

Product Number:	100200
Lot Number:	501649
Expiration:	12-Jun-2019
Platform(s):	HD-1 & SR-X



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

Component	Item Number	Lot Number
Bead Reagent	101089	820605
Detector Reagent	101091	821316
SBG Reagent	101101	819112
Sample Diluent	101099	818005
Calibrator Diluent	101095	816308
Calibrator Concentrate	101093	828403
RGP Reagent	103159	N/A ¹

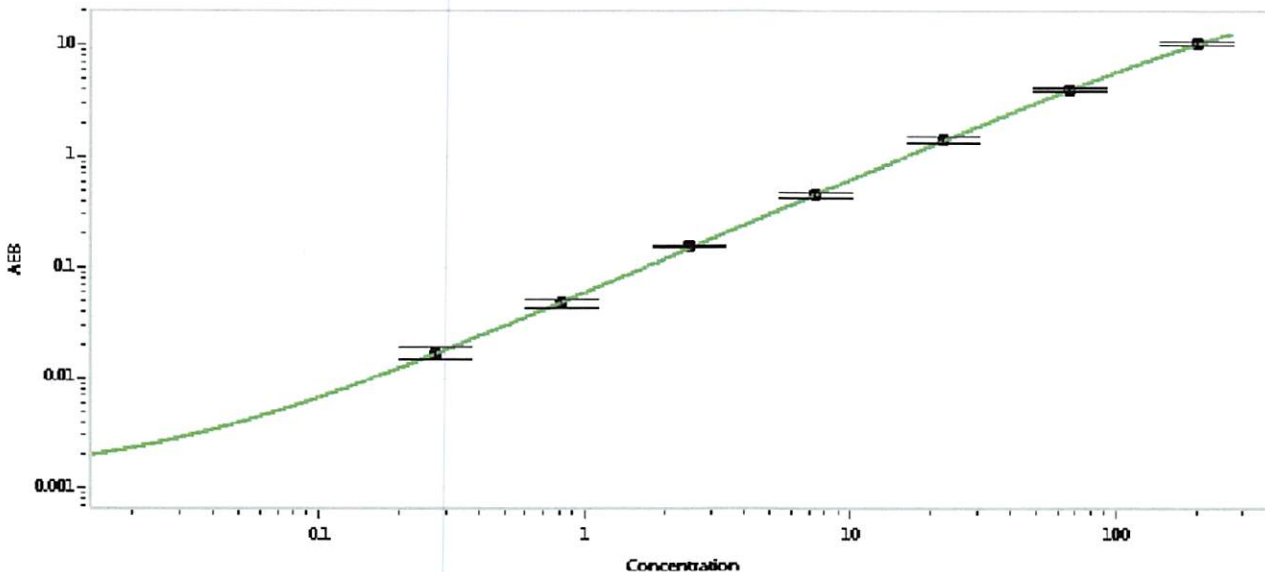
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-092215-03	8.19	6.99-11.4
Control 2	MF-092215-03	75.2	67.3-112
Panel 1	MF-012116-06	1.75	1.58-2.80
Panel 2	MF-093015-02	3.02	1.45-5.96
Panel 3	MF-093015-03	31.3	24.1-59.5

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

Calibrator Concentrate Value Assignment (pg/mL)	1,751
---	-------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.274
C	0.823
D	2.47
E	7.41
F	22.2
G	66.7
H	200

Review/Approval	Patrizia Stadler QC Assoc. Director	<i>P. Stadler</i> 08 Feb 2019
	Name Title	Signature/Date