

Product Number:	100988
Lot Number:	502357
Expiration:	18-Feb-2021
Platform(s):	HD-1/HD-X



Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	100974	012301
Detector Reagent	100976	011902
SBG Reagent	100987	011901
Sample Diluent	100985	005228
Calibrator Diluent	100980	004915
Calibrator Concentrate	100978	934402
RGP Reagent	103159	N/A ¹

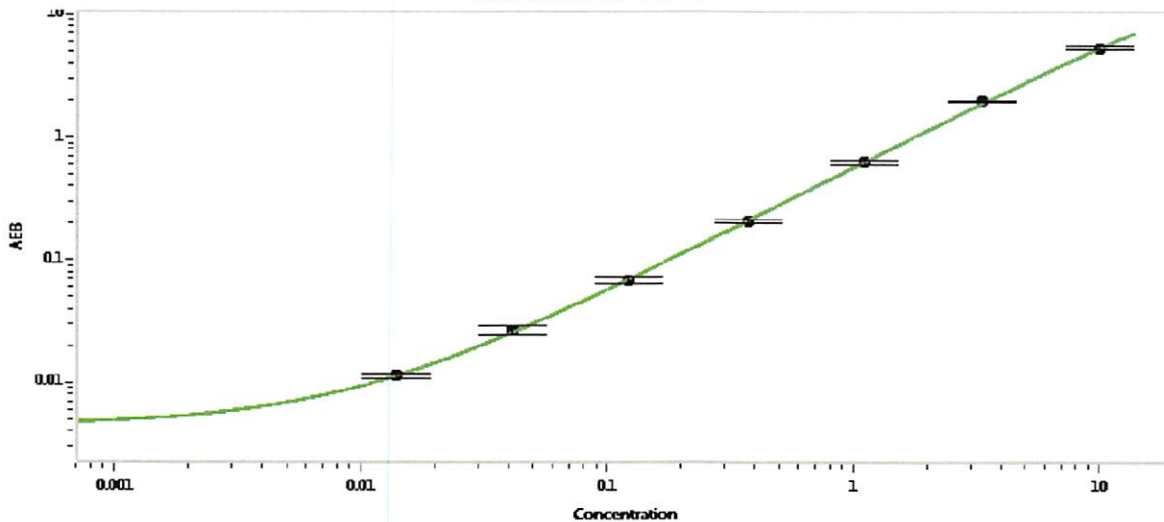
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-101816-02	0.131	0.068-0.161
Control 2	MF-101816-02	7.94	7.31-11.2
Panel 1	MF-101716-04	0.275	0.167-0.327
Panel 2	MF-101716-05	0.072	0.021-0.087
Panel 3	MF-101816-03	17.1	15.6-23.9

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

Calibrator Concentrate Value Assignment (pg/mL)	225
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.014
C	0.041
D	0.123
E	0.370
F	1.11
G	3.33
H	10.0

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 13 May 2020

Name

Title

Signature/Date