

Product Number:	102215
Lot Number:	501671
Expiration:	8-Feb-2020
Platform(s):	HD-1 & SR-X



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

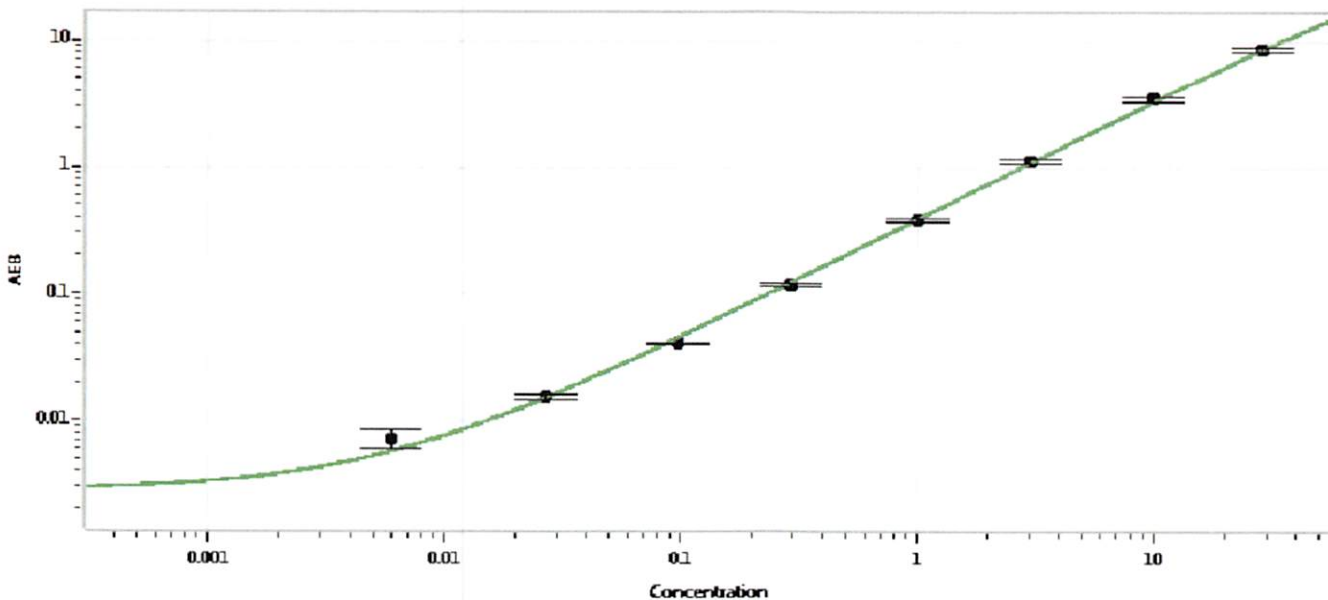
Component	Part Number	Lot Number
Bead Reagent	102211	903905
Detector Reagent	102212	904206
SBG Reagent	102213	903907
Sample Diluent	102214	903906
Calibrators	101310	904501
Controls	100702-100703	904502
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	904502	0.089	0.073-0.111
Control 2	904502	10.7	8.03-12.0
Panel 1	MF-092217-01	0.259	0.177-0.353
Panel 2	MF-092217-02	2.71	1.99-3.51
Panel 3	MF-101117-01	17.8	12.8-20.8

* 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	0.006
C	0.027
D	0.097
E	0.293
F	0.993
G	2.99
H	9.95
I	28.7

Review/Approval

Patrizia Stadler

Assoc. Director, QC

P. Stadler 25 Feb 2019

Name

Title

Signature/Date