

<b>Product Number:</b>	103293
<b>Lot Number:</b>	504558
<b>Expiration:</b>	3-Sep-2026
<b>Platform(s):</b>	SR-X and HD-1/HD-X



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

Component	Part Number	Lot Number	Control Lot Number
Bead Reagent	103281	508413	N/A
Detector Reagent	103283	506217	N/A
SBG Reagent	103285	506216	N/A
Sample Diluent Reagent	103286	506218	506218A
Calibrator Diluent Reagent	103288	506313	N/A
Calibrator Concentrate	103287	506416	N/A
SBG Diluent (SR-X)†	103158	N/A <sup>1</sup>	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>	N/A <sup>1</sup>

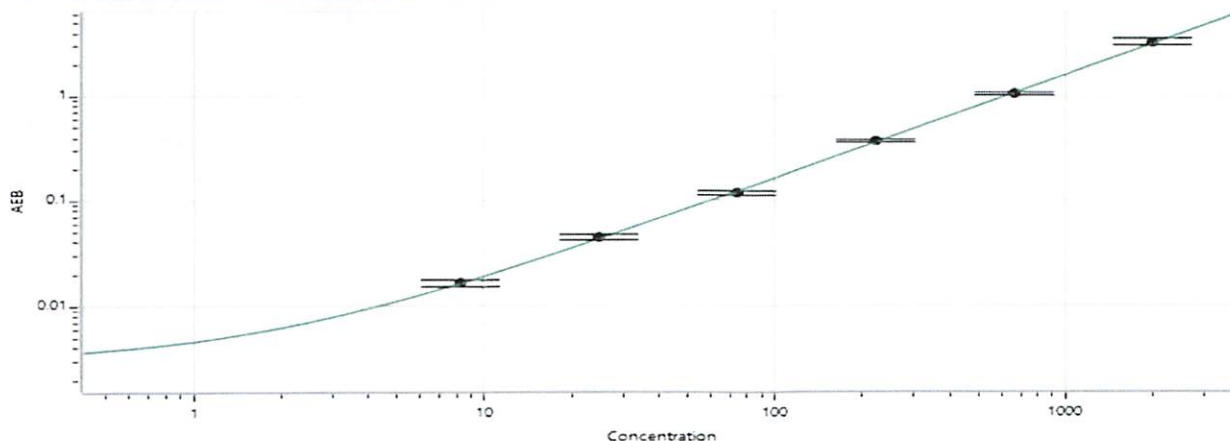
<sup>1</sup> Reagents are not Kit Lot Specific

† Specific to use on the SR-X platform.  
Refer to the Simoa TDP-43 Kit Instructions SR-X (KI-0132).

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	506416	187	150 - 250
Control 2	506416	1911	1500 - 2500

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

**Calibrator Concentrate Value Assignment (pg/mL)** 38510



**Calibrator Levels (pg/mL)**

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

Review/Approval

*Dhruvi Patel* *DOC Control-II, QA* *DP 02 Apr 2025*

Name

Title

Signature/Date