

Product Number:	104073
Lot Number:	503725
Expiration:	20-Apr-2023
Platform(s):	HD-1 / HD-X / SR-X



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

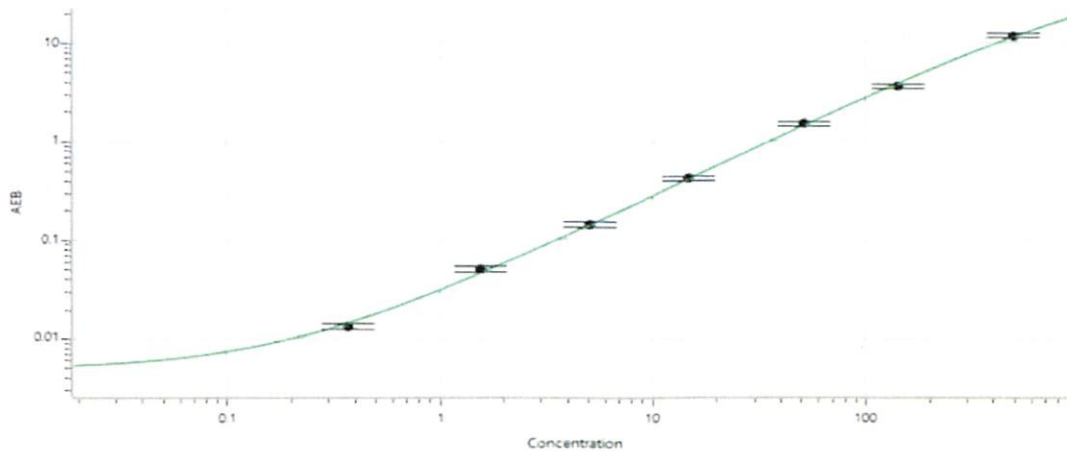
Component	Part Number	Lot Number
Bead Reagent	104067	213206
Detector Reagent	104068	212102
SBG Reagent	104069	212104
Sample Diluent	104071	212201
Calibrators	102255	227008
Control 1	102256	227009
Control 2	102257	227009
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	227009	9.67	7.98 - 12.0
Control 2	227009	369	295 - 442
Panel 1	224402	11.1	8.82 - 13.2
Panel 2	224402	8.56	6.19 - 9.75
Panel 3	224402	75.8	62.9 - 94.3
Panel 4	212511	243	186 - 366

* 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	0.373
C	1.56
D	5.06
E	14.8
F	51.1
G	142
H	494

Review/Approval Sonia Khillan QC Soni Khillan / 21 Nov 2022
 Name Title Signature/Date