

Product Number:	102860
Lot Number:	501777
Expiration:	3-Oct-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	102852	909316
Detector Reagent	102854	910302
SBG Reagent	102855	909305
Sample Diluent	102856	909304
Calibrator Diluent	102859	909318
Calibrator Concentrate	102857	81423
RGP Reagent	103159	N/A <sup>1</sup>

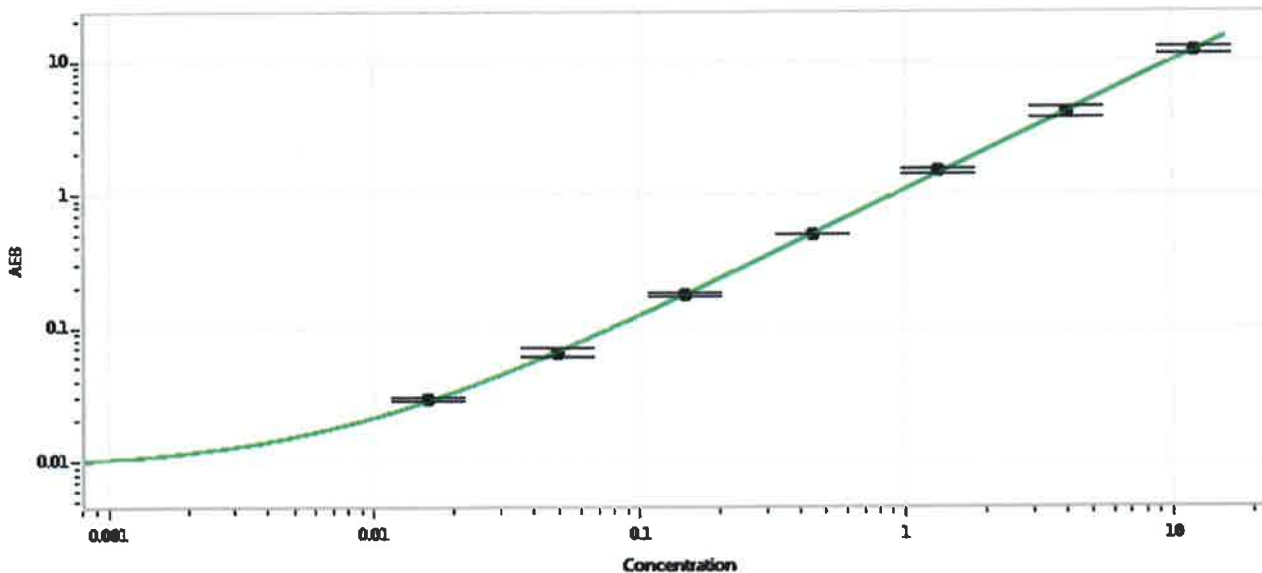
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-110817-02	0.284	0.248-0.372
Control 2	MF-110817-02	4.87	4.13-6.40
Panel 1	MF-100517-01	0.466	0.418-0.744
Panel 2	MF-100517-02	1.36	1.12-1.68
Panel 3	MF-100517-03	4.71	3.96-6.22

\* 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	QC Assoc. Director	<i>P. Stadler</i> 17 Apr 2019
	Name	Title	Signature/Date