

Product Number:	104073
Lot Number:	504012
Expiration:	22-Sep-2024
Platform(s):	HD-1 / HD-X / SR-X



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

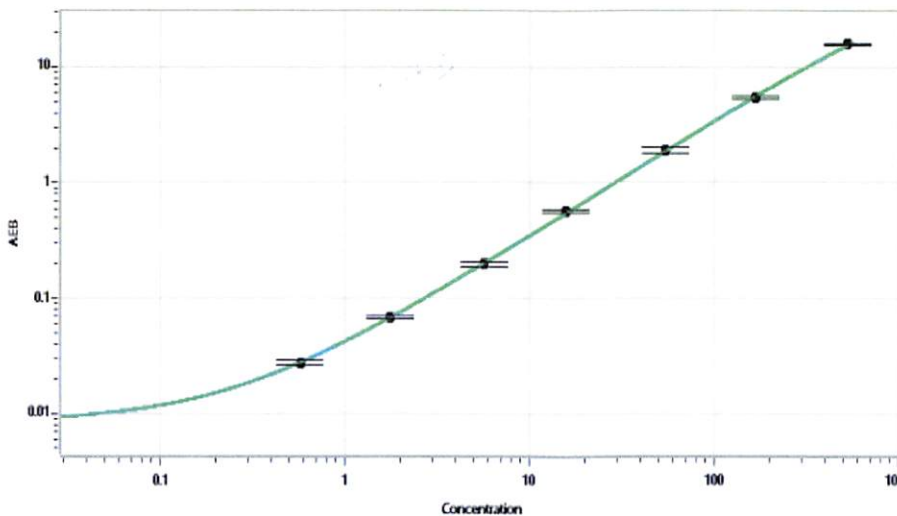
Component	Part Number	Lot Number
Bead Reagent	104067	328508
Detector Reagent	104068	327203
SBG Reagent	104069	330008
Sample Diluent	104071	326505
Calibrators	102255	327204
Control 1	102256	327205
Control 2	102257	327205
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	c. Range (pg/mL)
Control 1	327205	14.5	11.8 - 17.7
Control 2	327205	607	499 - 749
Panel 1	224402	9.57	7.86 - 11.8
Panel 2	224402	6.66	5.60 - 8.59
Panel 3	224402	68.2	56.1 - 84.1
Panel 4	301701	521	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	0.574
C	1.75
D	5.73
E	15.7
F	54.3
G	168
H	530

Review/Approval Sonia Khillan QC Sonia Khillan 09 NOV 2023
 Name Title Signature/Date