

Product Number:	100860
Lot Number:	503771
Expiration:	18-Jan-2024
Platform(s):	HD-1 & HD-X



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

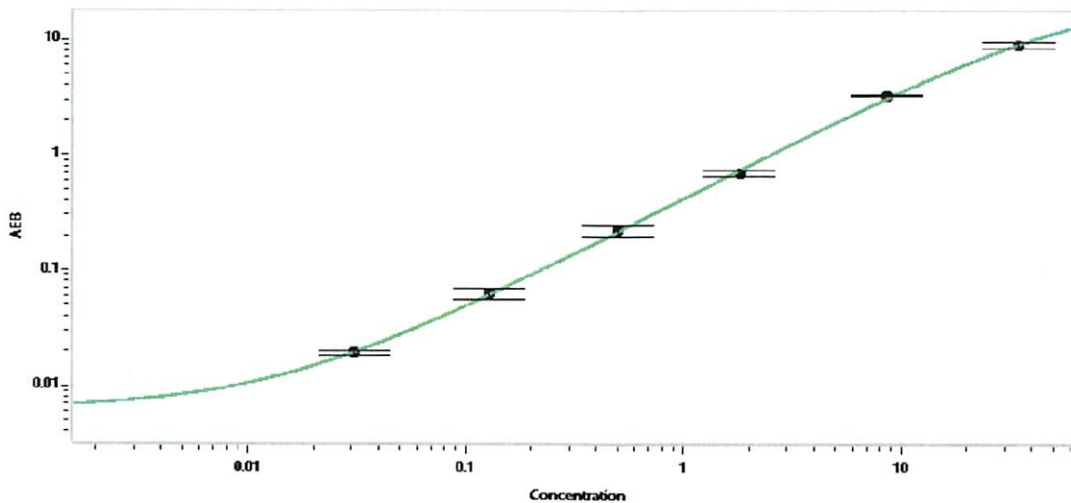
Component	Part Number	Lot Number
Bead Reagent	100865	308111
Detector Reagent	100862	303004
SBG Reagent	100868	303005
Sample Diluent	100870	301803
Calibrators	101111	307402
Control 1	100875	307403
Control 2	100876	307403
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	307403	5.86	3.52-7.51
Control 2	307403	0.116	0.081-0.129
Panel 1	306704	0.775	0.772-1.16
Panel 2	306704	4.90	2.98-7.27
Panel 3	306704	16.3	8.18-26.3

* 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	0.031
C	0.128
D	0.500
E	1.81
F	8.56
G	34.4

Review/Approval Matt Koubouris QC Supervisor [Signature] 30 Mar 2023
 Name Title Signature/Date