

Product Number:	103575
Lot Number:	504110
Expiration:	11-Dec-2024
Platform(s):	SR-X and HD-1/HD-X



Component	Part Number	Lot Number
Bead Reagent	103564	402921
Detector Reagent	103566	403007
SBG Reagent	103568	403008
Sample Diluent	103569	401619
Calibrator Diluent	103571	402217
Calibrator Concentrate	103570	403003
SBG Diluent (SRX)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

† Specific to use on the SRX platform.

Refer to the SNAP-25 Advantage SR-X Kit Instructions (KI-0149).

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

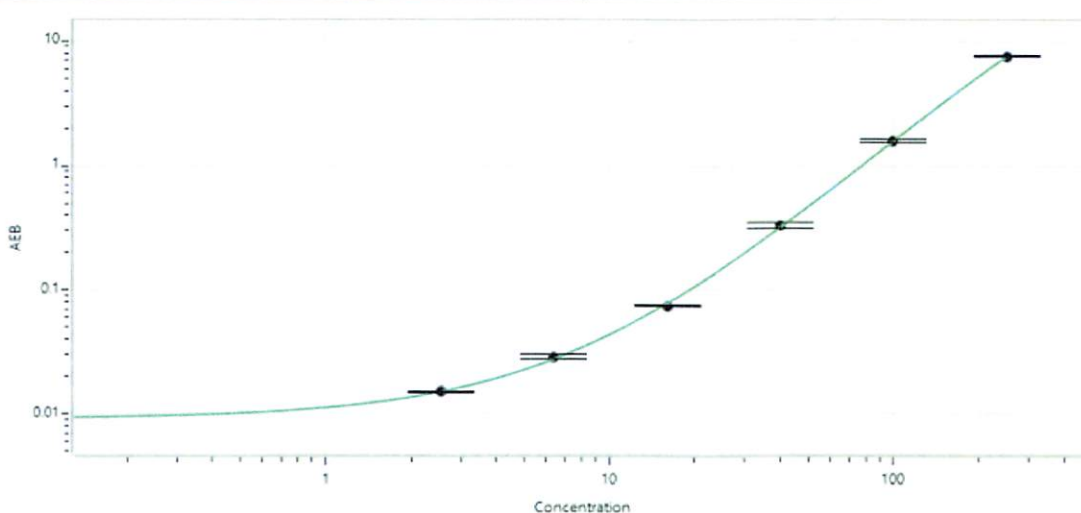
Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	403003	49.2	37.5 - 62.5
Control 2	403003	403	300 - 500
Panel 1	313805	36.2	29.4 - 46.9
Panel 2	N/A	N/A	N/A
Panel 3	313805	487	408 - 613

* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibrator Concentrate Value Assignment (pg/mL)	3819
--	------

Calibrator Levels (pg/mL)

A	0.000
B	2.56
C	6.40
D	16.0
E	40.0
F	100
G	250



Review/Approval

Cody Carlson
Name

QC Manager
Title

[Signature] 08 FEB 2024
Signature/Date