

Product Number:	101605
Lot Number:	502125
Expiration:	19-Nov-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	101607	932509
Detector Reagent	101606	932308
SBG Reagent	101608	932307
Sample Diluent	103200	932305
Calibrator Diluent	101610	932310
Calibrator Concentrate	100895	932507
RGP Reagent	103159	N/A <sup>1</sup>

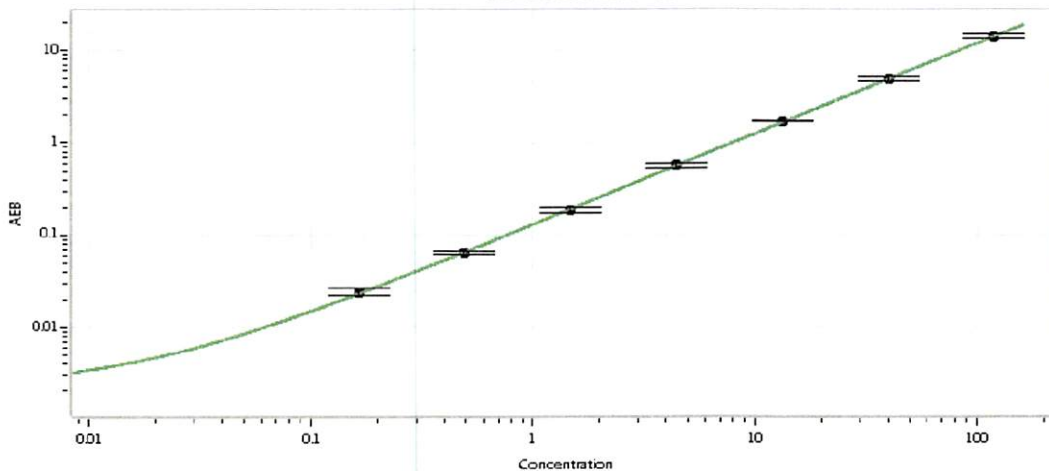
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	911406	2.55	1.27-2.97
Control 2	911406	147	80.9-189
Panel 1	829001	2.88	1.80-4.20
Panel 2	829002	5.95	0.938-7.09
Panel 3	829003	11.4	6.62-15.5
Panel 4	829004	52.5	41.1-61.6

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

Calibrator Concentrate Value Assignment (pg/mL)	3472
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.165
C	0.494
D	1.48
E	4.44
F	13.3
G	40.0
H	120

Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i>	<i>09 Dec 2019</i>
	Name	Title	Signature/Date	