

Product Number:	100794
Lot Number:	501531
Expiration:	19-Sep-2019
Platform(s):	HD-1 and SR-X

Component	Part Number	Lot Number
Bead Reagent	100781	833104
Detector Reagent	100782	833103
SBG Reagent	100783	833204
Sample Diluent	103202	826220
Calibrator Diluent	100793	829714
Calibrator Concentrate	100780	80254
SBG Diluent (SRX)†	103158	N/A ¹
RGP Reagent	103159	N/A ¹



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

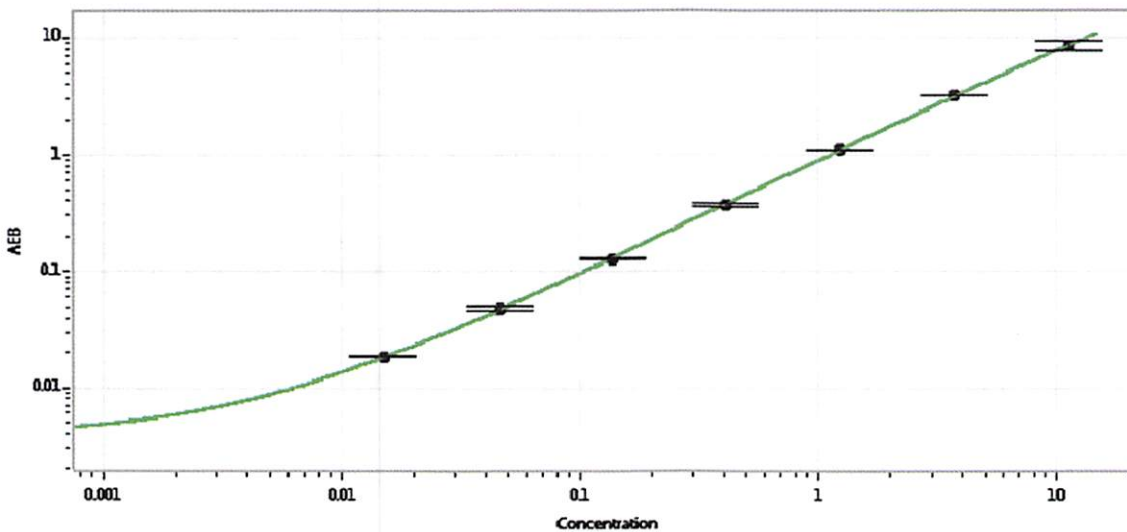
¹ Reagents are not Kit Lot Specific.

† Specific to use on the SR-X platform - refer to the SR-X Kit Instructions (KI-0085).

Release Materials*				* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Control 1	MF-062316-02	7.29	7.15-10.7	
Control 2	MF-062316-02	0.659	0.593-0.889	
Panel 1	MF-110316-05	3.90	3.68-5.52	
Panel 2	MF-110316-06	45.1	40.8-61.2	
Panel 3	MF-110316-07	3.43	3.25-4.88	

Calibrator Concentrate Value Assignment (pg/mL)	829
---	-----

Calibration Curve



Review/Approval

Sivasankari Murugan QC Supervisor
Name Title

[Signature] 11 DEC 2018
Signature/Date