

Product Number:	100196
Lot Number:	502115
Expiration:	4-Oct-2020
Platform(s):	HD-1/HD-X



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

Component	Part Number	Lot Number
Bead Reagent	100391	931906
Detector Reagent	100315	931803
SBG Reagent	100316	931804
Sample Diluent	100317	927716
Calibrator Diluent	100665	931909
Calibrator Concentrate	100664	931810
RGP Reagent	103159	N/A ¹

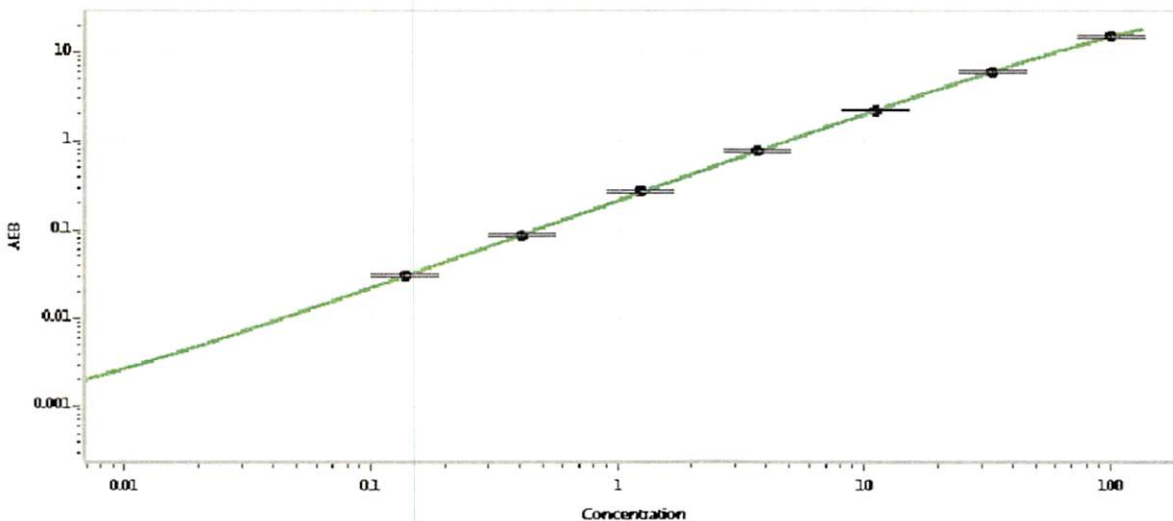
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	910002	8.01	7.14-10.7
Control 2	910002	87.9	73.6-110
Panel 1a	MF-062018-01	0.285	0.041-1.22
Panel 2	MF-062018-02	7.53	6.73-11.0
Panel 3	MF-062018-03	87.5	72.9-152

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

Calibrator Concentrate Value Assignment (pg/mL)	2096
---	------

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
Name

Assoc. Director, QC
Title

P. Stadler 09 Dec 2019
Signature/Date