

Product Number:	104073
Lot Number:	504141
Expiration Date:	28-Sep-25
Platform(s):	SR-X and HD-1/HD-X



Component	Part Number	Lot Number
Bead Reagent	104067	414902
Detector Reagent	104068	414903
SBG Reagent	104069	408810
Sample Diluent Reagent	104071	414905
Reference Calibrators	102255	415905 (A-G), 419404 (H)
Control 1	102256	415906
Control 2	102257	415906
RGP Reagent	103159	N/A ¹

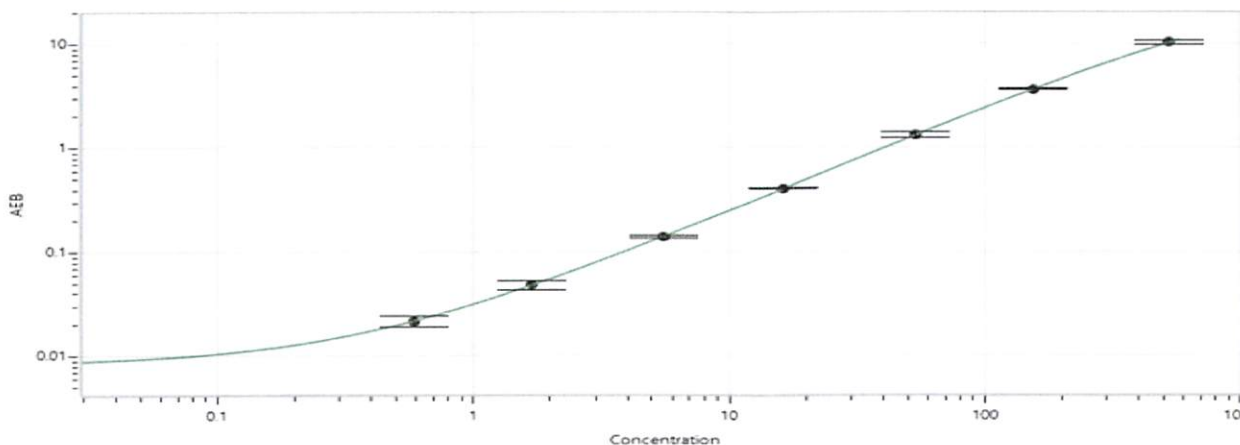
¹ Reagents are not Kit Lot Specific

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

Release Materials*			
NF-Light®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	415906	4.30	3.34 - 5.97
Control 2	415906	267	215 - 323

* 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	0.586
C	1.69
D	5.56
E	16.4
F	53.7
G	154
H	527

Review/Approval

Cody Carlson

QC Manager

Name

Title

[Signature] 05 Aug 2024
Signature/Date