

Product Number:	103293
Lot Number:	503660
Expiration:	18-Aug-2023
Platform(s):	HD-1/HD-x & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	103281	223801
Detector Reagent	103283	227211
SBG Reagent	103285	227212
Sample Diluent	103286	227214
Calibrators Diluent	103288	223010
Calibrator Concentrate	103287	233612
SBG Diluent (SR-X) †	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

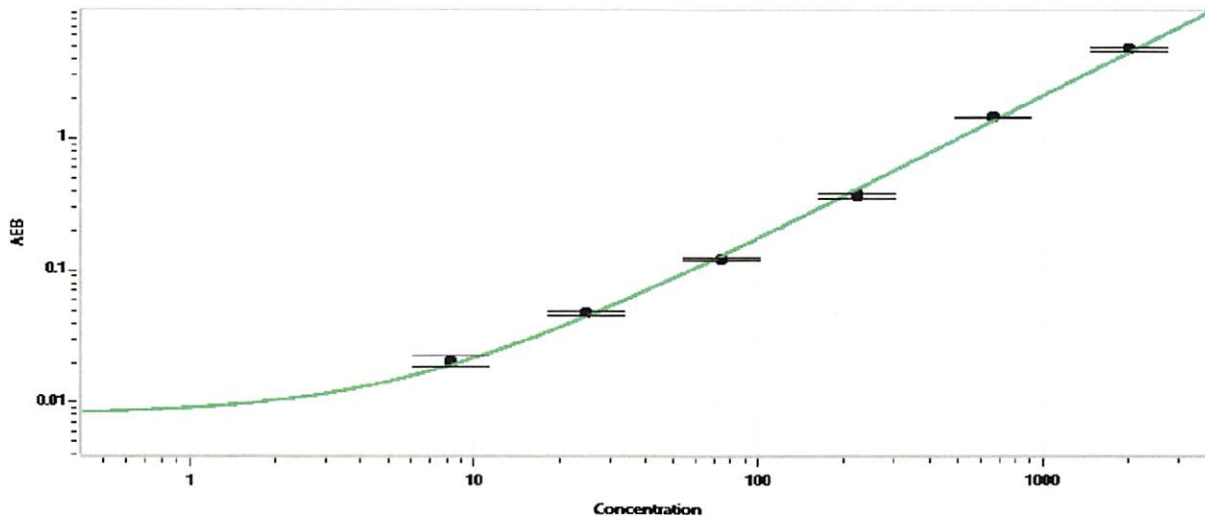
† Specific for use on the SR-X platform - refer to the Simoa™ TDP-43 SR-X Kit Instructions (KI-0132).

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	106304	81.4	70.8 - 127
Control 2	106304	695	510 - 1320
Panel 2	231407	31.4	18.9 - 71.0
Panel 3	232608	157	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)	22,287
---	--------

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: Sonia Khillan Name, QC Title, Sonia Khillan 1/29 DEC 2022 Signature/Date