

Product Number:	102860
Lot Number:	502301
Expiration:	23-Sep-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	008309
Detector Reagent	102854	008301
SBG Reagent	102855	008308
Sample Diluent	102856	008407
Calibrator Diluent	102859	008408
Calibrator Concentrate	102857	008305
RGP Reagent	103159	N/A ¹

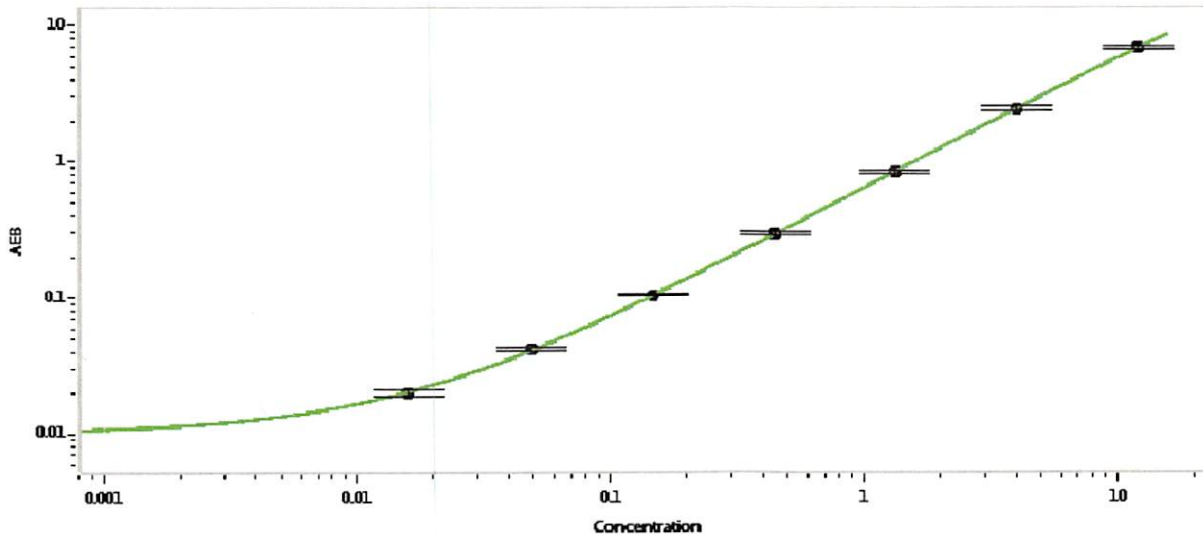
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-110817-02	0.294	0.247-0.371
Control 2	MF-110817-02	5.22	4.14-6.42
Panel 1	MF-100517-01	0.481	0.395-0.733
Panel 2	MF-100517-02	1.52	1.14-1.70
Panel 3	MF-100517-03	5.43	3.97-6.27

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

Calibrator Concentrate Value Assignment (pg/mL)	386
---	-----

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

Name

Title

P. Stadler

30 Mar 2020

Signature/Date