

Product Number:	101599
Lot Number:	502288
Expiration:	26-Feb-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	008306
Detector Reagent	102012	007703
SBG Reagent	101602	007805
Sample Diluent	101603	007804
Calibrator Diluent	102014	007806
Calibrator Concentrate	100778	007810
RGP Reagent	103159	N/A <sup>1</sup>

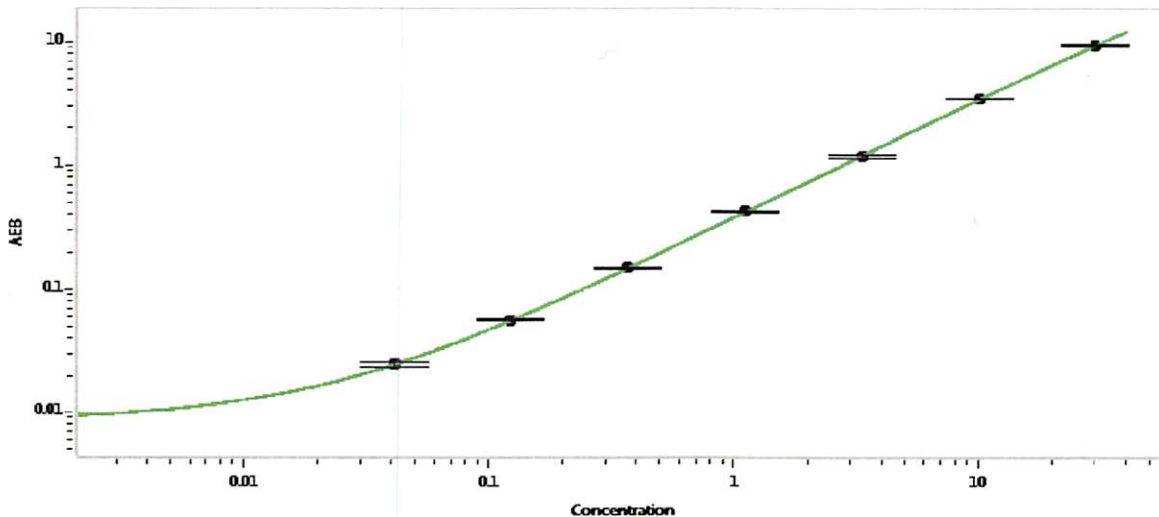
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-012518-02	0.383	0.352-0.537
Control 2	MF-012518-02	22.0	14.5-28.3
Panel 3	900903	8.01	6.49-9.73
Panel 4	900904	32.9	22.3-35.1

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

Calibrator Concentrate Value Assignment (pg/mL)	825
---	-----

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 Mar 2020

Name

Title

Signature/Date