

Product Number:	103186
Lot Number:	501814
Expiration:	3-Apr-2020
Platform(s):	HD-1 & SR-X



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

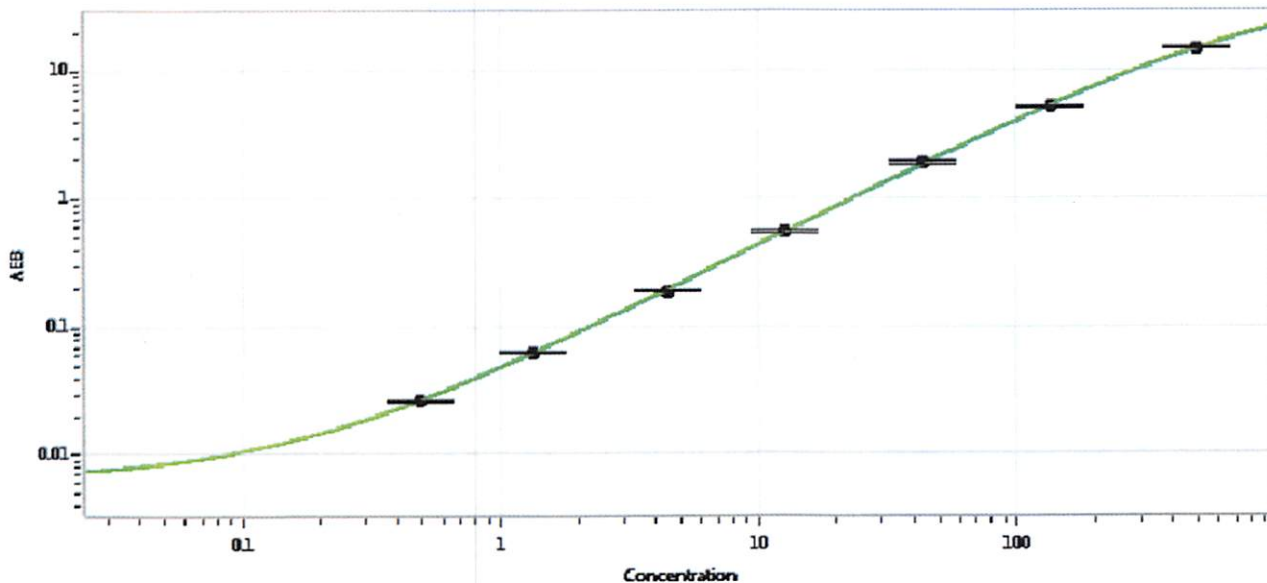
Component	Part Number	Lot Number
Bead Reagent	102246	909315
Detector Reagent	102248	910008
SBG Reagent	102250	909314
Sample Diluent	102252	909302
Reference Calibrators	102255	911901
Controls	102256 - 102257	911902
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	911902	3.99	3.04-4.81
Control 2	911902	153	122-184
Panel 1-A	823301	4.79	4.08-6.68
Panel 1-B	823302	5.39	4.90-7.61
Panel 2	823601	73.7	59.3-101
Panel 3	823602	270	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	0.485
C	1.33
D	4.43
E	12.7
F	43.6
G	136
H	498

Review/Approval: Patrizia Stadler QC Assoc. Director *P. Stolle* 06MAY2019
Name Title Signature/Date