

Product Number:	100196
Lot Number:	502250
Expiration:	14-Nov-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	005210
Detector Reagent	100315	005101
SBG Reagent	100316	004907
Sample Diluent	100317	933818
Calibrator Diluent	100665	931822
Calibrator Concentrate	100664	005203
RGP Reagent	103159	N/A ¹

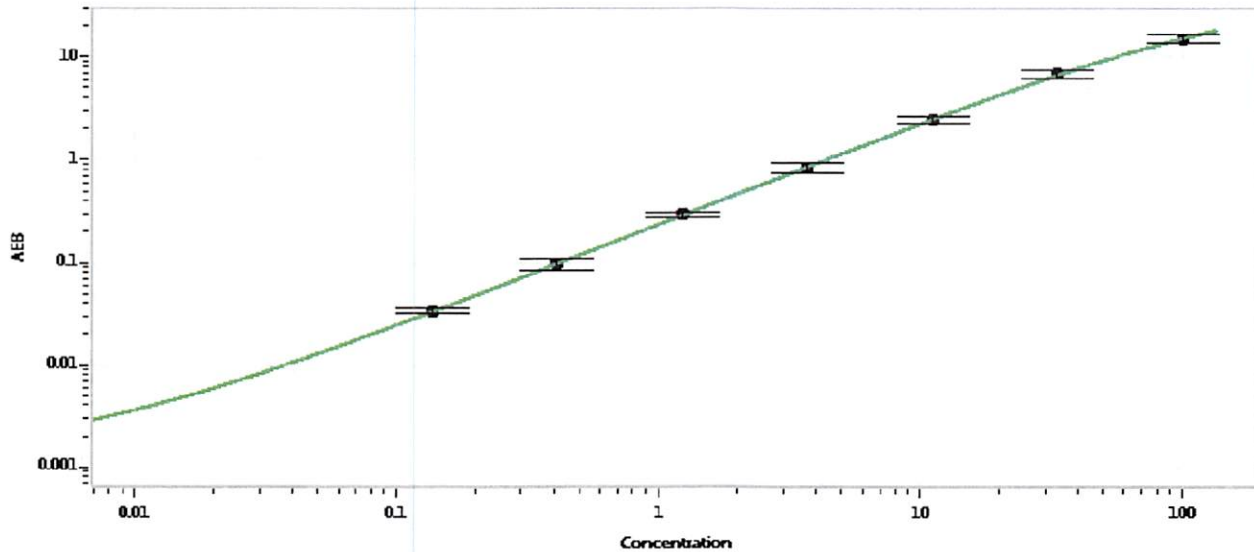
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	910002	9.40	7.19-10.8
Control 2	910002	100	74.4-112
Panel 1a	MF-062018-01	0.749	0.097-1.20
Panel 2	MF-062018-02	9.12	6.90-10.9
Panel 3	MF-062018-03	120	72.9-155

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

Calibrator Concentrate Value Assignment (pg/mL)	1747
---	------

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

Assoc. Director, QC

P. Stadler

28 Feb 2020

Name

Title

Signature/Date