

Product Number:	102860
Lot Number:	502004
Expiration:	22-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	102852	923804
Detector Reagent	102854	923807
SBG Reagent	102855	923806
Sample Diluent	102856	923412
Calibrator Diluent	102859	923413
Calibrator Concentrate	102857	81423
RGP Reagent	103159	N/A ¹

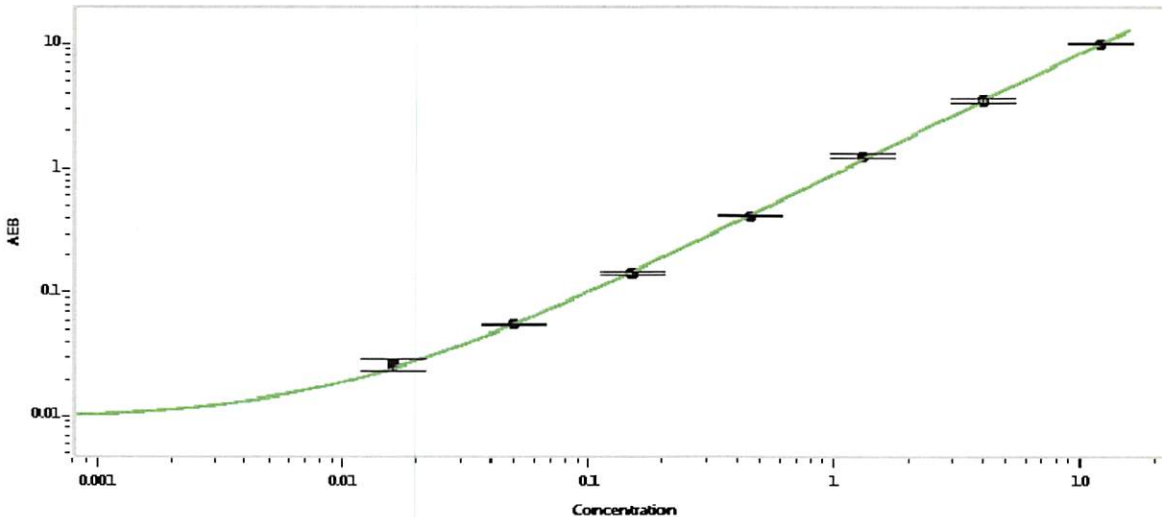
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-110817-02	0.263	0.229-0.379
Control 2	MF-110817-02	4.71	4.00-6.39
Panel 1	MF-100517-01	0.405	0.340-0.777
Panel 2	MF-100517-02	1.25	1.10-1.65
Panel 3	MF-100517-03	4.56	3.76-6.28

* 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 03 Sept 2019

Name

Title

Signature/Date