

Product Number:	104073
Lot Number:	504044
Expiration Date:	17-Jul-2025
Platform(s):	HD-X



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

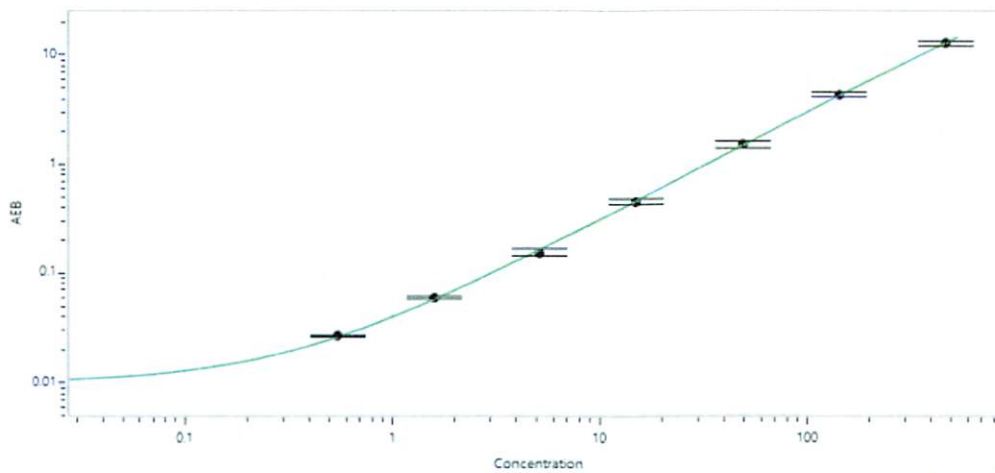
Component	Part Number	Lot Number
Bead Reagent	104067	401705
Detector Reagent	104068	401806
SBG Reagent	104069	401807
Sample Diluent	104071	401808
Calibrators	102255	402205
Control 1	102256	402206
Control 2	102257	402206
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> Reagents are not Kit Lot Specific

Release Materials* NF-Light®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	402206	12.2	10.4 - 16.3
Control 2	402206	482	440 - 660
Panel 1	224402	8.84	7.86 - 11.8
Panel 2	224402	6.11	5.60 - 8.59
Panel 3	224402	69.3	56.1 - 84.1
Panel 4	301701	519	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	0.544
C	1.59
D	5.12
E	15.0
F	48.9
G	143
H	469

Review/Approval

Cody Carlson

QC Manager

*Cody Carlson* 21 FEB 2024

Name

Title

Signature/Date