

Product Number:	100196
Lot Number:	501648
Expiration:	29-Jun-2019
Platform(s):	HD-1



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

Component	Part Number	Lot Number
Bead Reagent	100391	825612
Detector Reagent	100315	825513
SBG Reagent	100316	825514
Sample Diluent	100317	818011
Calibrator Diluent	100665	825515
Calibrator Concentrate	100664	72701
RGP Reagent	103159	N/A <sup>1</sup>

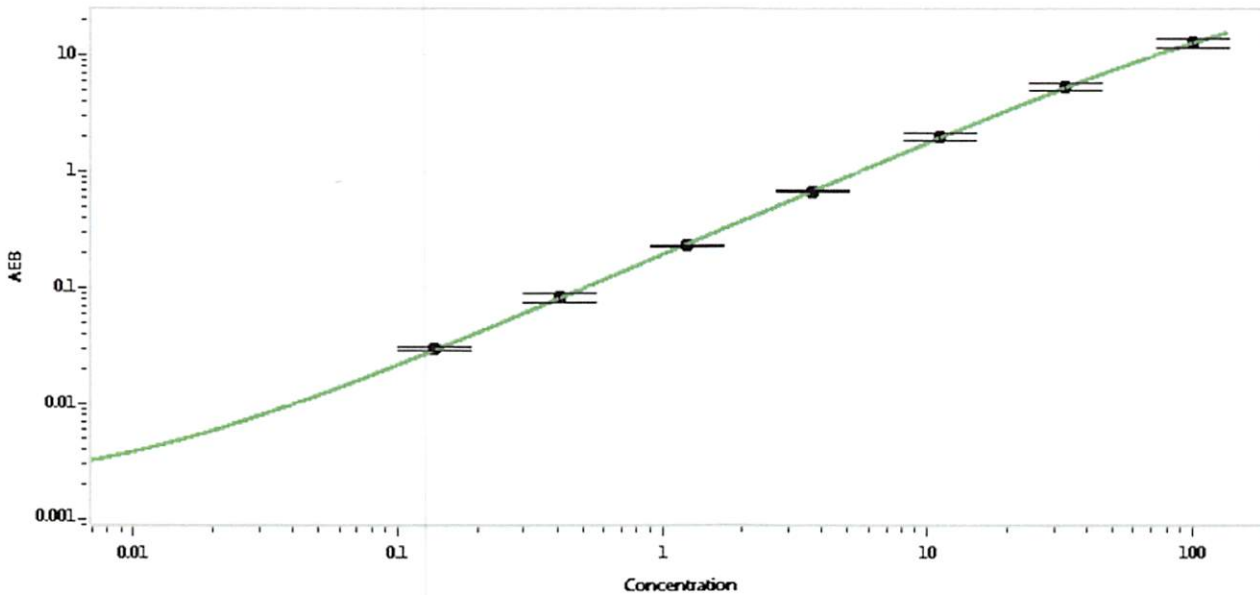
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	52521	97.4	75.6-118
Control 2	52521	9.07	7.72-11.6
Panel 1	MF-012116-05	1.02	0.871-1.73
Panel 2	MF-090815-18	6.73	4.45-9.13
Panel 3	MF-090815-19	83.6	53.9-94.9

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

Calibrator Concentrate Value Assignment (pg/mL)	2154
---	------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.137
C	0.412
D	1.23
E	3.70
F	11.1
G	33.3
H	100

Review/Approval

Patrizia Stadler QC Assoc. Director

*P. Stadler* 08 Feb 2019

Name

Title

Signature/Date