

Product Number:	101478
Lot Number:	501975
Expiration:	19-Jun-2020
Platform(s):	HD-1/HD-x & SR-X



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

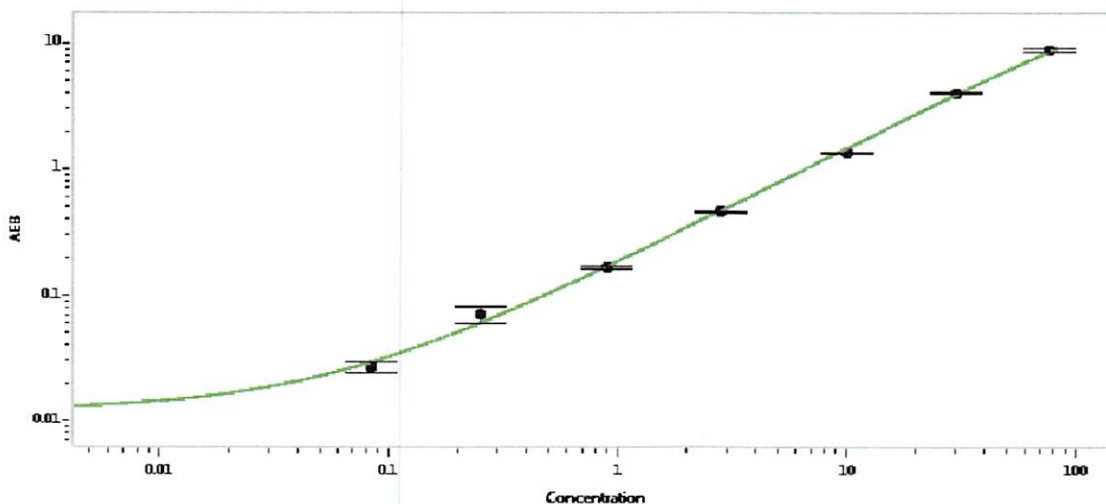
Component	Part Number	Lot Number
Bead Reagent	101480	919110
Detector Reagent	101481	917102
SBG Reagent	101482	917005
Sample Diluent	101483	917104
Calibrators	100758-100765	921201
Controls	100768 & 100769	921202
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	921202	2.45	2.16-3.24
Control 2	921202	61.6	51.2-76.8
Panel 1a	MF-051017-04	0.557	0.508-0.855
Panel 2	MF-042117-06	4.18	3.89-6.07
Panel 3	MF-042117-07	59.8	45.5-73.7

* 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	0.084
C	0.252
D	0.899
E	2.82
F	9.94
G	30.0
H	76.3

Review/Approval

Patrizia Stadler
Name

Assoc. Director, QC
Title

P. Stadler
Signature

09 Aug 2019
Date