

Product Number:	104073
Lot Number:	504172
Expiration Date:	7-Aug-24
Platform(s):	SR-X and HD-1/HD-X



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

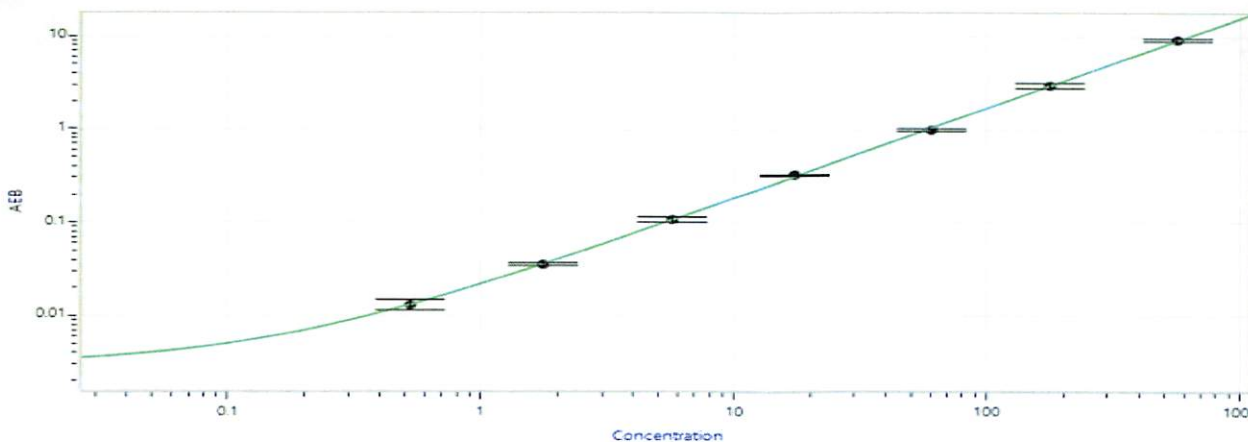
Component	Part Number	Lot Number	Control Lot Number
Bead Reagent	104067	321910	321910
Detector Reagent	104068	322002	322002
SBG Reagent	104069	322001	322001A
Sample Diluent	104071	322606	322606
Calibrators	102255	323302	323302
Control 1	102256	323303	323303
Control 2	102257	323303	323303
RGP Reagent	103159	N/A ¹	N/A ¹

¹ Reagents are not Kit Lot Specific

Release Materials* NF-Light®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	323303	13.0	10.8 - 17.1
Control 2	323303	632	514 - 771
Panel 1	224402	9.53	7.86 - 11.8
Panel 2	224402	6.23	5.60 - 8.59
Panel 3	224402	72.9	56.1 - 84.1
Panel 4	301701	582	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	0.525
C	1.75
D	5.72
E	17.3
F	59.8
G	176
H	564

Review/Approval

Cody Carlson

QC Manager

Name

Title

Cody Carlson 08 MAR 2024
Signature/Date