

<b>Product Number:</b>	103575
<b>Lot Number:</b>	504404
<b>Expiration:</b>	16-Jul-2025
<b>Platform(s):</b>	SR-X and HD-1/HD-X



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

Component	Part Number	Lot Number
Bead Reagent	103564	426706
Detector Reagent	103566	426909
SBG Reagent	103568	427703
Sample Diluent	103569	401623
Calibrator Diluent	103571	423224
Calibrator Concentrate	103570	427602
SBG Diluent (SR-X)†	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

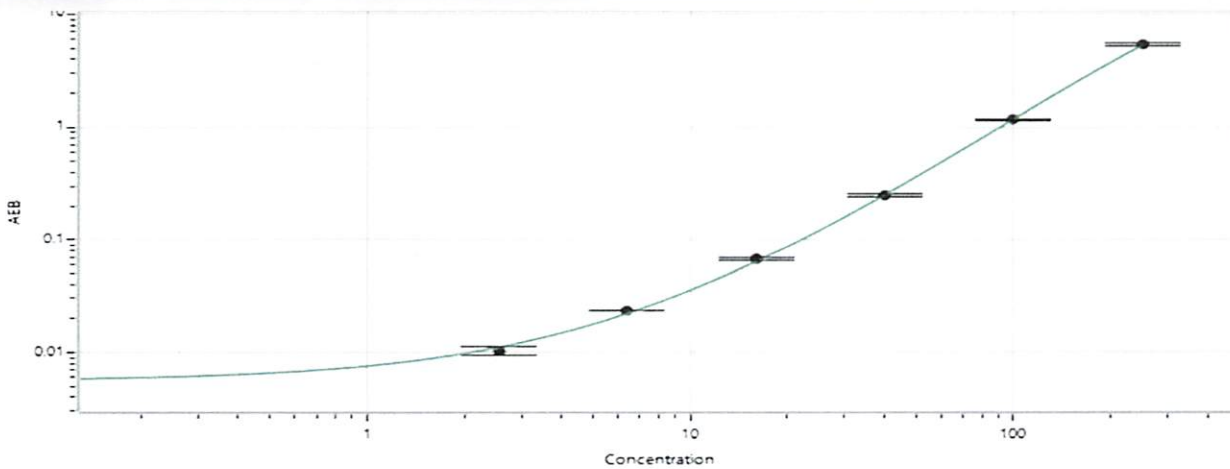
<sup>1</sup> Reagents are not Kit Lot Specific

† Specific to use on the SR-X platform.  
Refer to the SNAP-25 Advantage SR-X Kit Instructions (KI-0149).

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	427602	53.6	37.5 - 62.5
Control 2	427602	434	300 - 500

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

**Calibrator Concentrate Value Assignment (pg/mL)** 4923



**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

*Dipesh Patel* (Signature) *Doc control IT QA* (Title) *[Signature]* 27 Nov 2024 (Signature/Date)