

Product Number:	103186
Lot Number:	501892
Expiration:	16-Feb-2020
Platform(s):	HD-1/HD-x & SR-X



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

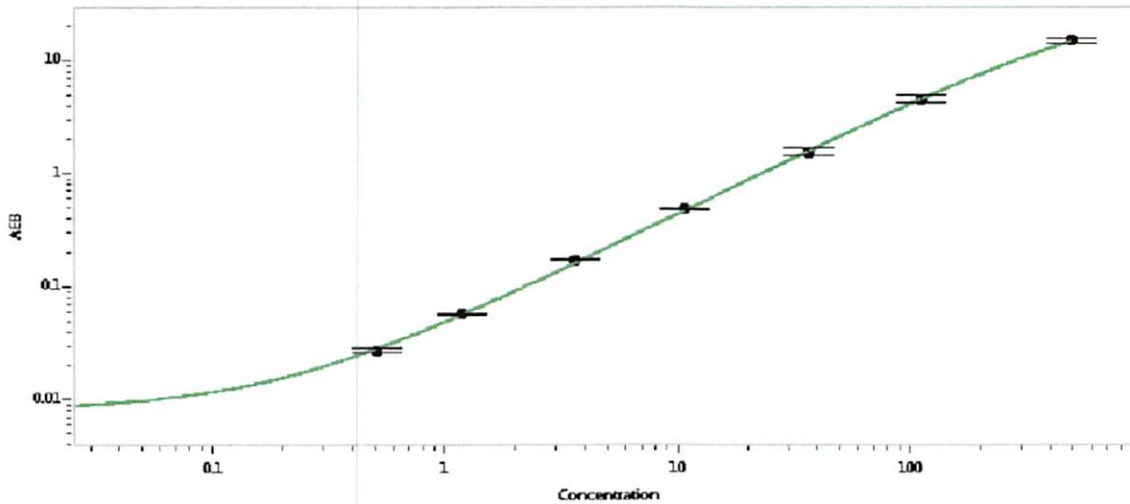
Component	Part Number	Lot Number
Bead Reagent	102246	911413
Detector Reagent	102248	910508
SBG Reagent	102250	910509
Sample Diluent	102252	904701
Calibrators	102255	913501
Controls	102256 - 102257	822806, 913601
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	822806	5.14	3.63-5.71
Control 2	913601	184	145-218
Panel 1a	823301	5.20	4.08-6.68
Panel 1b	823302	5.89	4.90-7.61
Panel 2	823601	77.4	59.3-101
Panel 3	823602	285	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 P. Stadler 08 Jul 2019
 Name Title Signature/Date