

Product Number:	103186
Lot Number:	501591
Expiration:	4-Jun-2019
Platform(s):	HD-1 & SR-X



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

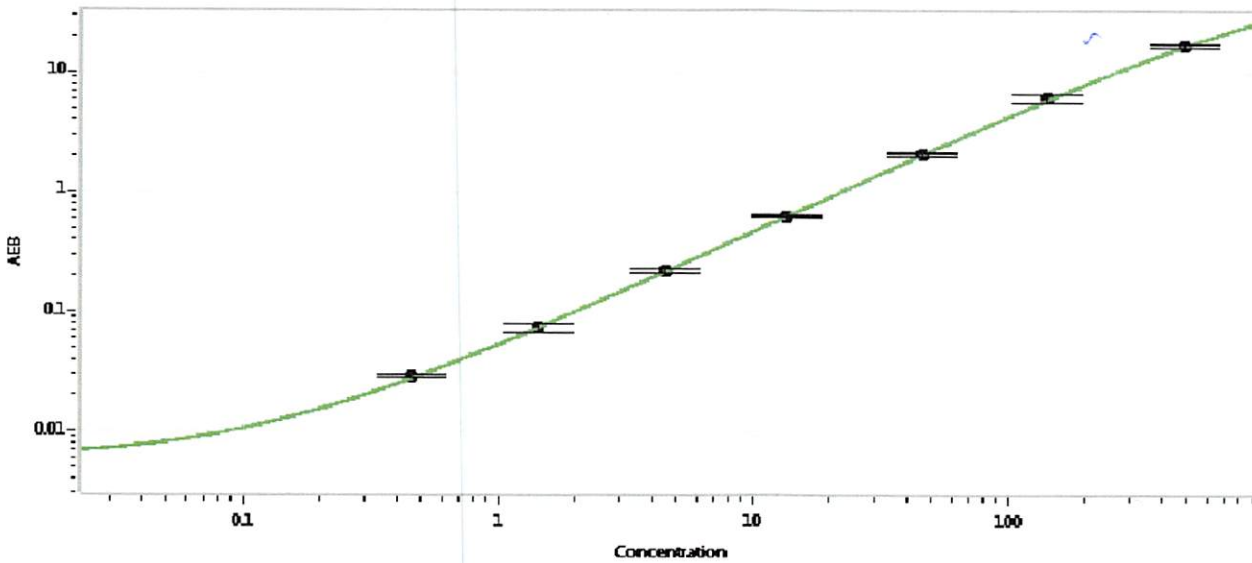
Component	Part Number	Lot Number
Bead Reagent	102246	815503
Detector Reagent	102248	824203
SBG Reagent	102250	833911
Sample Diluent	102252	834507
Reference Calibrators	102255	901401/833201
Controls	102256 - 102257	822806/830201
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	822806	4.90	3.63-5.71
Control 2	830201	159	125-187
Panel 1-A	823301	5.37	4.08-6.68
Panel 1-B	823302	6.36	4.90-7.61
Panel 2	823601	81.8	59.3-101
Panel 3	823602	314	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	0.458
C	1.44
D	4.48
E	13.5
F	45.9
G	142
H	488

Review/Approval

Patrizia Stadler  
Name

Associate Director,  
QC

Title

*P Stadler* 18 Jan 2019

Signature/Date