

Product Number:	101635
Lot Number:	501876
Expiration:	24-Jun-2020
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	917504
Detector Reagent	101636	917903
SBG Reagent	101638	917702
Sample Diluent	101639	917609
Calibrator Diluent	101640	917610
Calibrator Concentrate	100661	907709
RGP Reagent	103159	N/A ¹

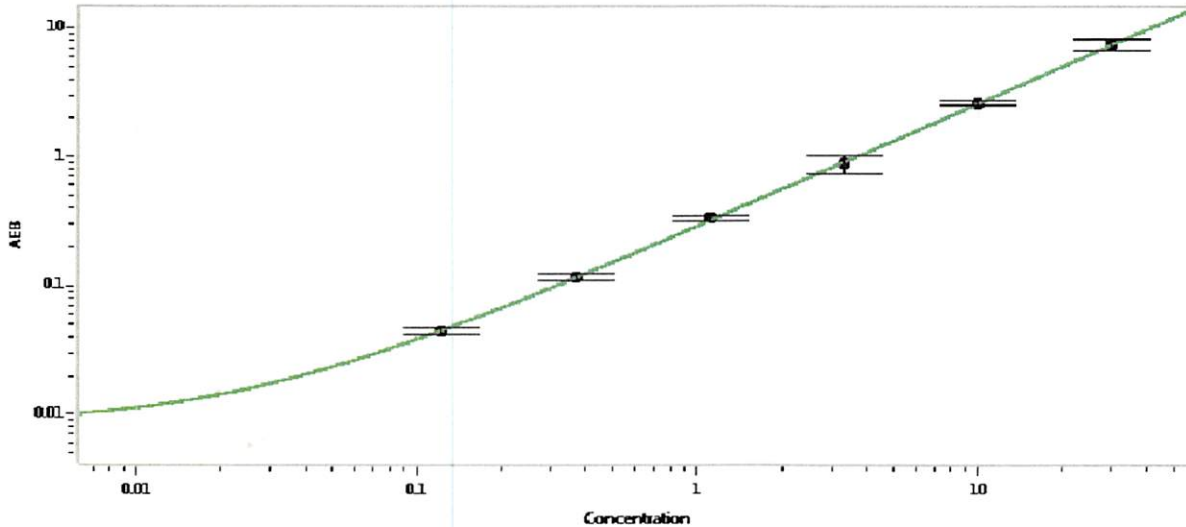
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	904406	0.662	0.554-0.831
Control 2	904406	70.6	54.0-81.0
Panel 1	MF-040618-02	0.562	0.413-0.655
Panel 2	MF-040618-03	5.49	4.21-6.31
Panel 3	MF-040618-04	60.8	44.4-66.6

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

Calibrator Concentrate Value Assignment (pg/mL)	1810
---	------

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
Name Title

P Stadler 09 Jul 2019
Signature/Date