

Product Number:	104073
Lot Number:	503957
Expiration:	5-Aug-2024
Platform(s):	HD-1 / HD-X / SR-X



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

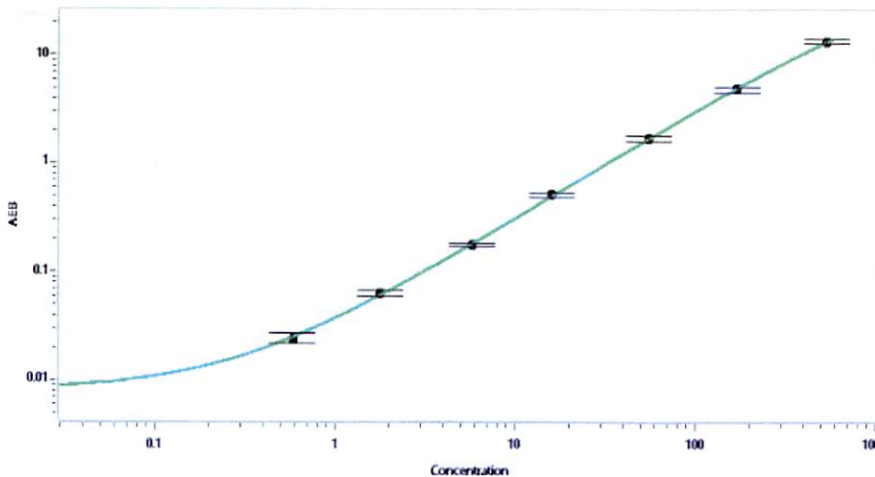
Component	Part Number	Lot Number
Bead Reagent	104067	327004
Detector Reagent	104068	327203
SBG Reagent	104069	327106
Sample Diluent	104071	326505
Calibrators	102255	327204
Control 1	102256	327205
Control 2	102257	327205
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	327205	15.1	12.1-18.1
Control 2	327205	631	483-725
Panel 1	224402	9.89	7.86-11.8
Panel 2	224402	6.84	5.60-8.59
Panel 3	224402	68.3	56.1-84.1
Panel 4	301701	544	462-719

* 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	0.585
C	1.78
D	5.84
E	16.0
F	55.3
G	171
H	540

Review/Approval Sonia Khillan QC Sonia Khillan 19 OCT 2023
 Name Title Signature/Date