

Product Number:	103180
Lot Number:	503135
Expiration:	2-Dec-2021
Platform(s):	HD-1/ HD-X



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

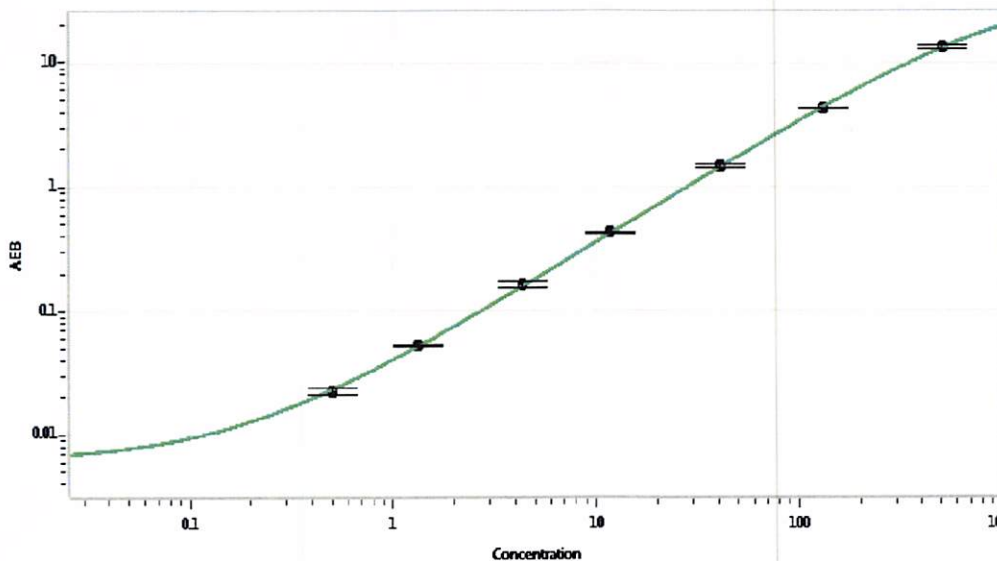
Component	Part Number	Lot Number
Bead Reagent	102246	115708
Detector Reagent	102248	115809
SBG Reagent	102250	115810
Sample Diluent	102252	115914
Calibrators	102255	933601
Control 1	102256	933901
Control 2	102257	933901
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	933901	3.77	2.60 - 4.70
Control 2	933901	151	122 - 183
Panel 1	008501	13.0	8.59 - 20.5
Panel 2	008501	10.2	7.39 - 15.1
Panel 3	008501	63.5	43.8 - 91.8
Panel 4	009401	204	128 - 302

* 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	0.501
C	1.32
D	4.32
E	11.8
F	41.1
G	131
H	506

Review/Approval Junpil Park QC Supervisor [Signature] 20 Sep 2021
Name Title Signature/Date