

Product Number:	101552
Lot Number:	503311
Expiration:	31-Jan-2023
Platform(s):	HD-1/HD-x & SR-X



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

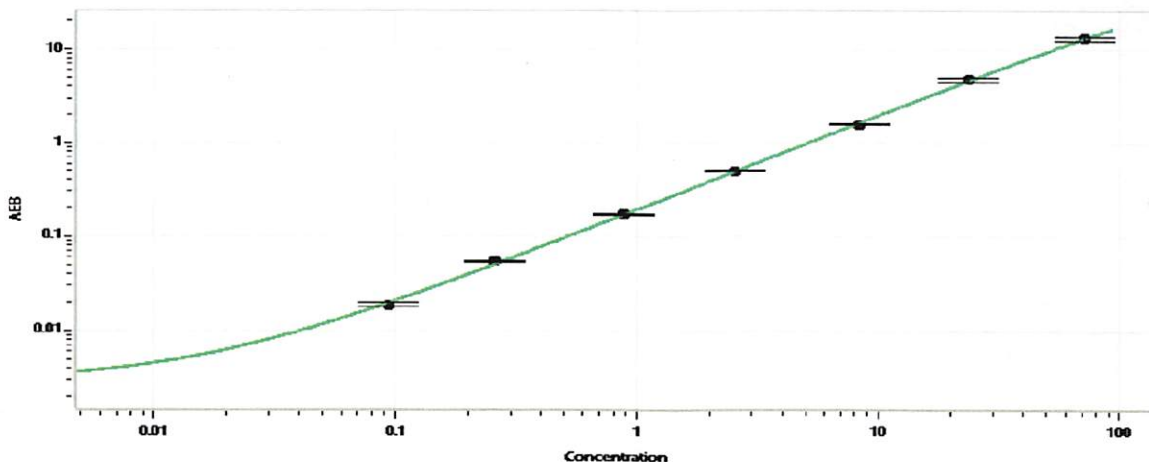
Component	Part Number	Lot Number
Bead Reagent	101554	203205
Detector Reagent	101553	203110
SBG Reagent	101555	203206
Sample Diluent	101556	203209
Calibrators	101559	203804
Control 1	101560	203805
Control 2	101561	203805
SBG Diluent (SR-X) †	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	203805	0.917	0.682-1.15
Control 2	203805	50.2	41.6-62.3
Panel 1	117205	5.01	3.49-5.23
Panel 2	117205	3.03	1.89-4.80
Panel 3	117205	22.8	17.8-26.7
Panel 4	119612	35.3	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	0.094
C	0.257
D	0.880
E	2.53
F	8.33
G	23.8
H	71.0

Review/Approval: Sonia Khillan Name: Sonia Khillan Title: QC Signature/Date: Sonia Khillan / 08NOV2022