

Product Number:	103293
Lot Number:	503818
Expiration:	1-Mar-2024
Platform(s):	HD-1/HD-x & SR-X



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

Component	Part Number	Lot Number	Control Number
Bead Reagent	103281	316707	316707A
Detector Reagent	103283	306806	N/A
SBG Reagent	103285	306807	N/A
Sample Diluent	103286	306118	N/A
Calibrators Diluent	103288	306009	N/A
Calibrator Concentrate	103287	310011	N/A
SBG Diluent (SR-X) †	103158	N/A	N/A
RGP Reagent	103159	N/A ¹	N/A ¹

¹ RGP is not Kit Lot Specific

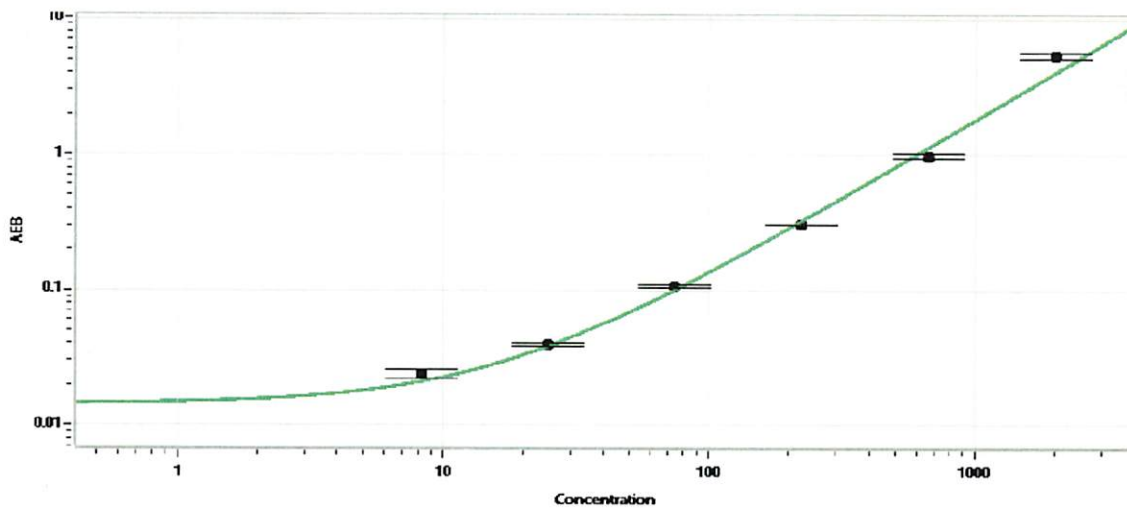
† Specific for use on the SR-X platform - refer to the Simoa™ TDP-43 SR-X Kit Instructions (KI-0132).

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	106304	109	70.8 - 127
Control 2	106304	1061	510 - 1320
Panel 2	231407	52.8	18.9 - 71.0
Panel 3	232608	204	122 - 239

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

Calibrator Concentrate Value Assignment (pg/mL)	15537
---	-------

Calibration Curve



Calibrator Levels (pg/mL)

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

Review/Approval

ROBERT SOMOGYI
Name

QC MANAGER
Title

[Signature] 04 AUG 2023
Signature/Date