

Product Number:	103186
Lot Number:	501603
Expiration:	15-Jan-2020
Platform(s):	HD-1 & SR-X



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

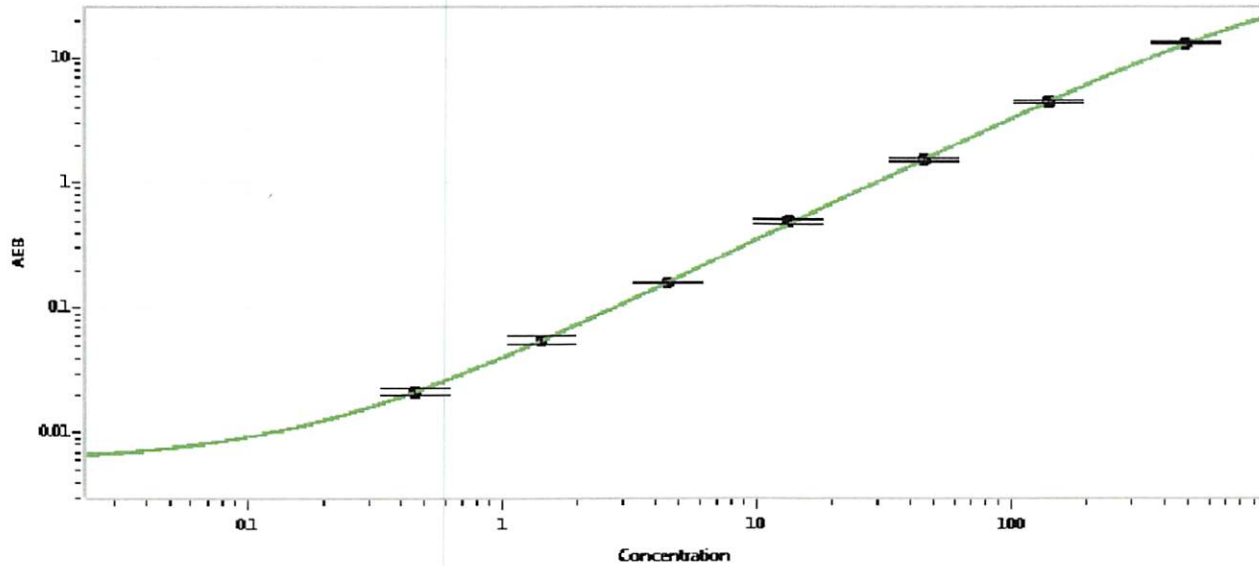
Component	Part Number	Lot Number
Bead Reagent	102246	901504
Detector Reagent	102248	901604
SBG Reagent	102250	901603
Sample Diluent	102252	901606
Reference Calibrators	102255	901401/833201
Controls	102256 - 102257	822806/830201
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	822806	4.85	3.63-5.71
Control 2	830201	172	125-187
Panel 1-A	823301	5.60	4.08-6.68
Panel 1-B	823302	7.08	4.90-7.61
Panel 2	823601	88.3	59.3-101
Panel 3	823602	350	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	Associate Director, QC	<i>P. Stadler</i> 26 Jan 2019
	Name	Title	Signature/Date