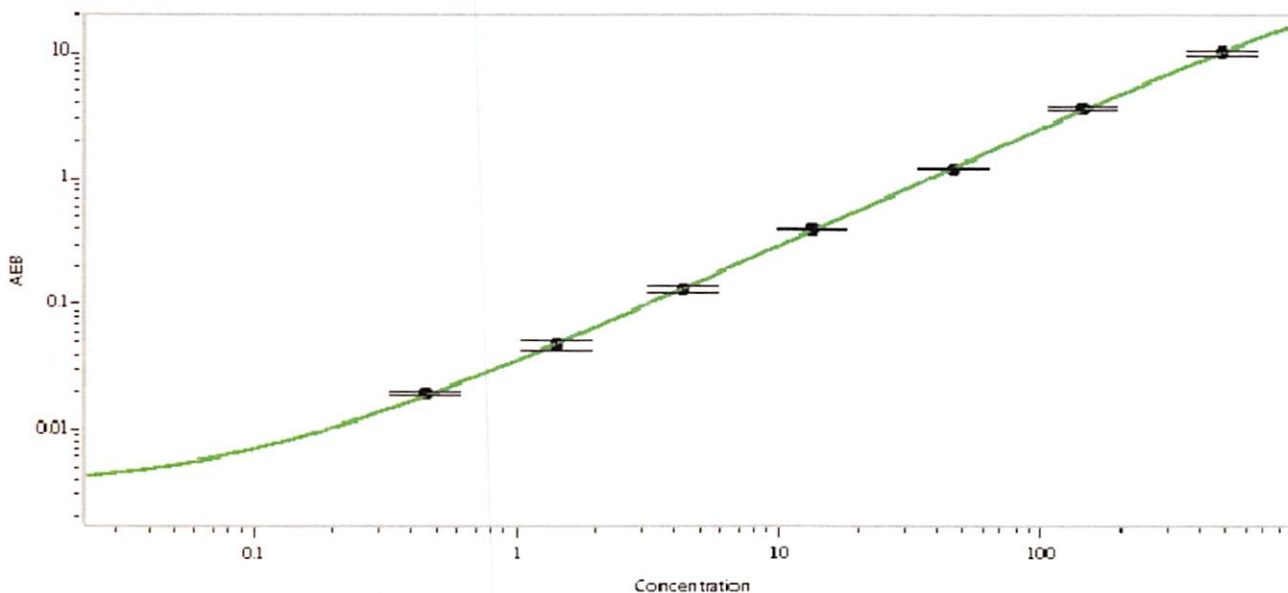
Component	Part Number	Lot Number
Bead Reagent	102246	007811
Detector Reagent	102248	007905
SBG Reagent	102250	007907
Sample Diluent	102252	002704
Calibrators	102255	008001
Controls	102256 - 102257	008002
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	008002	3.39	2.43-4.30
Control 2	008002	160	123-184
Panel 1	922601	15.6	14.1-21.1
Panel 2	922602	5.48	4.94-7.41
Panel 3	922603	37.1	28.4-42.6
Panel 4	922604	168	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	0.451
C	1.42
D	4.33
E	13.3
F	46.2
G	142
H	486

Review/Approval: Patrizia Stadler Name, Assoc. Director, QC Title, P. Stadler Signature, 31 Dec 2020 Date