

Product Number:	101635
Lot Number:	501895
Expiration:	29-Mar-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	910811
Detector Reagent	101636	909506
SBG Reagent	101638	912012
Sample Diluent	101639	908813
Calibrator Diluent	101640	908812
Calibrator Concentrate	100661	907709
RGP Reagent	103159	N/A ¹

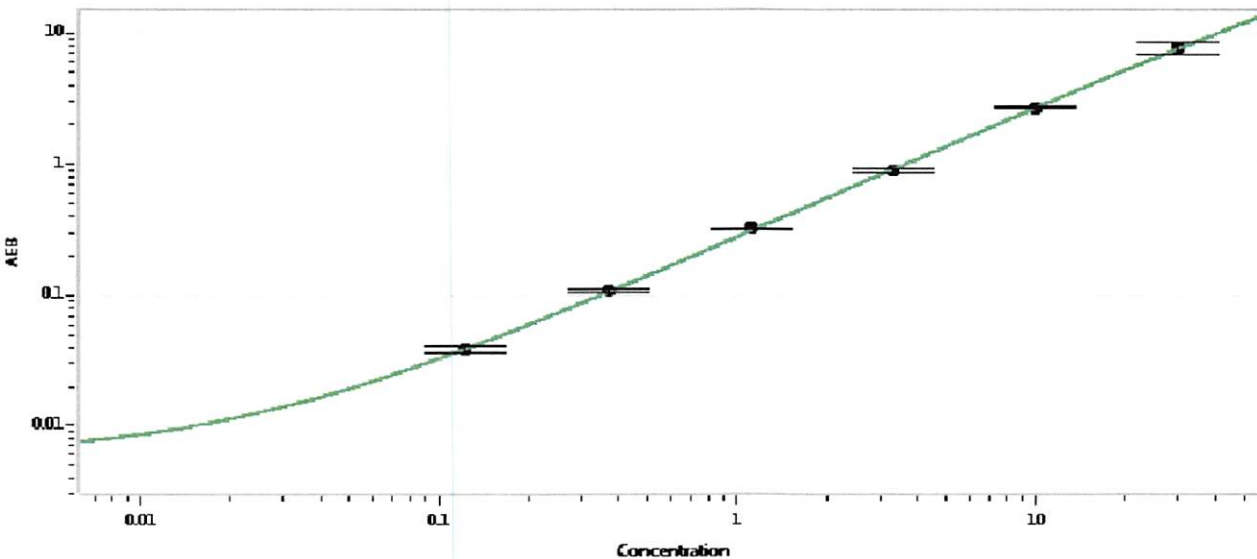
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	904406	0.734	0.558-0.837
Control 2	904406	71.6	53.6-80.4
Panel 1	MF-040618-02	0.54	0.407-0.654
Panel 2	MF-040618-03	5.66	4.18-6.27
Panel 3	MF-040618-04	54	43.8-65.7

* 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	<i>P. Stadler</i>	17 Jul 2019
	Name	Title	Signature/Date	