

Product Number:	103337
Lot Number:	501789
Expiration:	29-Mar-2020
Platform(s):	HD-1 & SR-X



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

Component	Item Number	Lot Number
Bead Reagent	103334	911906
Detector Reagent	101091	909811
SBG Reagent	101101	909812
Sample Diluent	103336	909813
Calibrator Diluent	101095	908817
Calibrator Concentrate	101093	906801
RGP Reagent	103159	N/A ¹

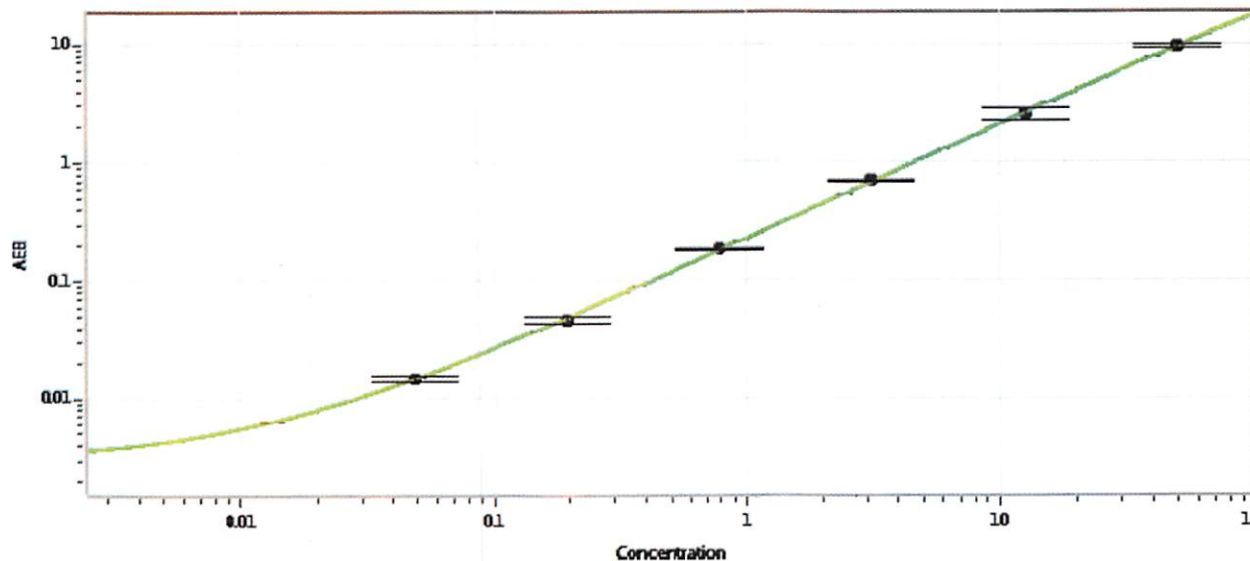
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-033018-02	21.1	12.6-22.1
Control 2	MF-033018-02	209	128-218
Panel 1	MF-033018-03	3.75	2.51-3.76
Panel 2	MF-033018-04	6.69	4.20-6.96
Panel 3	MF-033018-05	63.0	44.1-66.1

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

Calibrator Concentrate Value Assignment (pg/mL)	1,325
---	-------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.049
C	0.195
D	0.781
E	3.13
F	12.5
G	50.0

Review/Approval: Patrizia Stadler QC Assoc. Director
 Name: Patrizia Stadler Title: QC Assoc. Director Signature/Date: P. Stadler 06 MAY 2019