

Product Number:	100860
Lot Number:	503074
Expiration:	19-Jul-2022
Platform(s):	HD-1/ HD-X



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

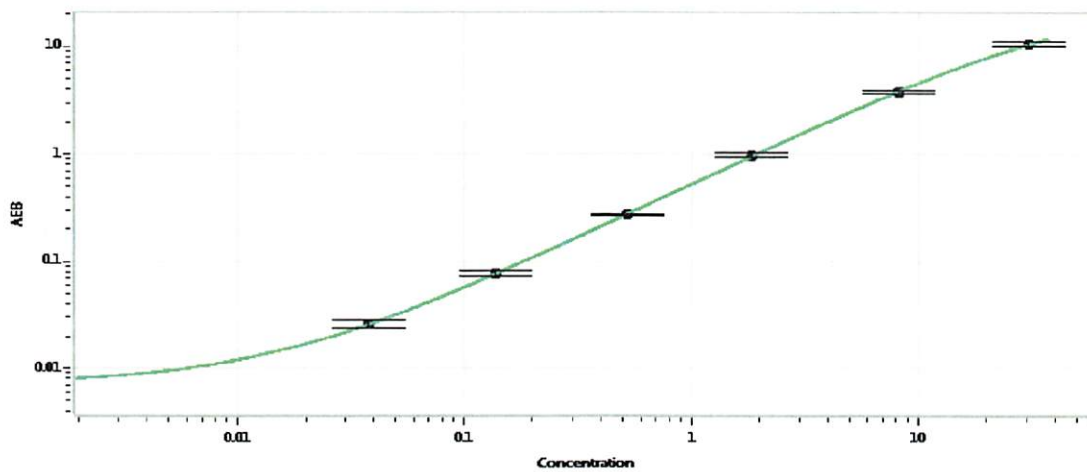
Component	Part Number	Lot Number
Bead Reagent	100865	120704
Detector Reagent	100862	120801
SBG Reagent	100868	120802
Sample Diluent	100870	120907
Calibrators	101111	120902
Control 1	100875	120903
Control 2	100876	120903
RGP Reagent	103159	N/A <sup>1</sup>

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

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	120903	5.87	4.90 - 7.35
Control 2	120903	0.106	0.055 - 0.192
Panel 1	115916	1.14	0.911 - 1.37
Panel 2	115916	4.55	3.25 - 9.53
Panel 3	115916	26.0	21.4 - 52.0

\* 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	0.038
C	0.137
D	0.521
E	1.85
F	8.17
G	30.6

Review/Approval: Shivani Goel Name: Shivani Goel QC Manager: Shivani Goel Title: QC Manager Signature/Date: 09 Mar 2022