

<b>Product Number:</b>	104073
<b>Lot Number:</b>	504450
<b>Expiration Date:</b>	24-Mar-26
<b>Platform(s):</b>	SR-X and HD-1/HD-X



Component	Part Number	Lot Number
Bead Reagent	104067	430914
Detector Reagent	104068	431001
SBG Reagent	104069	431002
Sample Diluent Reagent	104071	431003
Reference Calibrators	102255	430902
Control 1	102256	430903
Control 2	102257	430903
RGP Reagent	103159	N/A <sup>1</sup>

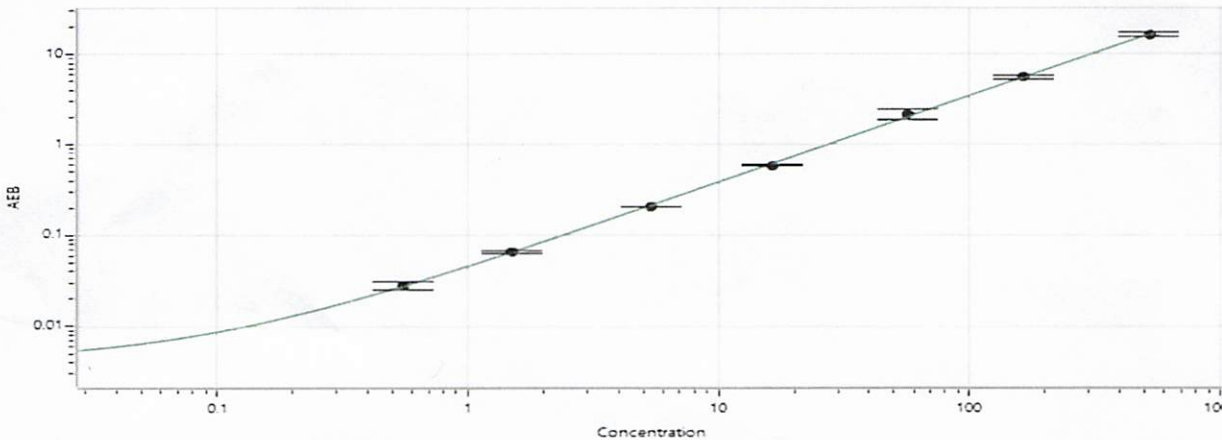
<sup>1</sup> Reagents are not Kit Lot Specific

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	430903	4.17	3.33 - 5.37
Control 2	430903	274	228 - 342

\* 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	0.554
C	1.49
D	5.34
E	16.3
F	57.0
G	164
H	520

Review/Approval

*Kyle Giroux* Senior Document Control Specialist, QA  
Name Title

*[Signature]* 03Dec2024  
Signature/Date