

Product Number:	103293
Lot Number:	503055
Expiration:	21-Mar-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	119709
Detector Reagent	103283	119804
SBG Reagent	103285	119802
Sample Diluent	103286	115223
Calibrators Diluent	103288	119415
Calibrator Concentrate	103287	114518
SBG Diluent (SR-X)†	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> RGP is not Kit Lot Specific

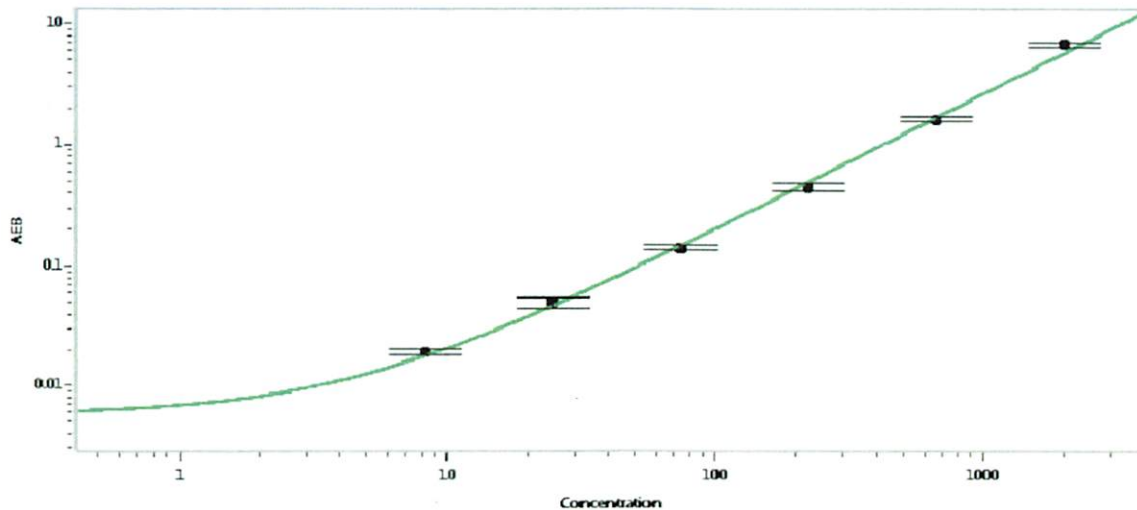
† Specific for use on the SR-X platform - refer to the Simoa™ TDP-43 SR-X Kit Instructions (KI-0132).

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	106304	115	70.8-127
Control 2	106304	1201	510-1320
Panel 1	108513	54.4	19.4-67.0
Panel 2	108513	73.3	31.0-82.3
Panel 3	108513	754	312-807
Panel 4	108513	5202	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)	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

23 Jul 2021