

Product Number:	102860
Lot Number:	502614
Expiration:	11-May-2021
Platform(s):	HD-1/HD-X



Data below represents results generated on the Simoa™ HD-X Analyzer.

Component	Part Number	Lot Number
Bead Reagent	102852	031608
Detector Reagent	102854	032504
SBG Reagent	102855	032206
Sample Diluent	102856	032211
Calibrator Diluent	102859	032212
Calibrator Concentrate	102857	002401
RGP Reagent	103159	N/A ¹

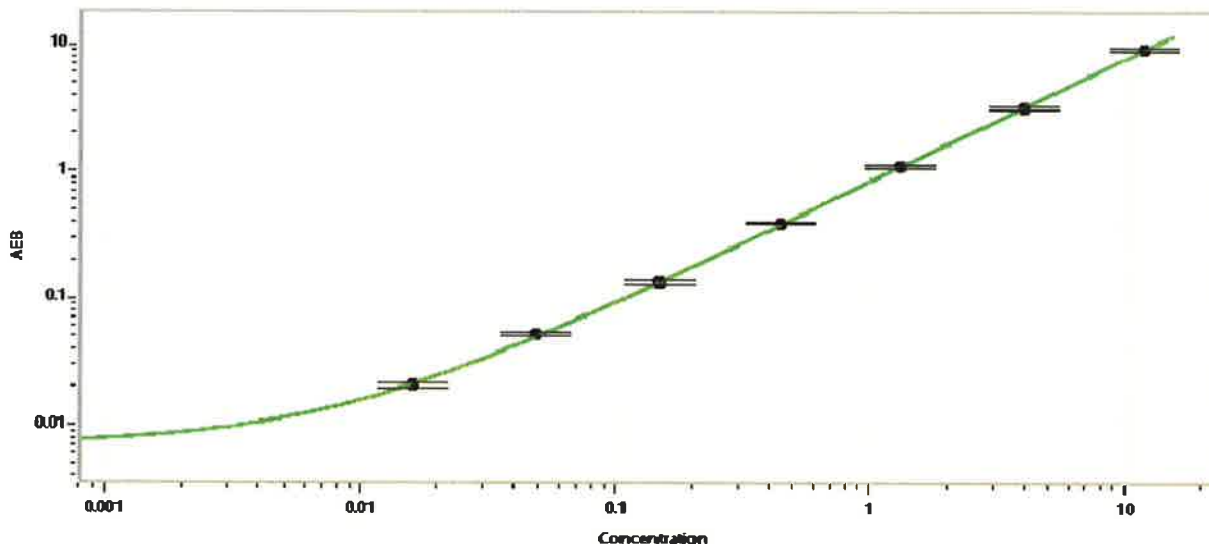
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-110817-02	0.306	0.247-0.371
Control 2	MF-110817-02	5.27	4.14-6.42
Panel 1	033802	0.153	0.114-0.258
Panel 2	033802	1.87	1.65-2.47
Panel 3	033802	9.54	7.87-11.8

* 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

Shivani Goel
Name

QC Manager
Title

Shivani Goel 09 Dec 2020
Signature/Date