

Product Number:	101599
Lot Number:	501614
Expiration:	10-Jan-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	901507
Detector Reagent	102012	902502
SBG Reagent	101602	901011
Sample Diluent	101603	901501
Calibrator Diluent	102014	901013
Calibrator Concentrate	100778	81562
RGP Reagent	103159	N/A ¹

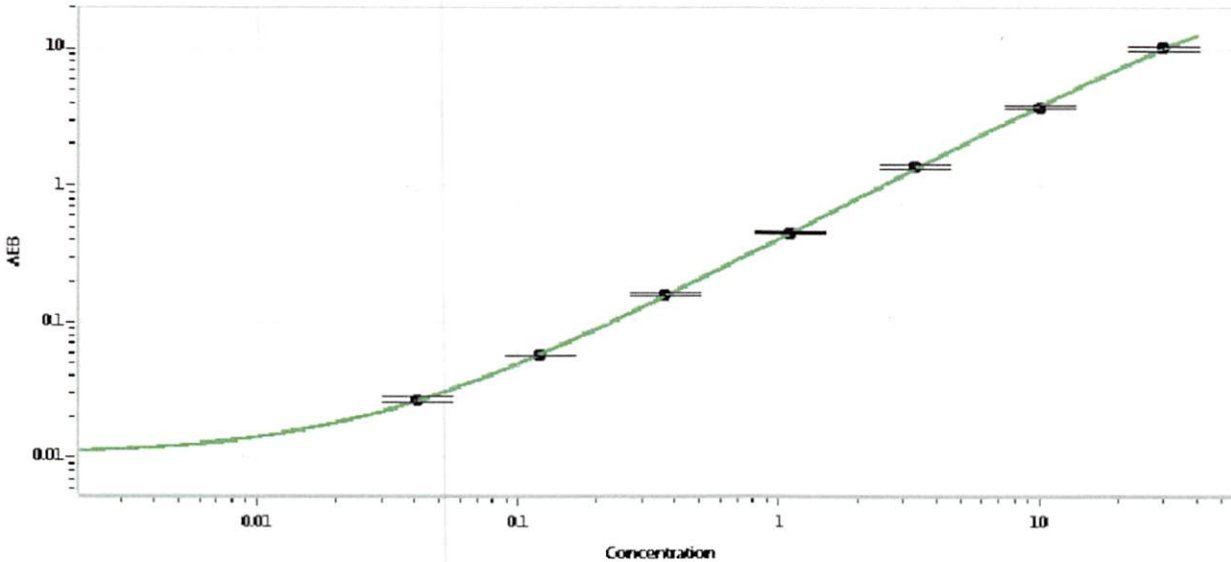
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-012518-02	0.448	0.363-0.545
Control 2	MF-012518-02	23.4	14.7-24.7
Panel 1	MF-120115-02	0.838	0.608-1.27
Panel 2	MF-071615-02	7.32	5.58-8.37
Panel 3	MF-071615-03	2.18	1.68-2.52

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

Calibrator Concentrate Value Assignment (pg/mL)	800
---	-----

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
Name

QC Assoc. Director
Title

P. Stadler 31 Jan 2019
Signature/Date