

Product Number:	101478
Lot Number:	502073
Expiration:	11-Oct-2020
Platform(s):	HD-1/HD-X & SR-X



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

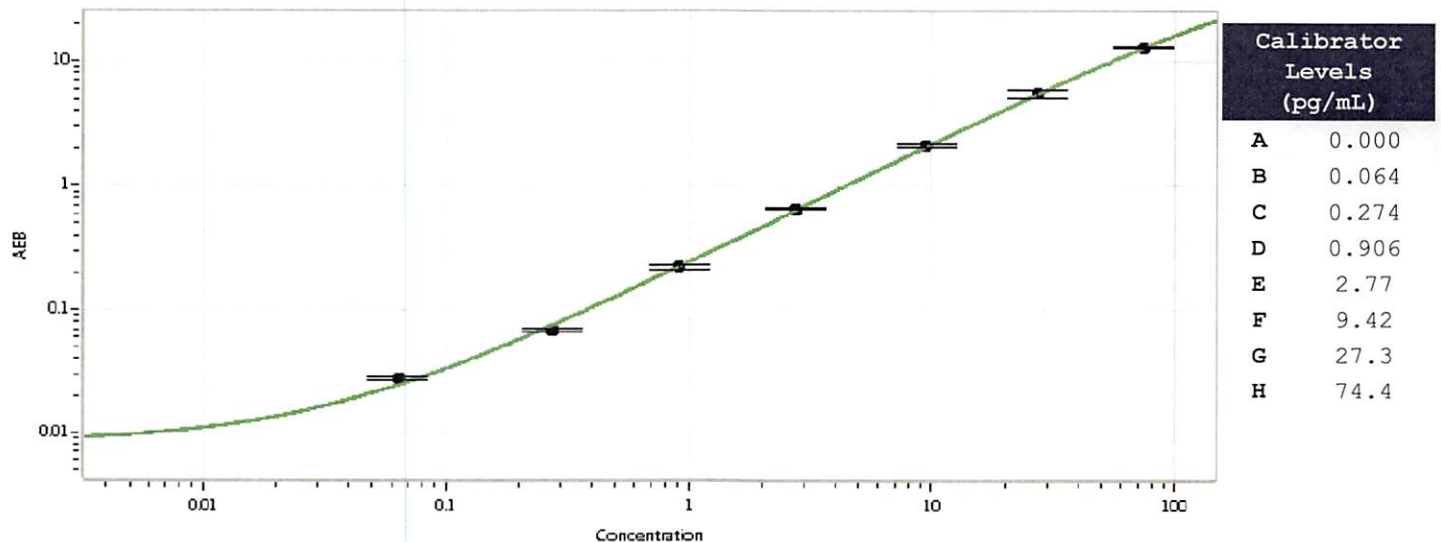
Component	Part Number	Lot Number
Bead Reagent	101480	928408
Detector Reagent	101481	928411
SBG Reagent	101482	928410
Sample Diluent	101483	928702
Calibrators	100758-100765	928901
Controls	100768 & 100769	928902
RGP Reagent	103159	N/A <sup>1</sup>

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

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	928902	2.73	2.18-3.27
Control 2	928902	57.0	47.4-71.1
Panel 1	MF-051017-04	0.547	0.508-0.855
Panel 2	MF-042117-06	4.09	3.89-6.07
Panel 3	MF-042117-07	51.6	45.5-73.7

\* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibration Curve



Review/Approval: Patrizia Stadler Assoc. Director, QC P. Stodler 06 NOV 2019  
Name Title Signature/Date