

Product Number:	101478
Lot Number:	502267
Expiration:	29-Jan-2021
Platform(s):	HD-1/HD-X & SR-X



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

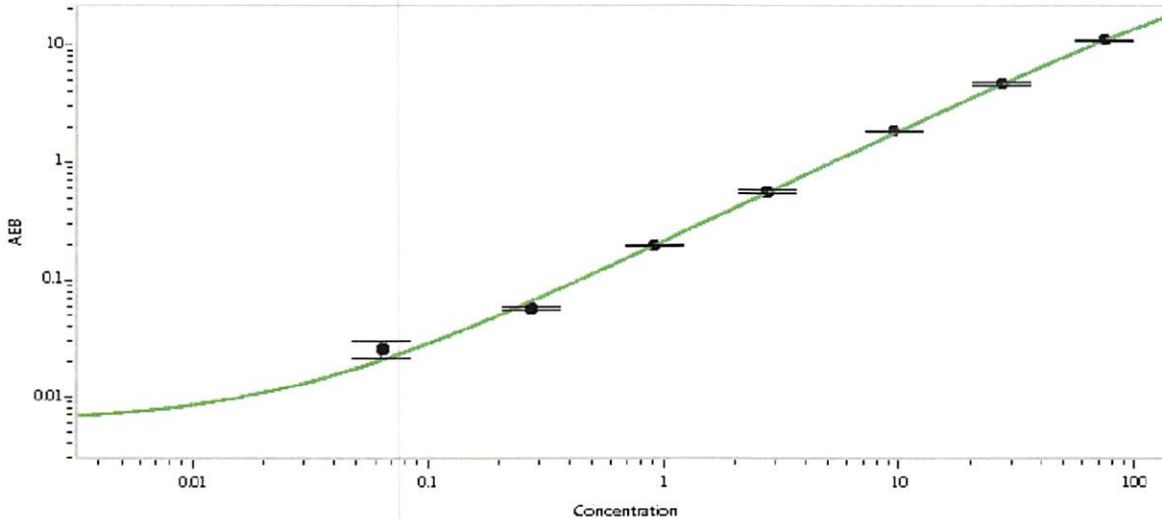
Component	Part Number	Lot Number
Bead Reagent	101480	006312
Detector Reagent	101481	006313
SBG Reagent	101482	006314
Sample Diluent	101483	006315
Calibrators	100758-100765	928901
Controls	100768 & 100769	928902
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	928902	2.61	2.16-3.24
Control 2	928902	60.2	47.5-71.3
Panel 1	MF-051017-04	0.577	0.508-0.855
Panel 2	MF-042117-06	4.04	3.89-6.07
Panel 3	MF-042117-07	47.7	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	0.064
C	0.274
D	0.906
E	2.77
F	9.42
G	27.3
H	74.4

Review/Approval

Patrizia Stadler

Assoc. Director, QC

P. Stadler

11 MAR 2020

Name

Title

Signature/Date