

Product Number:	103186
Lot Number:	501642
Expiration:	15-Jan-2020
Platform(s):	HD-1 & SR-X



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

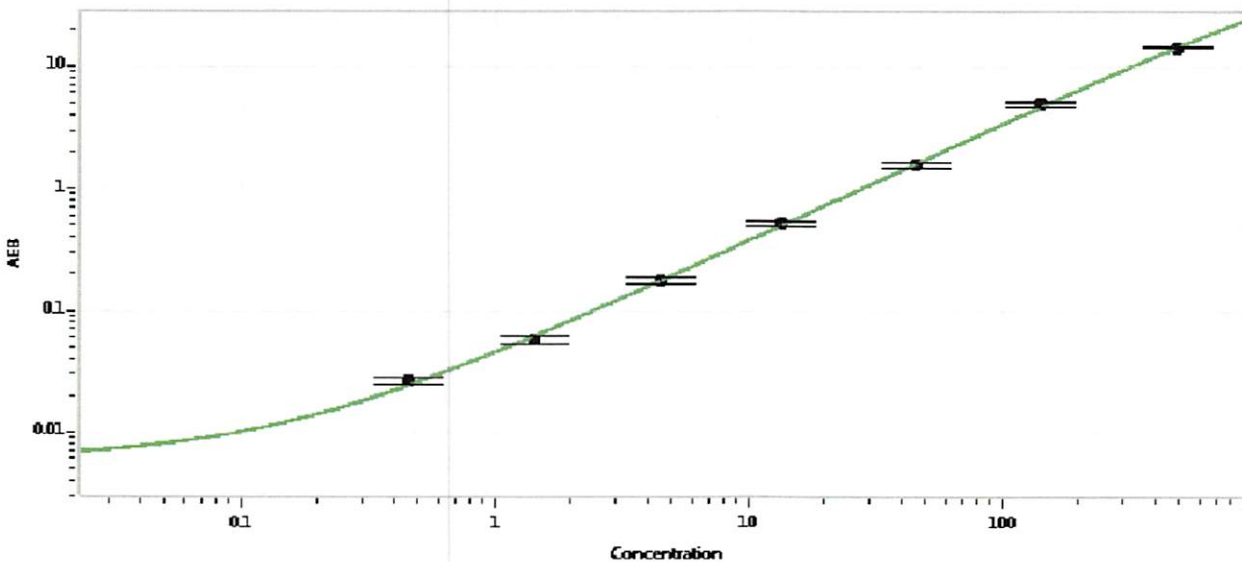
Component	Part Number	Lot Number
Bead Reagent	102246	901504
Detector Reagent	102248	901604
SBG Reagent	102250	901603
Sample Diluent	102252	901606
Reference Calibrators	102255	903501
Controls	102256 - 102257	903502
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	903502	4.76	3.85-5.77
Control 2	903502	209	157-236
Panel 1-A	823301	5.23	4.08-6.68
Panel 1-B	823302	6.29	4.90-7.61
Panel 2	823601	91.4	59.3-101
Panel 3	823602	351	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	0.568
C	1.50
D	5.05
E	14.1
F	48.1
G	147
H	487

Review/Approval	Patrizia Stadler	QC Assoc. Director	<i>P. Stadler</i>	07 Feb 2019
	Name	Title	Signature/Date	