

Product Number:	101672
Lot Number:	501599
Expiration:	19-Sep-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	101674	834809
Detector Reagent	101673	901802
SBG Reagent	101675	834503
Sample Diluent	101676	826223
Calibrator Diluent	101677	829716
Calibrator Concentrate	101734	73205
RGP Reagent	103159	N/A ¹

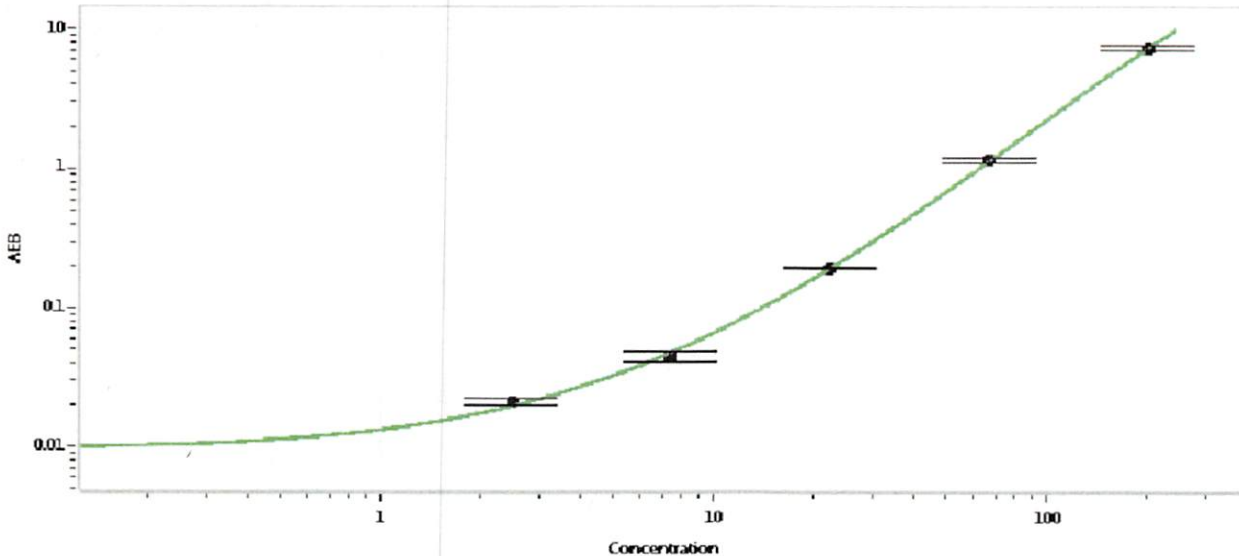
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-021816-02	21.2	17.1-28.3
Control 2	MF-021816-02	332	263-434
Panel 1	MF-040518-01	174	140-211
Panel 2	MF-040518-02	54.5	34.7-60.3

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

Calibrator Concentrate Value Assignment (pg/mL)	2711
---	------

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

Patrizia Stadler
Name
Associate Director, QC
Title

P. Stadler
Signature
22 Jan 2019
Date