

Product Number:	104292
Lot Number:	503887
Expiration:	22-Feb-2024
Platform(s):	HD-1 / HD-X



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

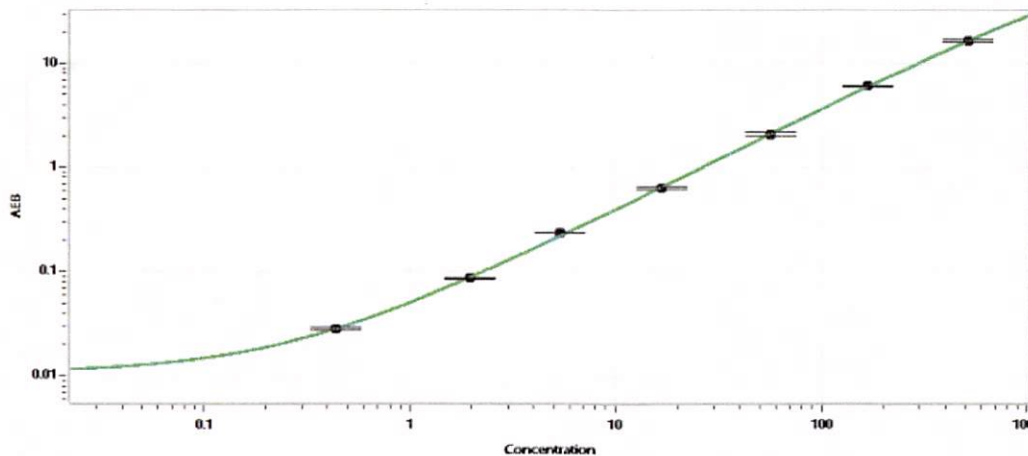
Component	Part Number	Lot Number
Bead Reagent	104067	307612
Detector Reagent	104068	308008
SBG Reagent	104069	308002
Sample Diluent	104071	308003
Calibrators	102255	308204
Control 1	102256	308205
Control 2	102257	308205
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	308205	5.59	4.75 - 7.12
Control 2	308205	263	208 - 312
Panel 1	224402	9.99	8.84 - 13.3
Panel 2	224402	7.24	6.29 - 9.65
Panel 3	224402	74.5	63.0 - 94.5
Panel 4	212511	278	175 - 376

* 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 0.440
C 1.95
D 5.35
E 16.6
F 57.0
G 169
H 518

Review/Approval Sonia Khillan AC Manager Sonia Khillan 24 MAY 2023
 Name Title Signature/Date