

Product Number:	103293
Lot Number:	503262
Expiration:	30-Jul-2022
Platform(s):	HD-1/HD-x & SR-X



Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	103281	129001
Detector Reagent	103283	129011
SBG Reagent	103285	129012
Sample Diluent	103286	121116
Calibrators Diluent	103288	128015
Calibrator Concentrate	103287	129601
SBG Diluent (SR-X) †	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

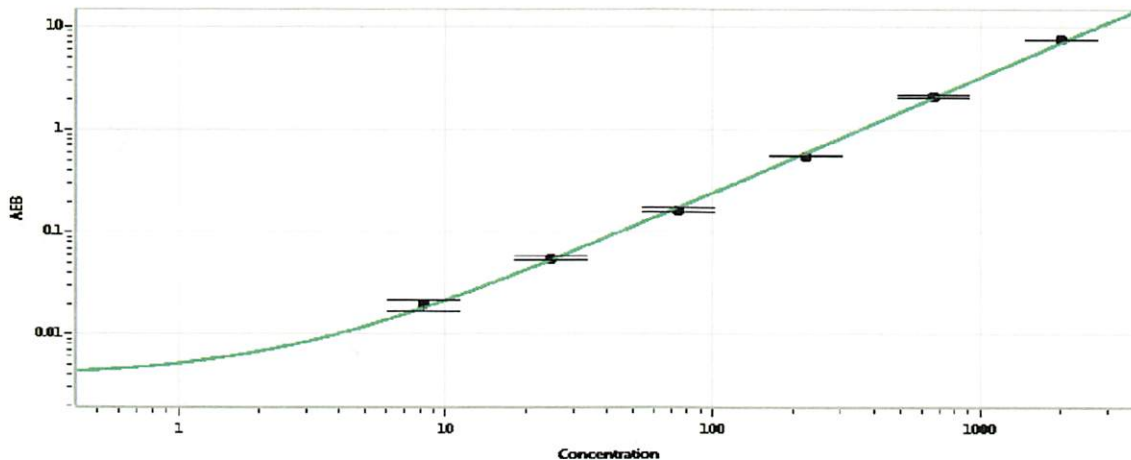
† Specific for use on the SR-X platform - refer to the Simoa™ TDP-43 SR-X Kit Instructions (KI-0132).

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	106304	90.0	70.8-127
Control 2	106304	936	510-1320
Panel 1	108513	41.2	19.4-67.0
Panel 2	108513	56.7	31.0-82.3
Panel 3	108513	529	312-807
Panel 4	108513	3926	1860-5614

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

Calibrator Concentrate Value Assignment (pg/mL)	36653
---	-------

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 Shivani Goel QC Shivani Goel 08Dec2021
Name Title Signature/Date