

Product Number:	103186
Lot Number:	501946
Expiration:	16-Feb-2020
Platform(s):	HD-1/ HD-X & SR-X



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

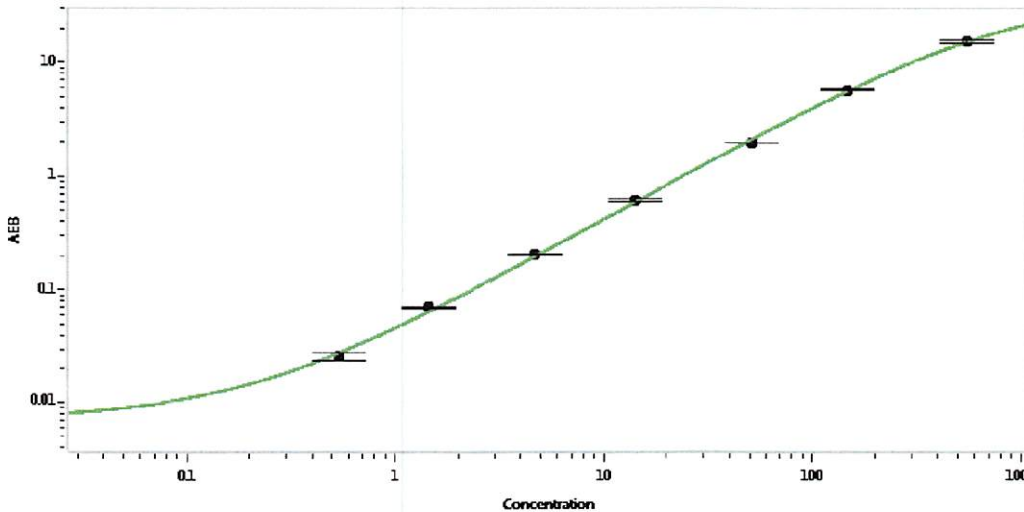
Component	Part Number	Lot Number
Bead Reagent	102246	906503
Detector Reagent	102248	904703
SBG Reagent	102250	904702
Sample Diluent	102252	904701
Calibrators	102255	910701
Controls	102256-102257	910702
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	910702	6.32	4.38-6.78
Control 2	910702	164	127-191
Panel 1	823301	5.29	4.08-6.68
Panel 2	823302	6.12	4.90-7.61
Panel 3	823601	76.8	59.3-101
Panel 4	823602	305	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	0.535
C	1.46
D	4.70
E	14.2
F	50.9
G	147
H	549

Review/Approval

Patrizia Stadler Assoc. Director, QC

Name

Title

Signature/Date