

Product Number:	104073
Lot Number:	504538
Expiration Date:	10-Sep-26
Platform(s):	SR-X and HD-1/HD-X

Component	Part Number	Lot Number
Bead Reagent	104067	506902
Detector Reagent	104068	507110
SBG Reagent	104069	506901
Sample Diluent Reagent	104071	507107
Reference Calibrators	102255	507305
Control 1	102256	507306
Control 2	102257	507306
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

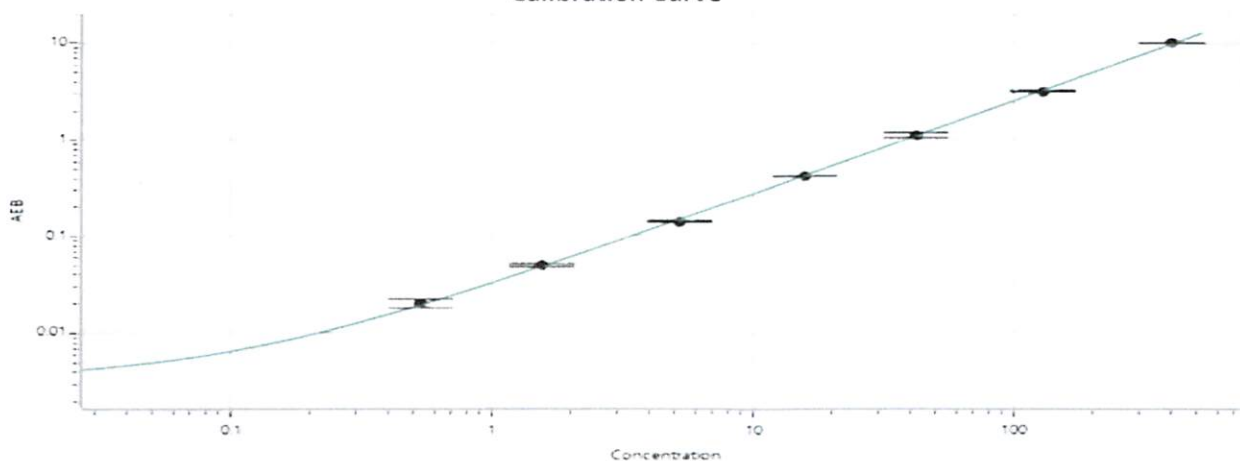


Data below represents results generated on the Simoa® HD-X Analyzer.

Release Materials*			
NF-Light®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	507306	4.33	3.62 - 5.48
Control 2	507306	218	174 - 268

* 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	0.535
C	1.56
D	5.22
E	15.7
F	42.3
G	128
H	398

Review/Approval

Dhruvi Patel

DOC Control II, QA

30 APR 2025

Name

Title

Signature/Date