

Product Number:	101580
Lot Number:	502460
Expiration:	2-Apr-2021
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	101582	018807
Detector Reagent	101581	016703
SBG Reagent	101583	015902
Sample Diluent	103201	015412
Calibrator Diluent	101585	015901
Calibrator Concentrate	100655	909204
RGP Reagent	103159	N/A ¹

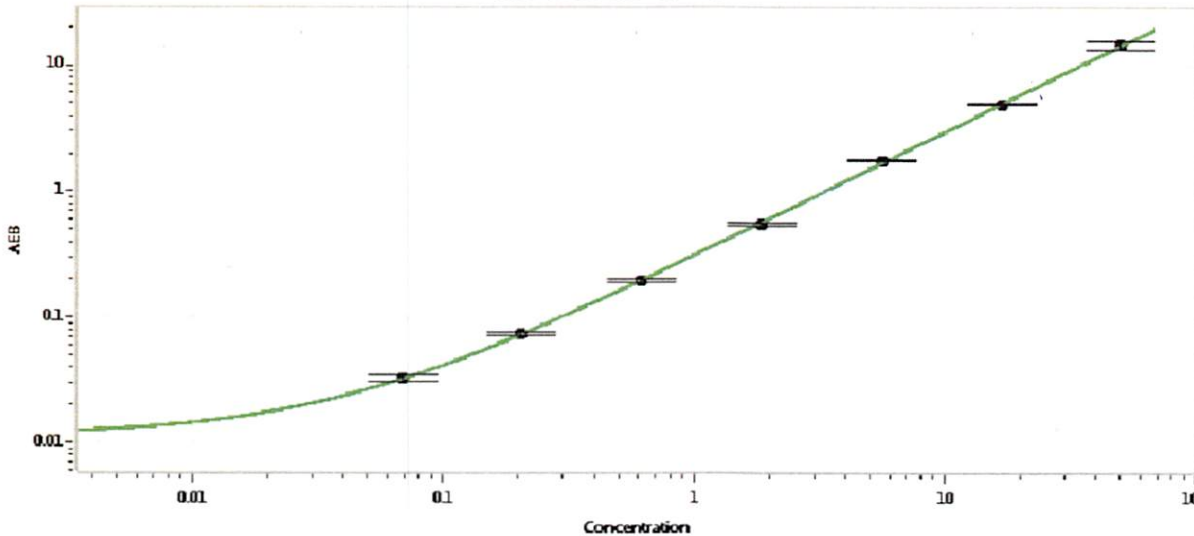
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-042817-01	2.73	2.08-3.12
Control 2	MF-051617-02	47.5	35.1-52.7
Panel 1	MF-051017-03	28.7	19.9-32.1
Panel 2	MF-032217-11	2.11	1.33-2.23
Panel 3	MF-033018-06	3.22	1.68-3.50

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

Calibrator Concentrate Value Assignment (pg/mL)	1065
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.069
C	0.206
D	0.617
E	1.85
F	5.56
G	16.7
H	50.0

Review/Approval

Shivani Goel

QC Manager

Shivani Goel 13 Jul 2020

Name

Title

Signature/Date