

Product Number:	101635
Lot Number:	502606
Expiration:	22-Oct-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	030903
Detector Reagent	101636	031124
SBG Reagent	101638	031117
Sample Diluent	101639	031005
Calibrator Diluent	101640	031008
Calibrator Concentrate	100661	031115
RGP Reagent	103159	N/A <sup>1</sup>

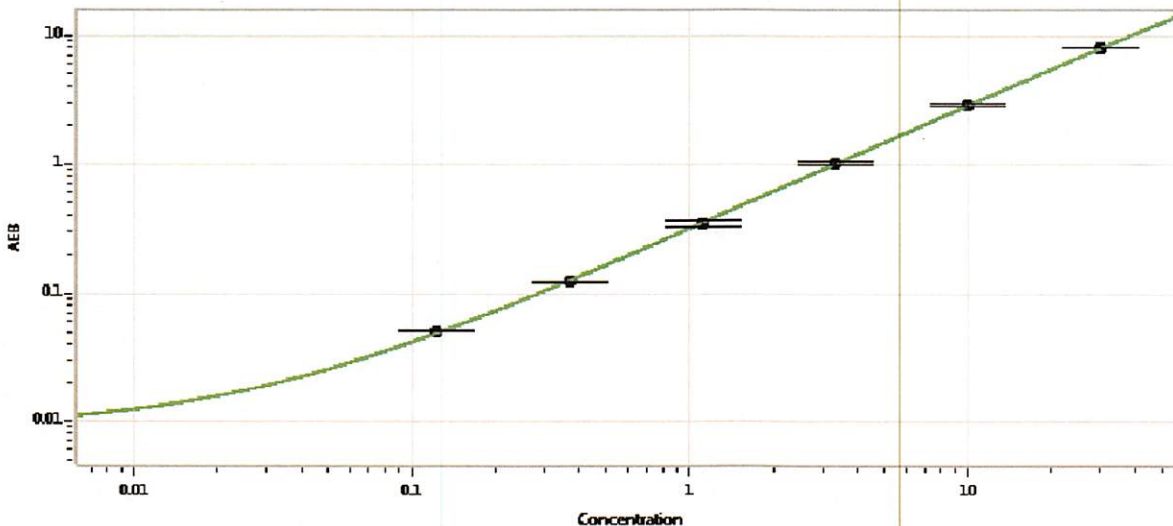
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	904406	0.553	0.551-0.827
Control 2	904406	54.8	54.7-82.0
Panel 1	MF-040618-02	0.459	0.434-0.651
Panel 2	MF-040618-03	4.32	4.19-6.28
Panel 3	MF-040618-04	43.6	43.5-65.3

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

Calibrator Concentrate Value Assignment (pg/mL)	1883
---	------

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
H	

Review/Approval

Patrizia Stadler Associate Director, QC

*P. Stadler* 18 Nov 2020

Name

Title

Signature/Date