

Product Number:	104073
Lot Number:	503892
Expiration:	30-Aug-2024
Platform(s):	HD-1 / HD-X / SR-X



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

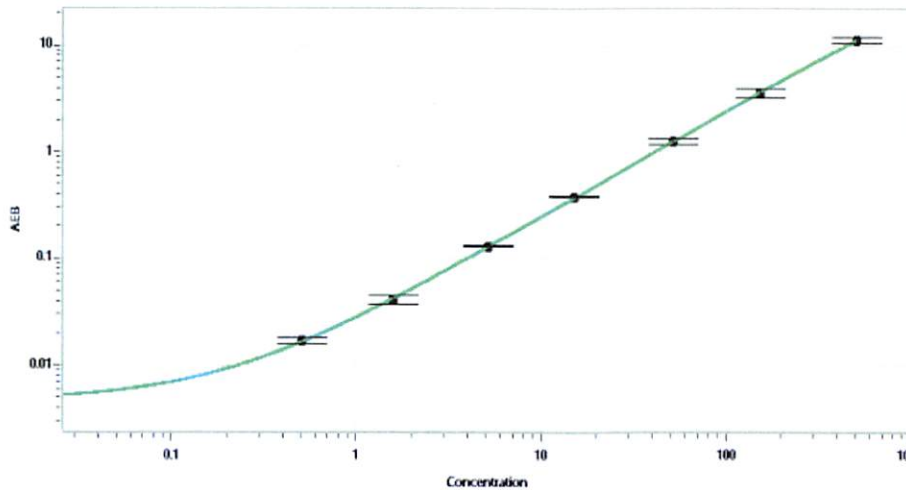
Component	Part Number	Lot Number
Bead Reagent	104067	324202
Detector Reagent	104068	324203
SBG Reagent	104069	324301
Sample Diluent	104071	324208
Calibrators	102255	324204
Control 1	102256	324205
Control 2	102257	324205
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Range (pg/mL)
Control 1	324205	13.5	11.0-16.5
Control 2	324205	572	472-731
Panel 1	224402	10.9	7.86-11.8
Panel 2	224402	7.90	5.60-8.59
Panel 3	224402	77.3	56.1-84.1
Panel 4	301701	594	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	0.514
C	1.58
D	5.21
E	15.1
F	51.9
G	153
H	506

Review/Approval ROBERT SOMOGYI QC, MANAGER [Signature] 27SEP2023
 Name Title Signature/Date