

Product Number:	101599
Lot Number:	501954
Expiration:	12-Jul-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	919703
Detector Reagent	102012	919606
SBG Reagent	101602	920707
Sample Diluent	101603	919305
Calibrator Diluent	102014	919008
Calibrator Concentrate	100778	919304
RGP Reagent	103159	N/A <sup>1</sup>

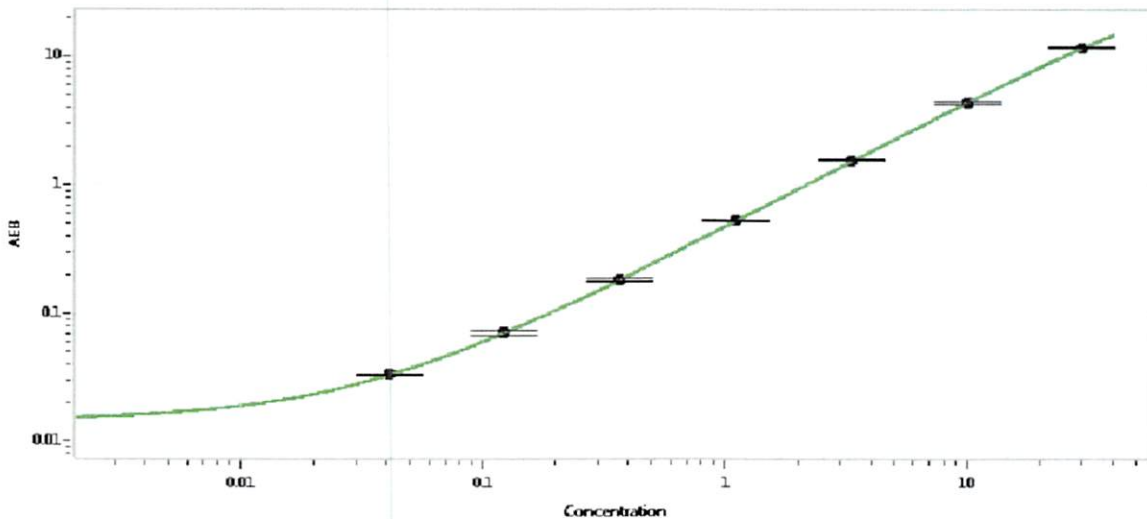
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-012518-02	0.377	0.365-0.548
Control 2	MF-012518-02	21.5	13.9-27.2
Panel 1	900901	2.00	0.392-3.05
Panel 2	900902	1.69	1.11-2.32
Panel 3	900903	8.54	6.42-9.63
Panel 4	900904	31.0	21.8-34.5

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

Calibrator Concentrate Value Assignment (pg/mL)	798
---	-----

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

Patrizia Stadler Assoc. Director, QC

*P. Stadler*

31 Jul 2019

Name

Title

Signature/Date