

Product Number:	103186
Lot Number:	503546
Expiration:	3-Aug-2023
Platform(s):	HD-1 / HD-X



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

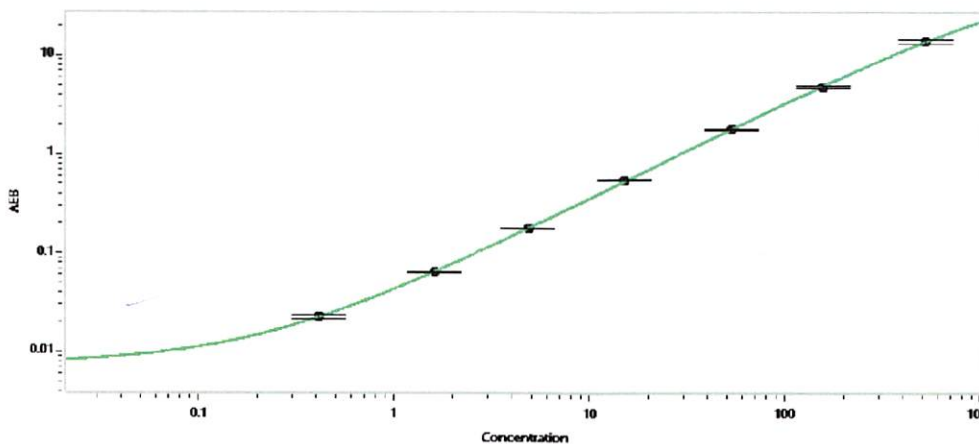
Component	Part Number	Lot Number
Bead Reagent	102246	222101
Detector Reagent	102248	222107
SBG Reagent	102250	222108
Sample Diluent	102252	222201
Calibrators	102255	227201
Control 1	102256	227202
Control 2	102257	227202
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	227202	9.63	7.61-11.5
Control 2	227202	383	317-476
Panel 1	224402	9.94	7.63-12.7
Panel 2	224402	7.71	5.89-9.40
Panel 3	224402	72.4	60.5-92.6
Panel 4	212511	255	150-356

* 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	0.415
C	1.60
D	4.83
E	15.1
F	52.6
G	155
H	517

Review/Approval

Matt Koubouris
Name

QC Supervisor
Title

[Signature] 20sep2022
Signature/Date