

Product Number:	103293
Lot Number:	501511
Expiration:	19-Mar-2019
Platform(s):	HD-1 & SR-X



Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	103281	832406
Detector Reagent	103283	832305
SBG Reagent	103285	832306
Sample Diluent	103286	826219
Calibrator Diluent	103288	829713
Calibrator Concentrate	103287	833803
SBG Diluent (SR-X) †	103158	N/A ¹
RGP Reagent	103159	N/A ¹

¹ Reagents are not Kit Lot Specific

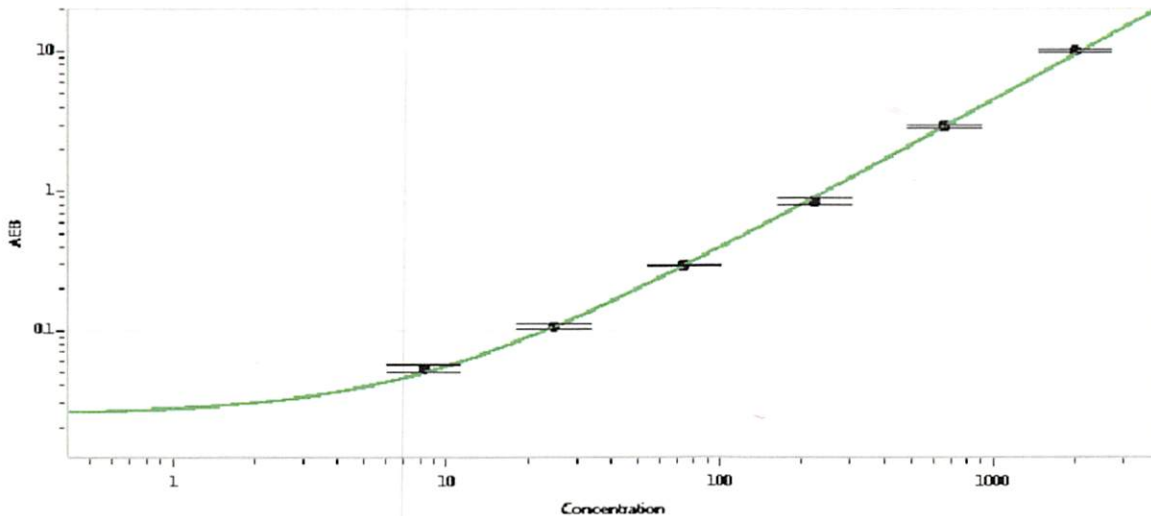
† Specific for use on the SR-X platform - refer to the Simoa™ TDP-43 SR-X Kit Instructions (KI-0132).

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	834401	197	166-249
Control 2	830401	1732	1294-1941
Panel 1	831201	50.7	36.3-70.1
Panel 2	831202	68.0	45.5-94.8
Panel 3	832006	297	196-335
Panel 4	832007	1567	1096-1758

* 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

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature]
Signature/Date