

Product Number:	103293
Lot Number:	502847
Expiration:	26-Jan-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	105303
Detector Reagent	103283	106407
SBG Reagent	103285	106502
Sample Diluent	103286	104619
Calibrators Diluent	103288	104726
Calibrator Concentrate	103287	106308
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

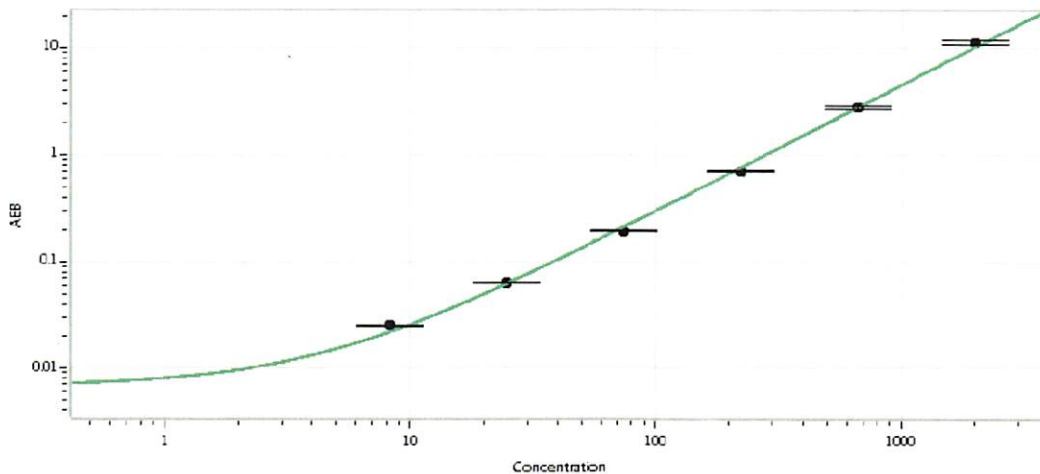
† Specific for use on the SR-X platform - refer to the Simoa™ TDP-43 SR-X Kit Instructions (KI-0132).

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	106304	96.5	77.6-118
Control 2	106304	854	561-1122
Panel 1	108513	37.8	22.5-61.8
Panel 2	108513	47.7	35.2-72.4
Panel 3	108513	488	415-624
Panel 4	108513	3232	2073-4705

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

Calibrator Concentrate Value Assignment (pg/mL)	26695
---	-------

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

20 Apr 2021