

Product Number:	104073
Lot Number:	503847
Expiration:	3-Apr-2024
Platform(s):	HD-1 / HD-X / SR-X



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

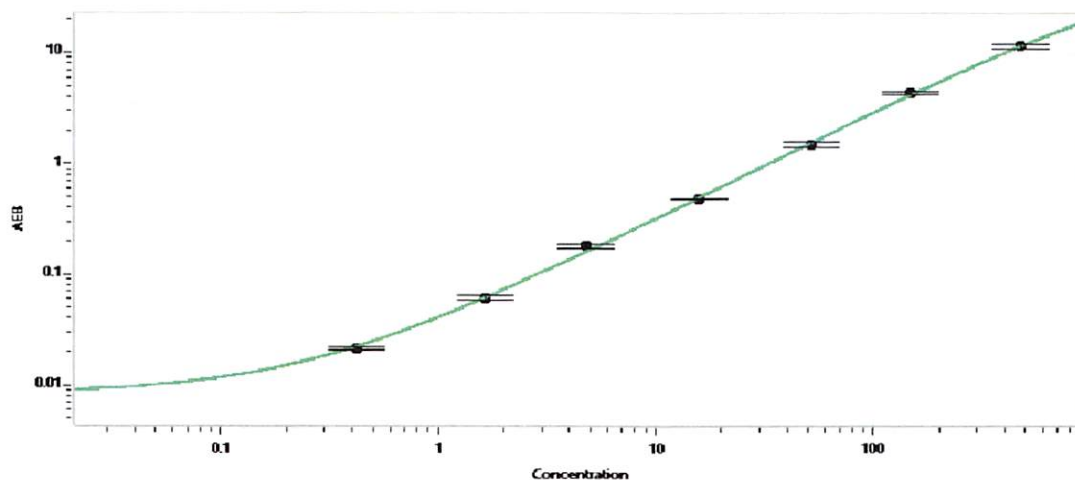
Component	Part Number	Lot Number
Bead Reagent	104067	309305
Detector Reagent	104068	309404
SBG Reagent	104069	309405
Sample Diluent	104071	309406
Calibrators	102255	309302
Control 1	102256	309303
Control 2	102257	309303
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	309303	5.44	3.76 - 6.65
Control 2	309303	248	185 - 277
Panel 1	224402	10.6	8.84 - 13.3
Panel 2	224402	7.08	6.29 - 9.65
Panel 3	224402	79.0	63.0 - 94.5
Panel 4	301701	626	519 - 808

* 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	0.421
C	1.65
D	4.76
E	15.7
F	51.2
G	148
H	474

Review/Approval

Sonia Khillan

Name

QC

Title

Sonia Kella 10MAY2023

Signature/Date