

Product Number:	104073
Lot Number:	503501
Expiration:	7-Jul-2023
Platform(s):	HD-1 / HD-X / SR-X



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

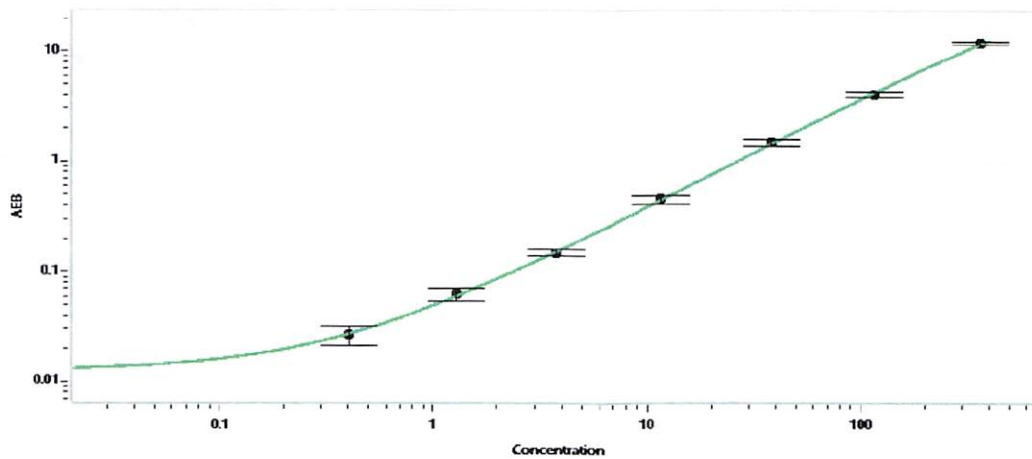
Component	Part Number	Lot Number
Bead Reagent	104067	221610
Detector Reagent	104068	218803
SBG Reagent	104069	218804
Sample Diluent	104071	218805
Calibrators	102255	214302
Control 1	102256	214303
Control 2	102257	214303
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	214303	4.34	3.16-4.75
Control 2	214303	156	126-190
Panel 1	224402	10.3	7.80-12.5
Panel 2	224402	7.01	5.71-9.38
Panel 3	224402	72.6	60.3-91.5
Panel 4	212511	326	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	0.403
C	1.28
D	3.80
E	11.6
F	38.1
G	115
H	365

Review/Approval Sonia Khillan QC Manager Sonia Khillan / 18 Oct 2022  
 Name Title Signature/Date