

Product Number:	102860
Lot Number:	502104
Expiration:	23-Mar-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	102852	926612
Detector Reagent	102854	930813
SBG Reagent	102855	930812
Sample Diluent	102856	930801
Calibrator Diluent	102859	930802
Calibrator Concentrate	102857	81423
RGP Reagent	103159	N/A <sup>1</sup>

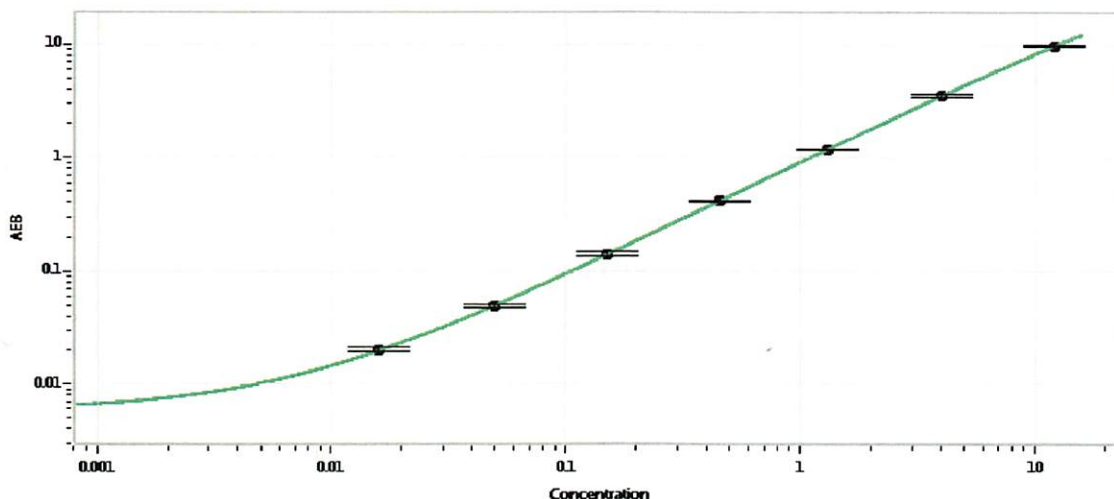
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-110817-02	0.300	0.247-0.370
Control 2	MF-110817-02	4.95	4.13-6.32
Panel 1	MF-100517-01	0.479	0.393-0.743
Panel 2	MF-100517-02	1.37	1.12-1.67
Panel 3	MF-100517-03	4.80	3.96-6.14

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

Calibrator Concentrate Value Assignment (pg/mL)	217
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.016
C	0.049
D	0.148
E	0.444
F	1.33
G	4.00
H	12.0

Review/Approval

Patrizia Stadler Assoc. Director, QC

*P. Stadler*

11 NOV 2019

Name

Title

Signature/Date