

Product Number:	101478
Lot Number:	501647
Expiration:	19-Oct-2019
Platform(s):	HD-1 & SR-X



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

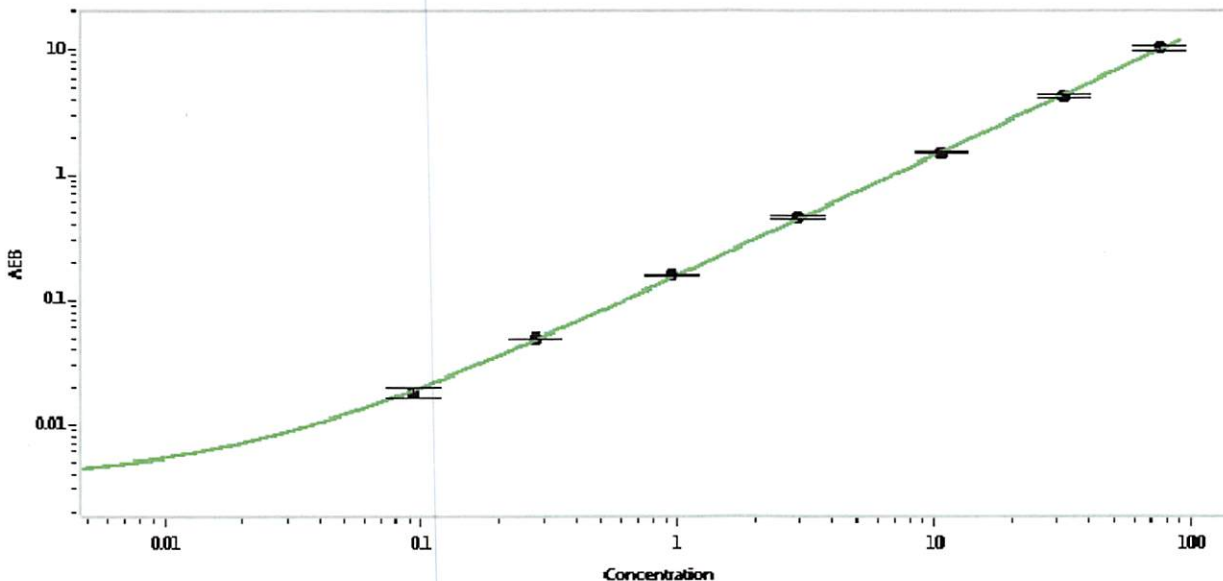
Component	Part Number	Lot Number
Bead Reagent	101480	831217
Detector Reagent	101481	829812
SBG Reagent	101482	829211
Sample Diluent	101483	829609
Calibrators	100758-100765	833801
Controls	100768 & 100769	833802
RGP Reagent	103159	N/A <sup>1</sup>

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

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	833802	2.94	2.28-3.42
Control 2	833802	63.9	49.8-74.6
Panel 1	MF-051017-04	0.638	0.508-0.855
Panel 2	MF-042117-06	4.33	3.89-6.07
Panel 3	MF-042117-07	56.3	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	0.093
C	0.278
D	0.948
E	2.95
F	10.6
G	32.1
H	75.1

Review/Approval: Patrizia Stadler QC Assoc. Director P. Stadler 08 Feb 2019

Name Title Signature/Date