

Product Number:	101599
Lot Number:	501675
Expiration:	13-Feb-2020
Platform(s):	HD-1 & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	101601	905208
Detector Reagent	102012	905106
SBG Reagent	101602	905107
Sample Diluent	101603	905105
Calibrator Diluent	102014	904413
Calibrator Concentrate	100778	904508
RGP Reagent	103159	N/A <sup>1</sup>

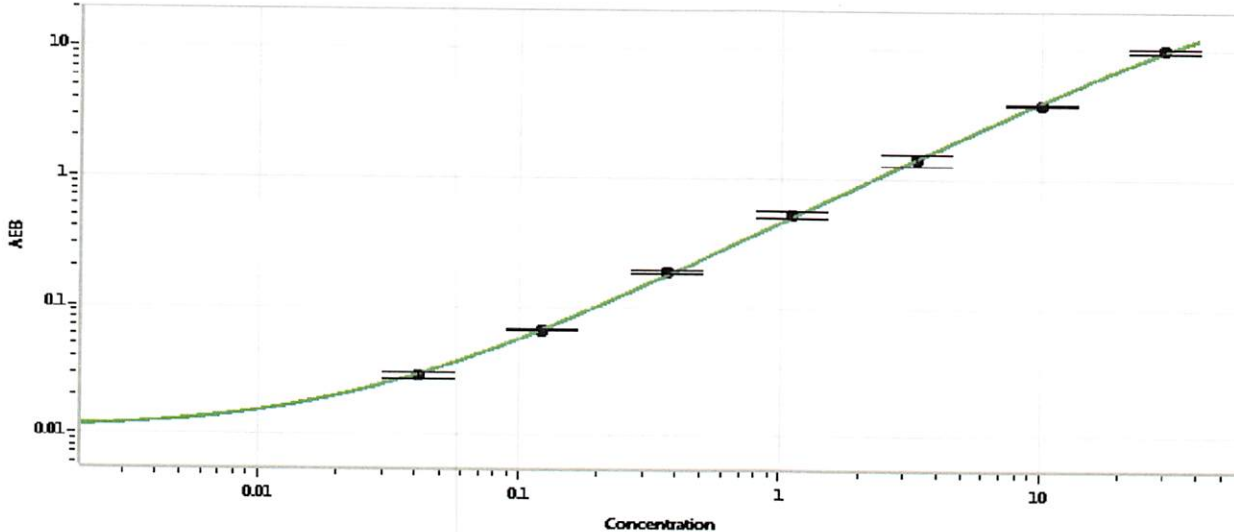
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-012518-02	0.376	0.363-0.545
Control 2	MF-012518-02	18.4	15.1-23.9
Panel 1	900901	1.25	0.963-1.97
Panel 2	900902	1.12	1.08-2.40
Panel 3	900903	7.72	6.39-6.58
Panel 4	900904	32.2	21.9-33.2

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

Calibrator Concentrate Value Assignment (pg/mL)	1006
---	------

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*

07/16/2019

Name

Title

Signature/Date