

Product Number:	100860
Lot Number:	502421
Expiration:	8-May-2021
Platform(s):	HD-1/ HD-X



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

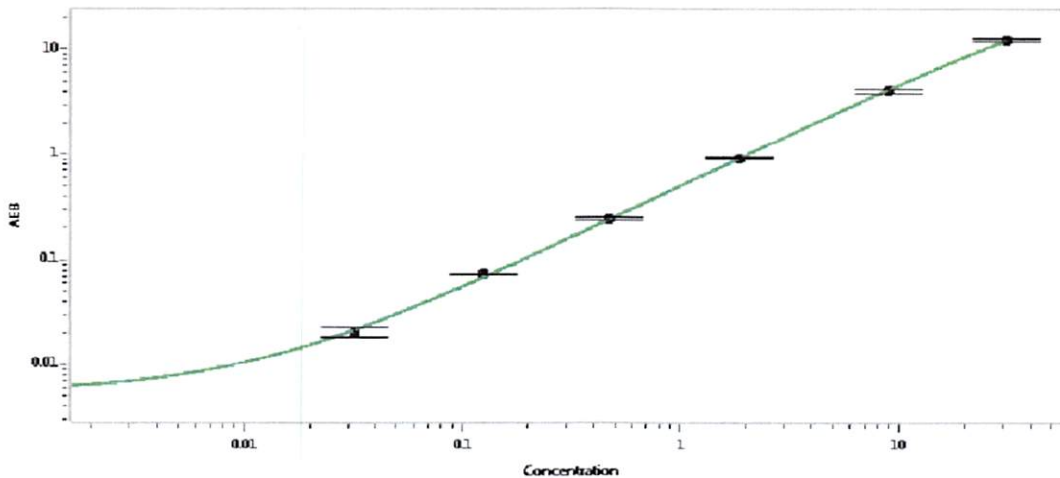
Component	Part Number	Lot Number
Bead Reagent	100865	016106
Detector Reagent	100862	016107
SBG Reagent	100868	012910
Sample Diluent	100870	013205
Calibrators	101111	016901
Controls	100875 & 100876	016902
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	016902	6.53	4.99-7.48
Control 2	016902	0.097	0.076-0.118
Panel 1	908501	1.25	0.631-2.10
Panel 2	908502	9.67	4.38-12.5
Panel 3	908503	32.4	12.1-47.2

* 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	0.032
C	0.126
D	0.470
E	1.85
F	8.88
G	30.9

Review/Approval _____

Shivani Goel

QC Manager

Shivani Goel 26 Jun 2020

Name

Title

Signature/Date