

Product Number:	104073
Lot Number:	504510
Expiration Date:	20-Aug-26
Platform(s):	SR-X and HD-1/HD-X



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

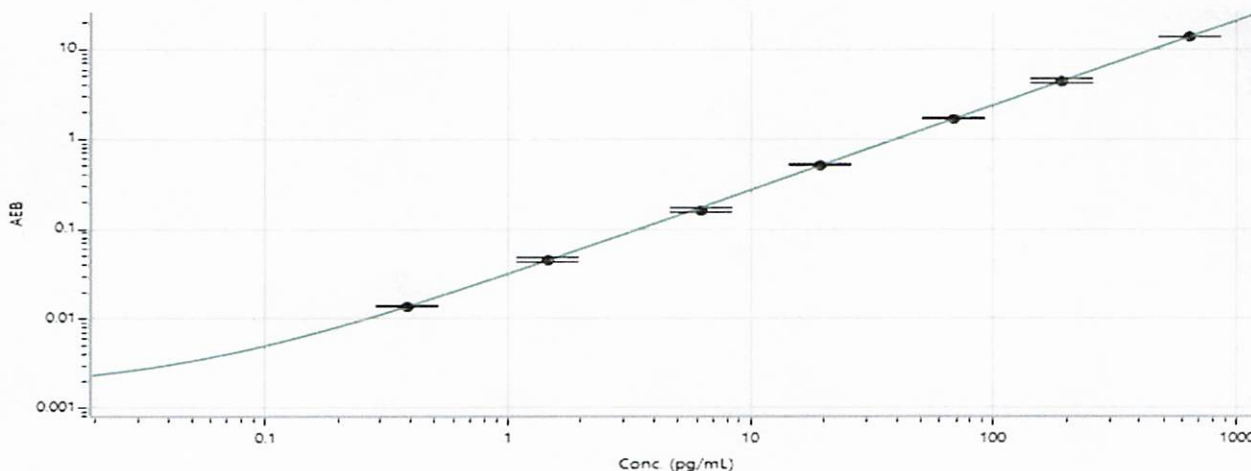
Component	Part Number	Lot Number	Lot Number
Bead Reagent	104067	505107	N/A
Detector Reagent	104068	505207	N/A
SBG Reagent	104069	505106	N/A
Sample Diluent Reagent	104071	505206	N/A
Reference Calibrators	102255	504211	N/A
Control 1	102256	504212	504212A
Control 2	102257	504212	504212A
RGP Reagent	103159	N/A ¹	N/A ²

¹ Reagents are not Kit Lot Specific

Release Materials*			
NF-Light®			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	504212A	4.50	3.43 - 6.33
Control 2	504212A	337	242 - 362

* 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	0.383
C	1.46
D	6.19
E	19.2
F	67.7
G	189
H	639

Review/Approval

Kyle Giroux
Name

Senior Document Control Specialist, QA
Title

[Signature] 13 Mar 2025
Signature/Date