

Product Number:	101635
Lot Number:	502198
Expiration:	14-Jan-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	101637	001411
Detector Reagent	101636	001404
SBG Reagent	101638	001405
Sample Diluent	101639	001406
Calibrator Diluent	101640	001412
Calibrator Concentrate	100661	908402
RGP Reagent	103159	N/A ¹

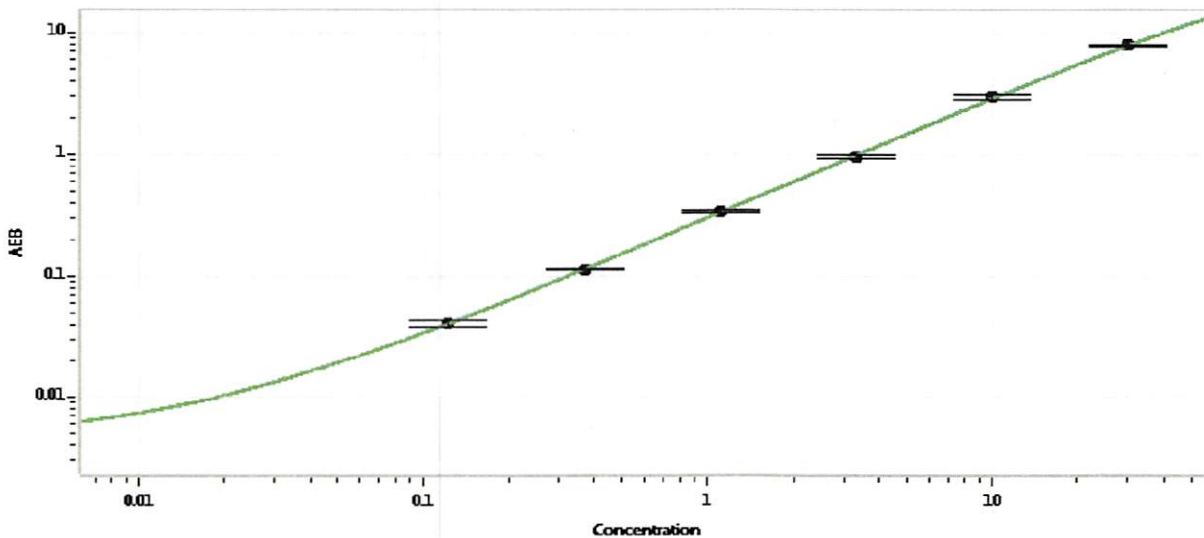
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	904406	0.645	0.550-0.825
Control 2	904406	69.1	54.1-81.2
Panel 1	MF-040618-02	0.546	0.414-0.657
Panel 2	MF-040618-03	4.92	4.18-6.26
Panel 3	MF-040618-04	52.1	44.1-66.2

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

Calibrator Concentrate Value Assignment (pg/mL)	1760
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.123
C	0.370
D	1.11
E	3.33
F	10.0
G	30.0

Review/Approval

Patrizia Stadler Assoc. Director, QC *P. Stadler* 29 Jan 2020
 Name Title Signature/Date