

Product Number:	103293
Lot Number:	502063
Expiration:	17-Sep-2019
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	928017
Detector Reagent	103283	928201
SBG Reagent	103285	927707
Sample Diluent	103286	927708
Calibrators Diluent	103288	926407
Calibrator Concentrate	103287	833803
SBG Diluent (SR-X)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

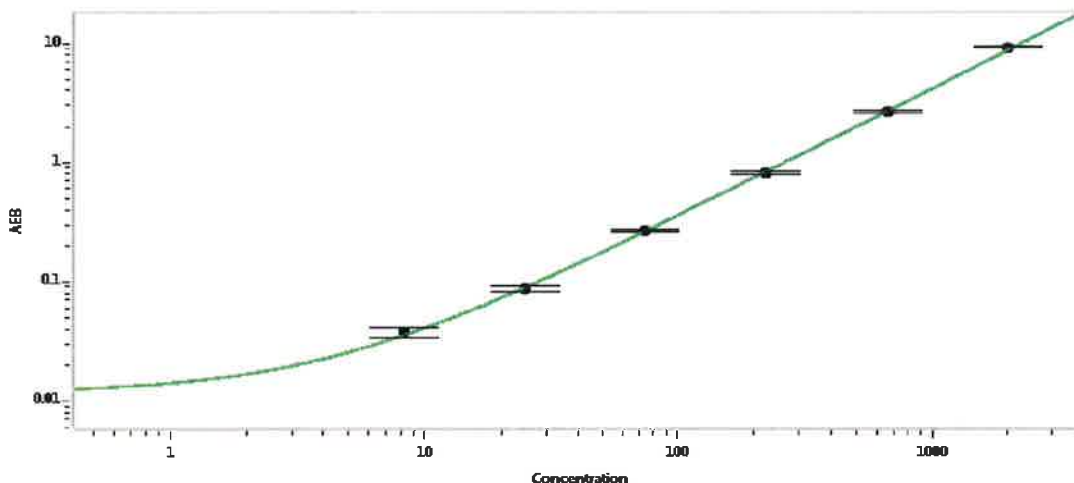
† Specific for use on the SR-X platform - refer to the Simoa™ TDP-43 SR-X Kit Instructions (KI-0132).

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	834401	220	168-251
Control 2	830401	1791	1314-1971
Panel 1	831201	38.6	30.6-72.3
Panel 2	831202	69.9	46.6-93.6
Panel 3	832006	293	200-339
Panel 4	832007	1441	1125-1734

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

Calibrator Concentrate Value Assignment (pg/mL) 34,946 pg/mL

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

Patrizia Stadler Assoc. Director, QC
Name Title

P. Stadler 21 Oct 2019
Signature/Date