

Product Number:	101599
Lot Number:	502194
Expiration:	13-Jan-2021
Platform(s):	HD-1/HD-X & SR-X



Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	101601	001305
Detector Reagent	102012	001314
SBG Reagent	101602	001315
Sample Diluent	101603	001313
Calibrator Diluent	102014	001312
Calibrator Concentrate	100778	001414
RGP Reagent	103159	N/A ¹

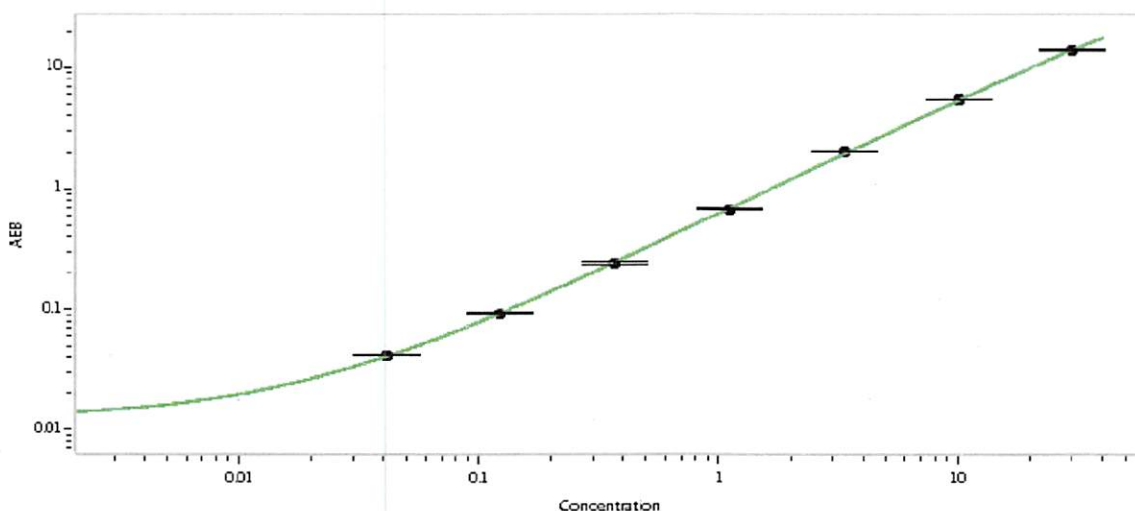
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-012518-02	0.358	0.352-0.537
Control 2	MF-012518-02	21.4	14.5-28.3
Panel 1	900901	1.00	0.723-1.90
Panel 3	900903	7.67	6.32-9.49
Panel 4	900904	28.5	22.3-33.4

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

Calibrator Concentrate Value Assignment (pg/mL)	706
---	-----

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* 31 Jan 2020
Name Title Signature/Date