

Product Number:	103293
Lot Number:	502929
Expiration:	21-Apr-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	112707
Detector Reagent	103283	112715
SBG Reagent	103285	112716
Sample Diluent	103286	111118
Calibrators Diluent	103288	111725
Calibrator Concentrate	103287	114518
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

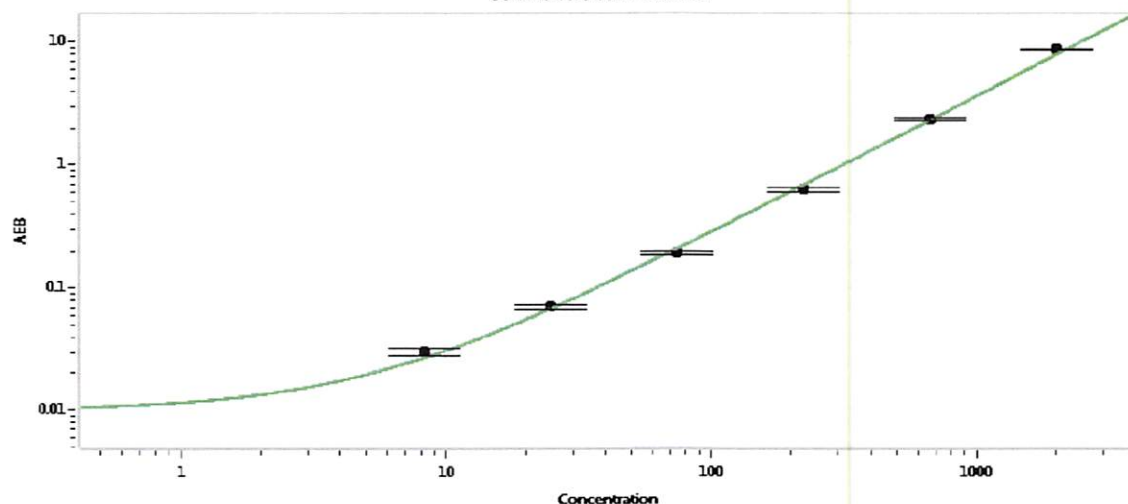
† Specific for use on the SR-X platform - refer to the Simoa™ TDP-43 SR-X Kit Instructions (KI-0132).

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	106304	86.4	73.6 - 118
Control 2	106304	923	593 - 1123
Panel 1	108513	36.1	20.8 - 60.8
Panel 2	108513	49.8	35.7 - 70.5
Panel 3	108513	471	408 - 632
Panel 4	108513	3287	2151 - 4586

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

Calibrator Concentrate Value Assignment (pg/mL)	22901
---	-------

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

Junpil Park
Name

QC Supervisor
Title

Signature/Date

09 Jun 2021