

Product Number:	101672
Lot Number:	501990
Expiration:	14-Mar-2020
Platform(s):	HD-1/HD-X



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

Component	Part Number	Lot Number
Bead Reagent	101674	911907
Detector Reagent	101673	912109
SBG Reagent	101675	912110
Sample Diluent	101676	907316
Calibrator Diluent	101677	910310
Calibrator Concentrate	101734	911207
RGP Reagent	103159	N/A ¹

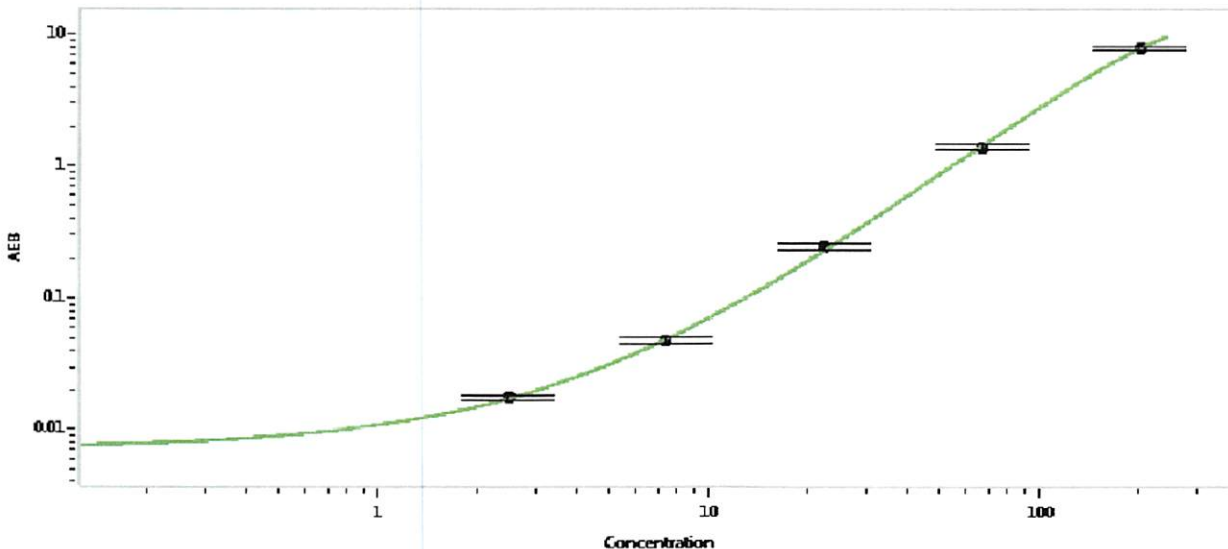
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-021816-02	24.1	17.1-28.3
Control 2	MF-021816-02	343	263-434
Panel 1	MF-040518-01	195	138-207
Panel 2	MF-040518-02	40.4	33.8-60.0

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

Calibrator Concentrate Value Assignment (pg/mL)	5847
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	2.47
C	7.41
D	22.2
E	66.7
F	200

Review/Approval

Daniel Ritchie Senior QC Analyst
Name Title

Signature/Date

*Signing on behalf of Patrizia Stadler