

Product Number:	101552
Lot Number:	503536
Expiration:	6-Jul-2023
Platform(s):	HD-1/HD-x & SR-X



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

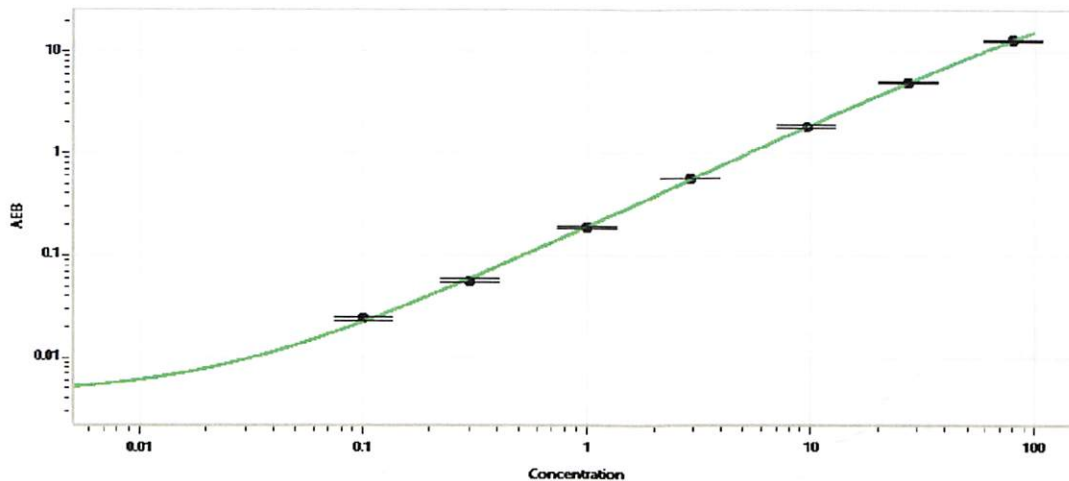
Component	Part Number	Lot Number
Bead Reagent	101554	218701
Detector Reagent	101553	218806
SBG Reagent	101555	218817
Sample Diluent	101556	218812
Calibrators	101559	218904
Control 1	101560	218905
Control 2	101561	218905
SBG Diluent (SR-X) †	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	218905	0.971	0.803-1.22
Control 2	218905	52.3	45.5-68.4
Panel 1	117205	3.64	3.49-5.23
Panel 2	117205	2.49	1.89-4.80
Panel 3	117205	20.0	17.8-26.7
Panel 4	119612	29.6	23.3-41.8

* 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	0.101
C	0.301
D	1.00
E	2.88
F	9.60
G	27.3
H	79.5

Review/Approval

Matt Koulouris

Name

QC Supervisor

Title

[Signature] 22 Nov 2022

Signature/Date