

Product Number:	101599
Lot Number:	501989
Expiration:	27-Feb-2020
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	101601	909503
Detector Reagent	102012	910601
SBG Reagent	101602	909815
Sample Diluent	101603	910210
Calibrator Diluent	102014	905826
Calibrator Concentrate	100778	904508
RGP Reagent	103159	N/A ¹

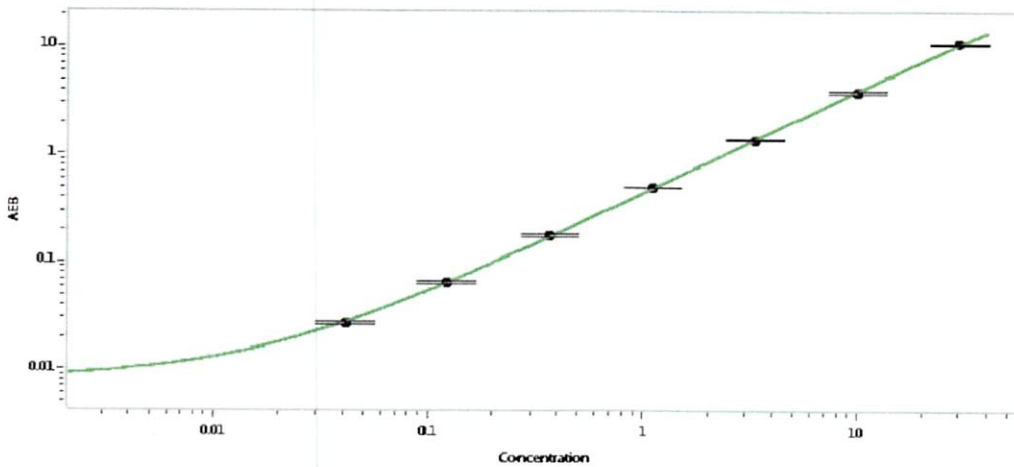
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-012518-02	0.452	0.365-0.548
Control 2	MF-012518-02	24.8	13.9-27.2
Panel 1	900901	0.958	0.744-2.05
Panel 2	900902	0.720	0.506-2.61
Panel 3	900903	9.40	6.48-9.90
Panel 4	900904	33.6	21.0-35.9

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

Calibrator Concentrate Value Assignment (pg/mL)	1006
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.041
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval: Daniel Ritchie Senior QC Analyst
 Name: Daniel Ritchie Title: Senior QC Analyst Signature/Date: [Signature] / 22 Aug 2019

*Signing on behalf of Patrizia Stadler