

Product Number:	101622
Lot Number:	502134
Expiration:	19-Aug-2020
Platform(s):	HD-1/HD-x & SR-X



Data below represents results generated on the Simoa™ HD-X Analyzer.

Component	Part Number	Lot Number
Bead Reagent	101624	923803
Detector Reagent	101623	923110
SBG Reagent	101625	923204
Sample Diluent	103203	927721
Calibrator Diluent	101627	930819
Calibrator Concentrate	100652	906001
RGP Reagent	103159	N/A <sup>1</sup>

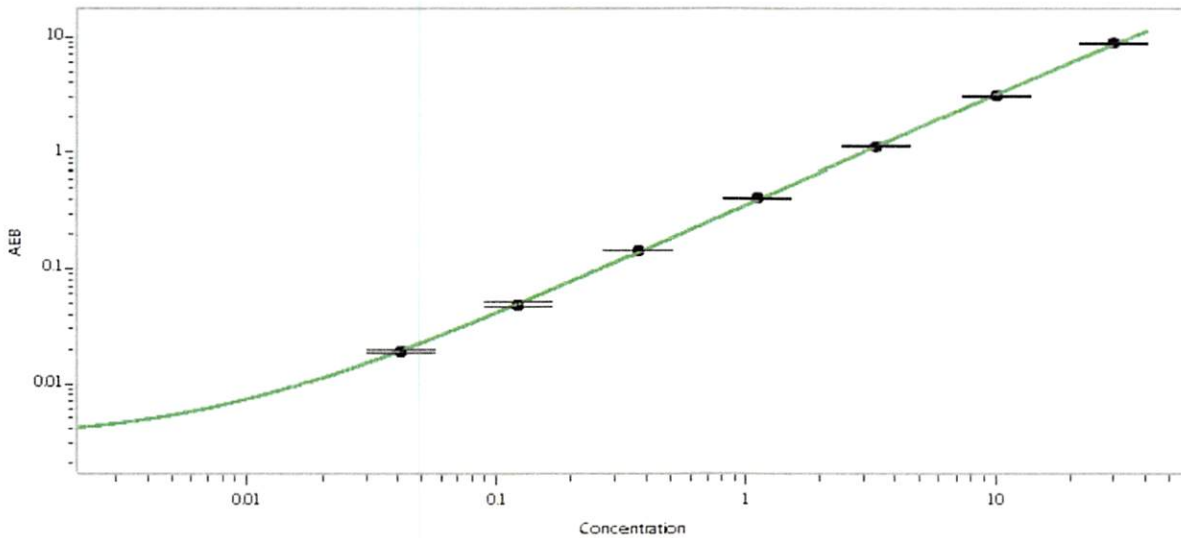
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	901702	1.29	0.944-1.42
Control 2	901702	21.3	15.4-23.1
Panel 1	MF-050918-01	55.7	41.6-67.5
Panel 2	MF-050918-02	8.37	5.03-8.84
Panel 3	MF-050918-03	1.15	0.644-1.24

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

Calibrator Concentrate Value Assignment (pg/mL)	864
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.041
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval: Patrizia Stadler Assoc. Director, QC P. Stadler 19 Dec 2019  
Name Title Signature/Date