

Product Number:	100860
Lot Number:	503166
Expiration:	3-Oct-2022
Platform(s):	HD-1 & HD-X



Data below represents results generated on the Simoa® HD-X Analyzer.

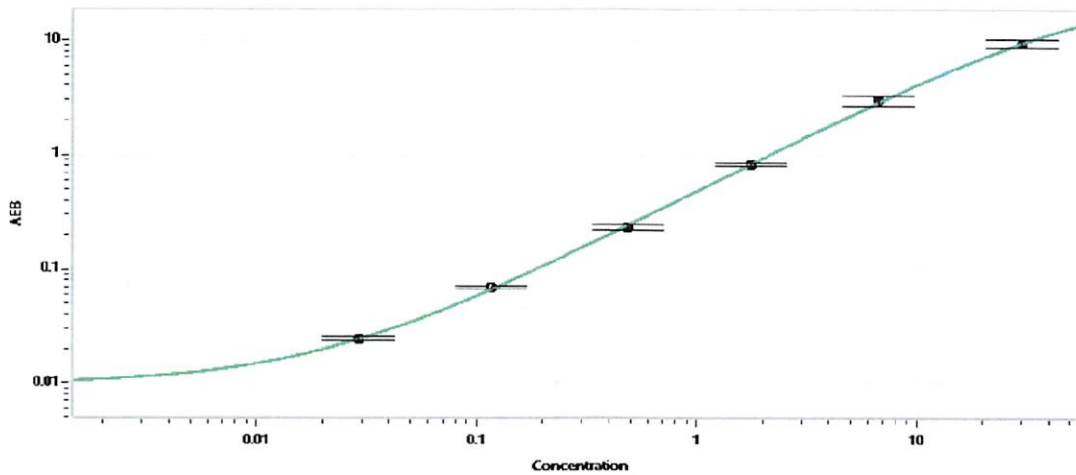
Component	Part Number	Lot Number
Bead Reagent	100865	127605
Detector Reagent	100862	127708
SBG Reagent	100868	210807
Sample Diluent	100870	127710
Calibrators	101111	130705
Control 1	100875	130706
Control 2	100876	130706
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	130706	4.63	2.99 - 7.18
Control 2	130706	0.092	0.071 - 0.126
Panel 1	115916	1.00	0.830 - 1.35
Panel 2	115916	4.11	1.78 - 9.41
Panel 3	115916	39.0	22.1 - 52.6

* 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 0.029
C 0.117
D 0.484
E 1.77
F 6.68
G 29.9

Review/Approval	<u>Matt Smart</u>	QC	<u>Matt Smart</u> 13 Jun 2022
	Name	Title	Signature/Date