

<b>Product Number:</b>	103293
<b>Lot Number:</b>	504209
<b>Expiration:</b>	8-Sep-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	103281	409205
Detector Reagent	103283	409305
SBG Reagent	103285	409306
Sample Diluent	103286	406817
Calibrator Diluent	103288	408516
Calibrator Concentrate	103287	409601
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

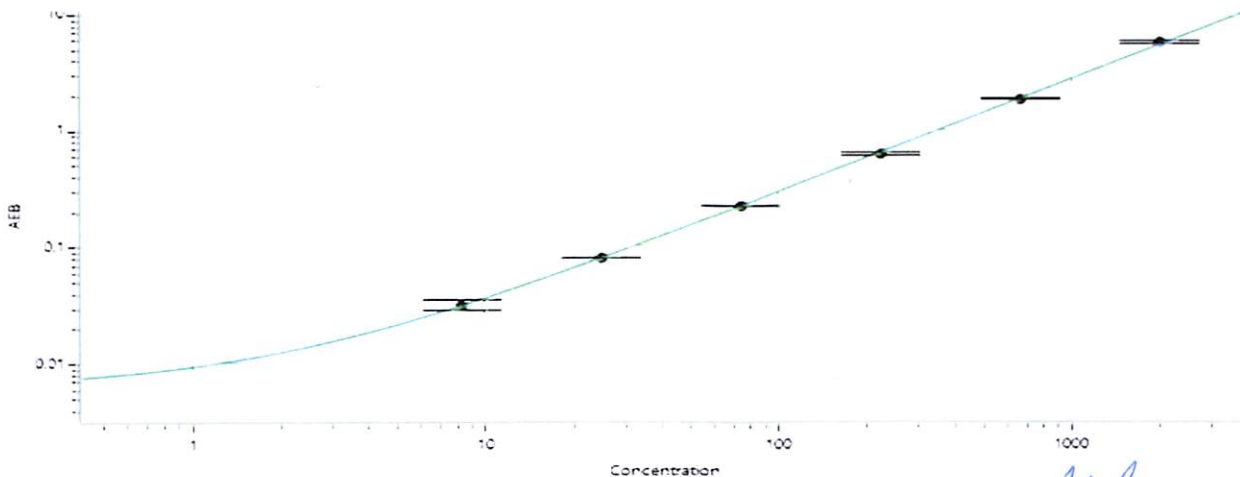
† Specific to use on the SRX platform.

Refer to the Simoa TDP-43 Kit Instructions SR-X (KI-0132).

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	409601	1988	1500 - 2500
Control 2	409601	208	150 - 250
Panel 1	406803	56.1	43.2 - 109
Panel 2	406803	95.9	72.4 - 187
Panel 3	406803	192	166 - 325
Panel 4	406803	3483	3000 - 4653

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

<b>Calibrator Concentrate Value Assignment (pg/mL)</b>	38548
--	-------



**Calibrator Levels (pg/mL)**

A	0.000
B	8.23
C	24.7
D	74.1
E	222
F	667
G	2000

Review/Approval

Cody Carlson

QC Manager

Name

Title

*[Signature]* 12 APR 2024

Signature/Date