

Product Number:	103293
Lot Number:	503756
Expiration:	28-Oct-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	234012
Detector Reagent	103283	234106
SBG Reagent	103285	234107
Sample Diluent	103286	230114
Calibrators Diluent	103288	231916
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

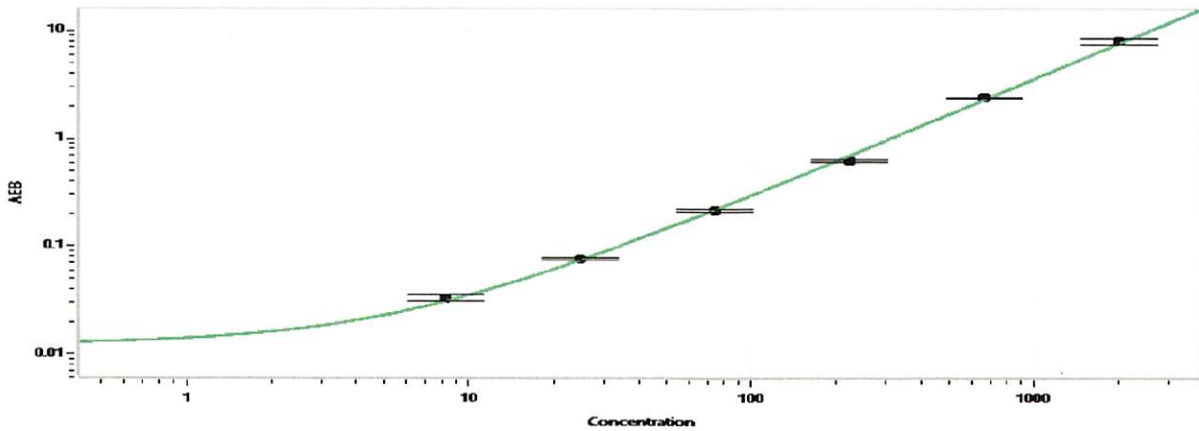
† Specific for use on the SR-X platform - refer to the Simoa™ TDP-43 SR-X Kit Instructions (KI-0132).

Release Materials*			
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1	106304	99.0	70.8 - 127
Control 2	106304	1216	510 - 1320
Panel 2	231407	53.0	18.9 - 71.0
Panel 3	232608	190	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 12/6/22 1/29 DEC 2022
Signature/Date