

Product Number:	101478
Lot Number:	501909
Expiration:	19-Jun-2020
Platform(s):	HD-1/HD-x & SR-X



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

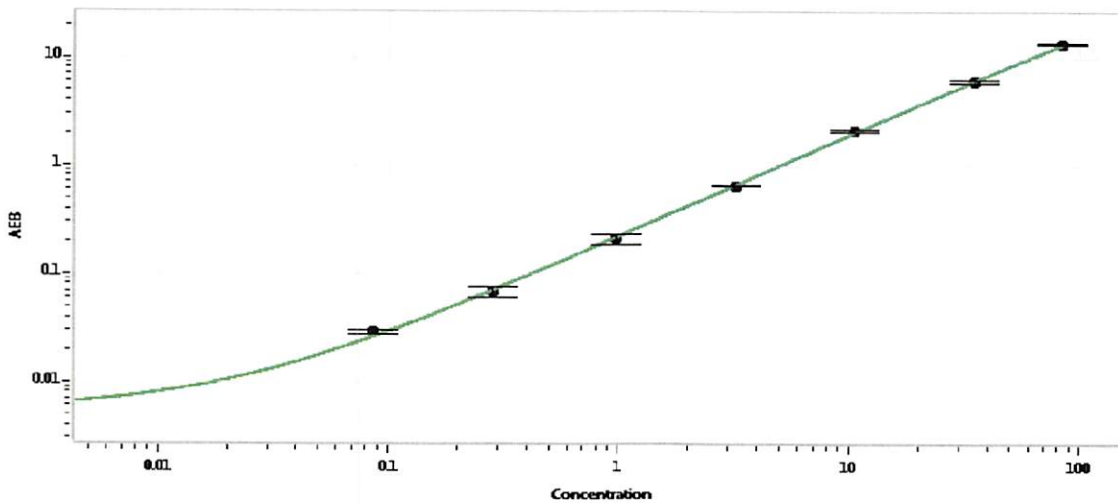
Component	Part Number	Lot Number
Bead Reagent	101480	919110
Detector Reagent	101481	917102
SBG Reagent	101482	917005
Sample Diluent	101483	917104
Calibrators	100758-100765	911301 / 912101
Controls	100768 & 100769	912102
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	912102	3.11	2.34-3.60
Control 2	912102	72.6	54.9-82.3
Panel 1a	MF-051017-04	0.670	0.508-0.855
Panel 2	MF-042117-06	4.88	3.89-6.07
Panel 3	MF-042117-07	60.1	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	0.086
C	0.288
D	0.980
E	3.23
F	10.6
G	34.6
H	82.9

Review/Approval: Patrizia Stadler Name, Assoc. Director, QC Title, P Stadler Signature, 17 Jul 2019 Date