

Product Number:	104073
Lot Number:	504139
Expiration Date:	13-Sep-25
Platform(s):	SR-X and HD-1/HD-X



Component	Part Number	Lot Number
Bead Reagent	104067	407342
Detector Reagent	104068	407406
SBG Reagent	104069	407407
Sample Diluent	104071	407604
Calibrators	102255	407504
Control 1	102256	407505
Control 2	102257	407505
RGP Reagent	103159	N/A ¹

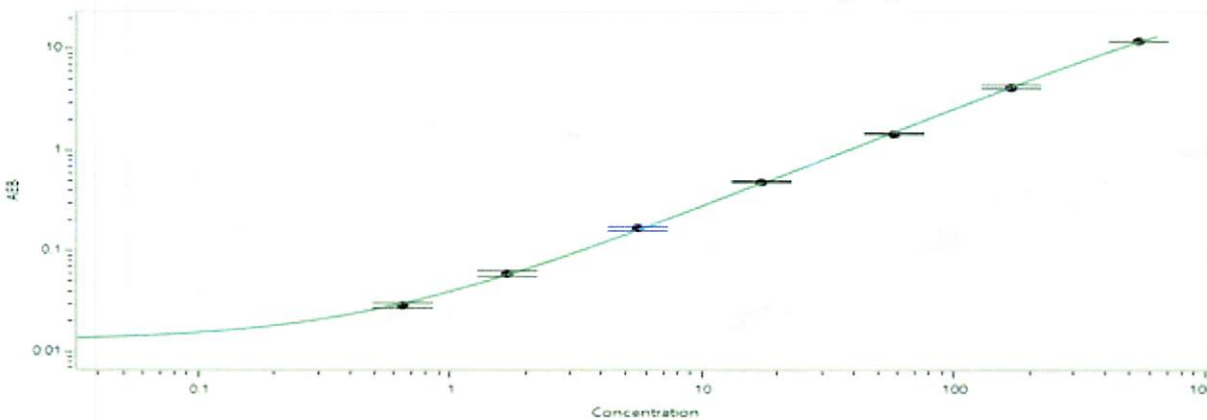
¹ Reagents are not Kit Lot Specific

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	407505	4.44	2.69 - 6.08
Control 2	407505	281	180 - 314
Panel 1	409602	10.6	6.53 - 11.4
Panel 2	409602	6.29	4.61 - 6.92
Panel 3	409602	102	57.3 - 107
Panel 4	412201	357	198 - 400

* 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	0.650
C	1.69
D	5.53
E	17.1
F	57.6
G	168
H	537

Review/Approval

ROBERT SOMOGYI QC MANAGER
Name Title

[Signature] 14 MAY 2024
Signature/Date