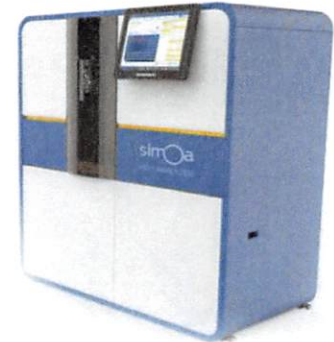


Product Number:	101599
Lot Number:	501988
Expiration:	10-Jan-2020
Platform(s):	HD-1/HD-X & SR-X



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	904508
RGP Reagent	103159	N/A ¹

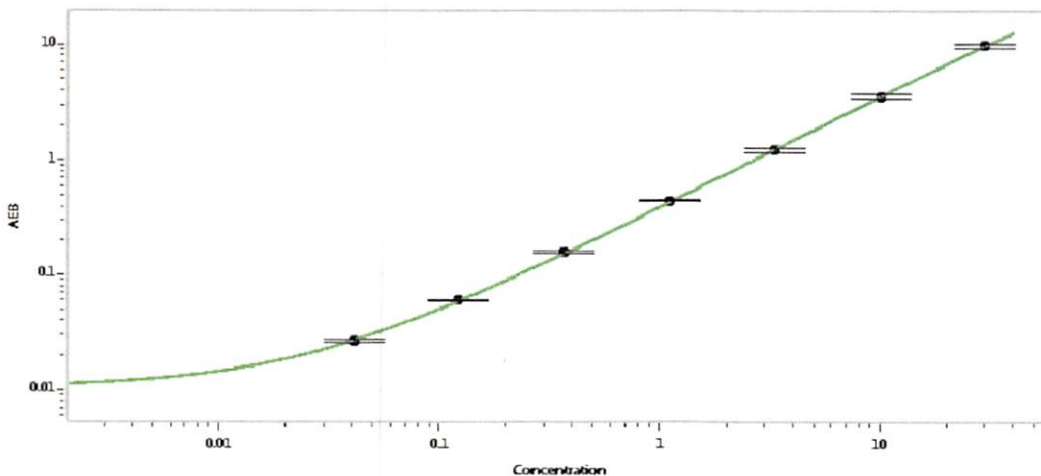
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-012518-02	0.438	0.365-0.548
Control 2	MF-012518-02	23.8	13.9-27.2
Panel 1	900901	1.29	0.392-3.05
Panel 2	900902	1.68	1.11-2.32
Panel 3	900903	8.59	6.42-9.63
Panel 4	900904	31.9	21.8-34.5

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

Calibrator Concentrate Value Assignment (pg/mL)	1006
---	------

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

Daniel Ritchie Senior QC Analyst

[Handwritten Signature] / 23 Aug 2019

Name

Title

Signature/Date

*Signing on behalf of Patrizia Stadler