

Product Number:	103186
Lot Number:	501942
Expiration:	16-Feb-2020
Platform(s):	HD-1/ HD-X & SR-X



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

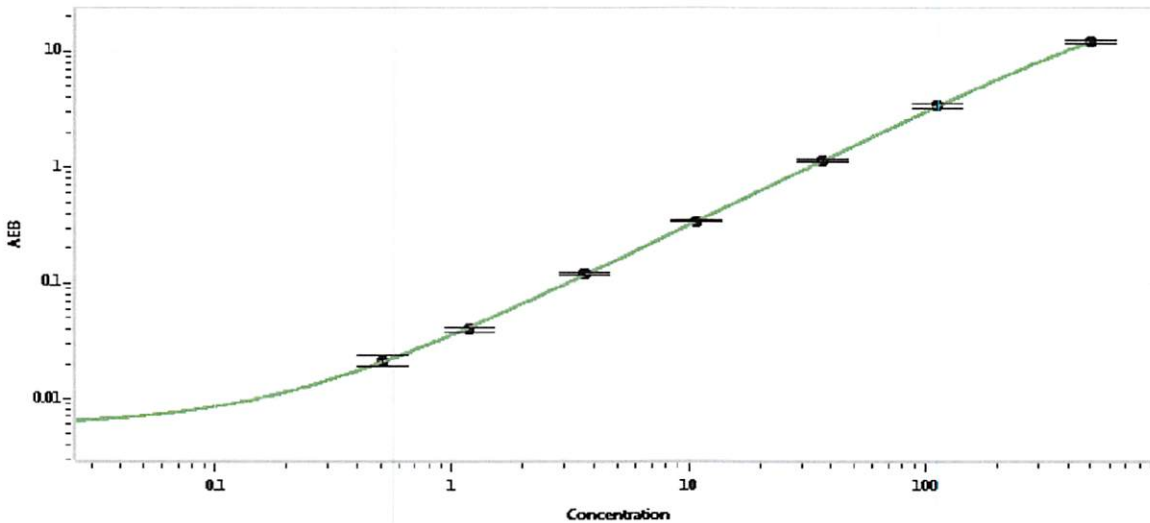
Component	Part Number	Lot Number
Bead Reagent	102246	906503
Detector Reagent	102248	904703
SBG Reagent	102250	904702
Sample Diluent	102252	904701
Reference Calibrators	102255	913501
Controls	102256 - 102257	913601
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	913601	4.70	3.56-5.35
Control 2	913601	196	145-218
Panel 1	823301	4.95	4.08-6.68
Panel 2	823302	6.19	4.90-7.61
Panel 3	823601	80.4	59.3-101
Panel 4	823602	313	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	0.509
C	1.18
D	3.63
E	10.8
F	36.7
G	112
H	493

Review/Approval

Patrizia Stadler Assoc. Director, QC
Name Title

P. Stadler

18 Sep 2019

Signature/Date