

Product Number:	102860
Lot Number:	502203
Expiration:	21-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	102852	002214
Detector Reagent	102854	002107
SBG Reagent	102855	002108
Sample Diluent	102856	002105
Calibrator Diluent	102859	002106
Calibrator Concentrate	102857	002401
RGP Reagent	103159	N/A ¹

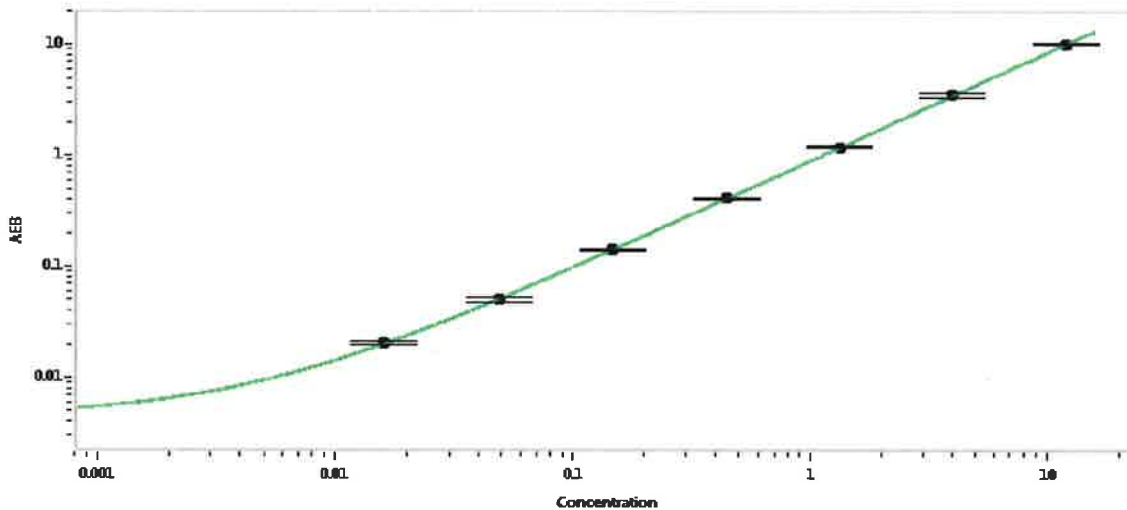
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-110817-02	0.311	0.247-0.371
Control 2	MF-110817-02	5.69	4.14-6.42
Panel 1	MF-100517-01	0.536	0.395-0.733
Panel 2	MF-100517-02	1.62	1.14-1.70
Panel 3	MF-100517-03	5.64	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)	240
---	-----

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 3 Feb 2020

Name

Title

Signature/Date