

Product Number:	103293
Lot Number:	504539
Expiration:	3-Sep-2026
Platform(s):	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	506501	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 ¹	N/A ¹
RGP Reagent	103159	N/A ¹	N/A ¹

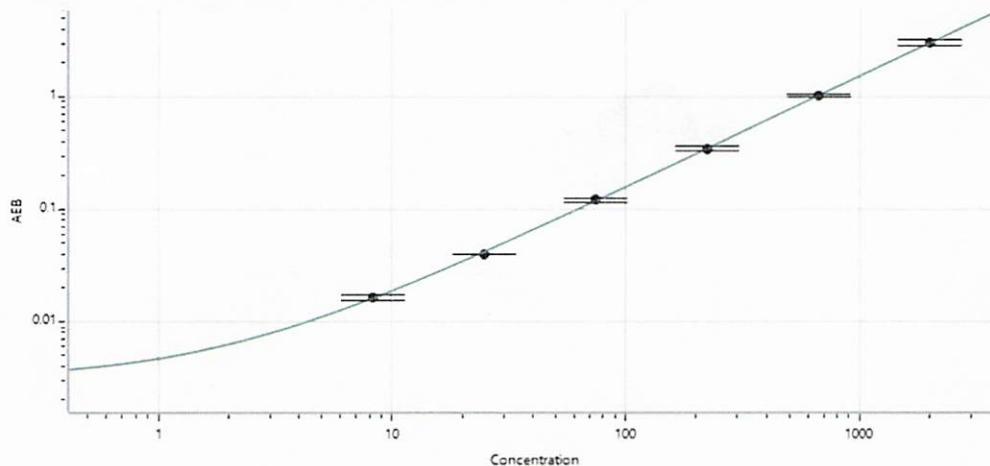
¹ Reagents are not Kit Lot Specific

† 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	191	150 - 250
Control 2	506416	1993	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

Kyle Giroux Senior Document Control Specialist QA *Kyle Giroux* 25 Mar 2025
Name Title Signature/Date