

Product Number:	101478
Lot Number:	503453
Expiration:	30-Apr-2023
Platform(s):	HD-1/HD-X & SR-X



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

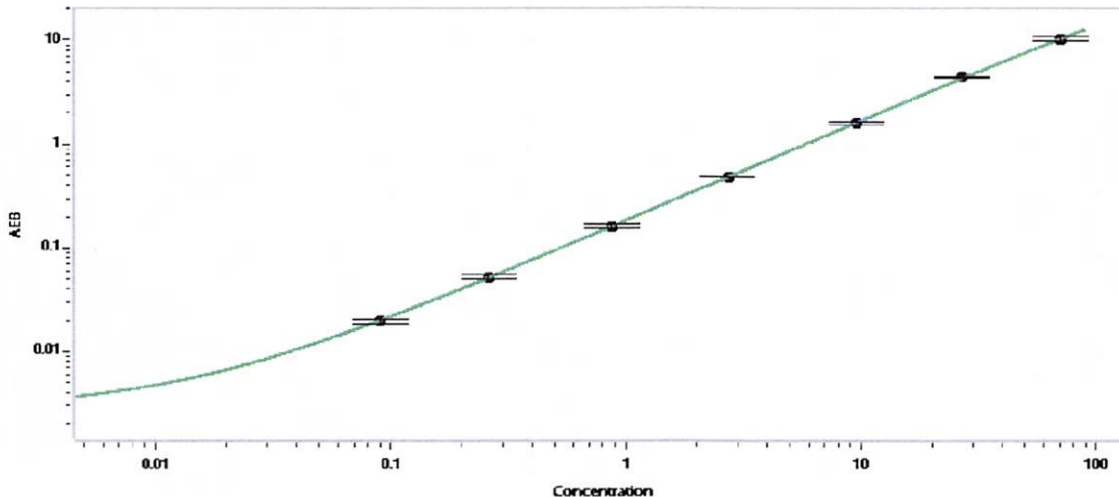
Component	Part Number	Lot Number
Bead Reagent	101480	212005
Detector Reagent	101481	212301
SBG Reagent	101482	212302
Sample Diluent	101483	212303
Calibrators	100758-100765	213502
Control 1	100768	213503
Control 2	100769	213503
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	213503	2.51	2.09-3.14
Control 2	213503	54.4	45.9-68.9
Panel 1	213804	0.828	0.441-1.26
Panel 2	122912	4.74	3.74-5.80
Panel 3	122912	46.1	37.9-69.4

* 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	0.090
C	0.260
D	0.878
E	2.71
F	9.56
G	26.9
H	70.0

Review/Approval Sonia Khillan QC Manager Sonia Khillan / 25 JUL 2023
 Name Title Signature/Date