

Product Number:	102215
Lot Number:	503952
Expiration:	25-Sep-2024
Platform(s):	HD-1/HD-X & SR-X



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

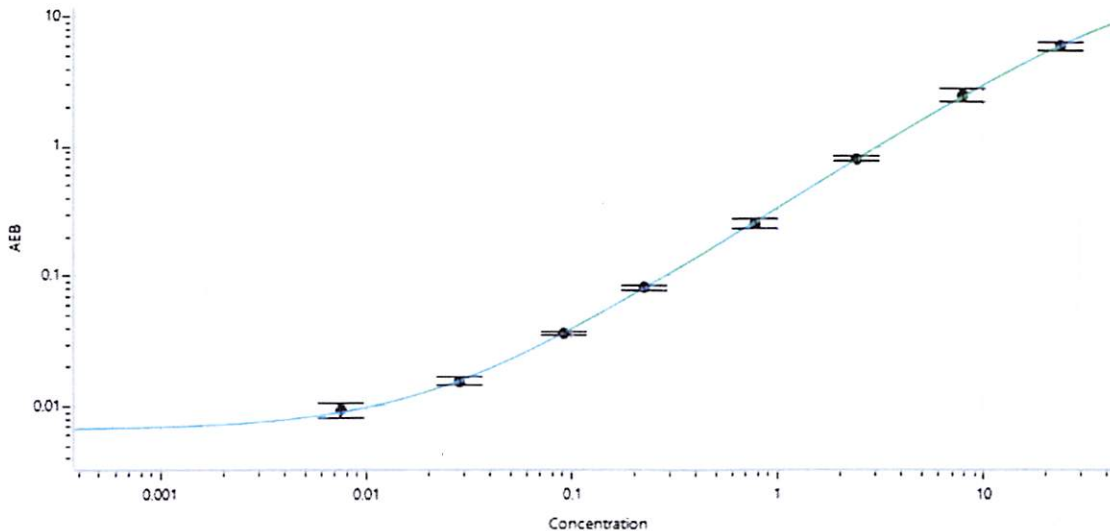
Component	Part Number	Lot Number
Bead Reagent	104297	326803
Detector Reagent	104338	327709
SBG Reagent	104340	326915
Sample Diluent	104342	326916
Calibrators	104343	326801
Control 1	104344	326802
Control 2	104345	326802
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	326802	0.089	0.070 - 0.112
Control 2	326802	7.72	6.15 - 10.3
Panel 1	019601	0.178	0.127 - 0.243
Panel 2	019601	1.79	1.56 - 2.34
Panel 3	019601	11.6	9.44 - 14.2

* 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	0.008
C	0.028
D	0.091
E	0.223
F	0.773
G	2.42
H	7.86
I	23.5

Review/Approval

Sonia Khillan

QC Manager

Sonia Khillan 16 Oct 2023

Name

Title

Signature/Date