

Product Number:	103186
Lot Number:	501548
Expiration:	2-May-2019
Platform(s):	HD-1 & SR-X



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

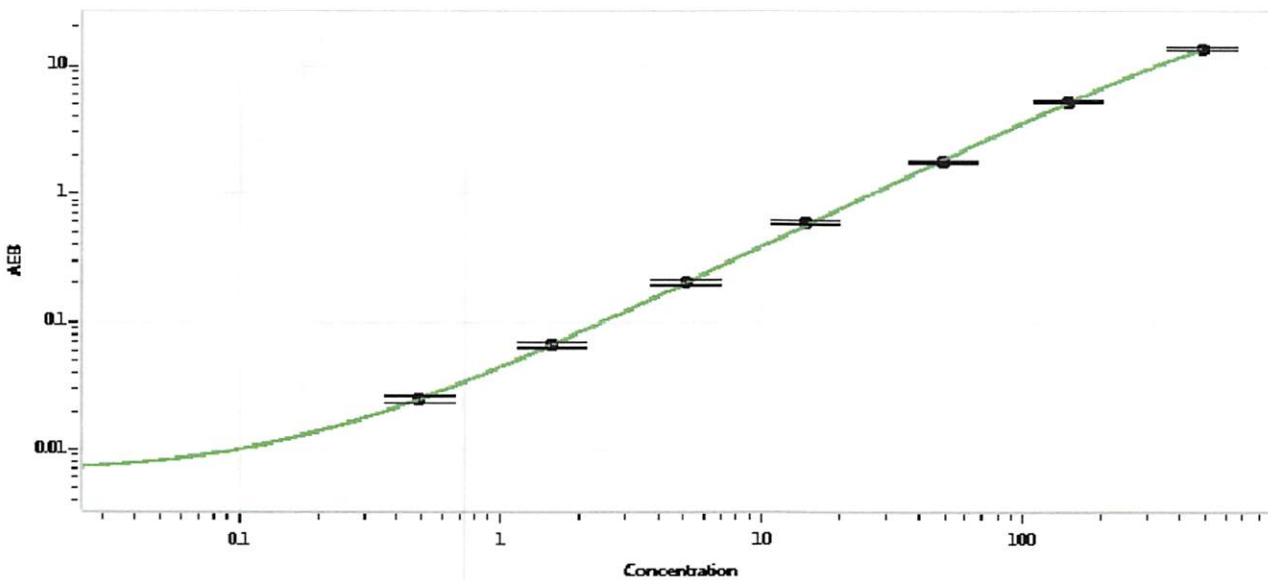
Component	Part Number	Lot Number
Bead Reagent	102246	815503
Detector Reagent	102248	814505
SBG Reagent	102250	812211
Sample Diluent	102252	812212
Reference Calibrators	102255	831901
Controls	102256 - 102257	822806 830201
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	822806	5.04	3.63-5.71
Control 2	830201	167	125-187
Panel 1-A	823301	5.54	4.08-6.68
Panel 1-B	823302	6.56	4.90-7.61
Panel 2	823601	87.6	59.3-101
Panel 3	823602	342	236-402

* 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	0.494
C	1.58
D	5.10
E	14.8
F	49.3
G	150
H	484

Review/Approval	Patrizia Stadler	Associate Director, QC	<i>P. Stadler</i>	14 Jan 2019
	Name	Title	Signature/Date	