

Product Number:	101664
Lot Number:	501927
Expiration:	4-Apr-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	101666	909409
Detector Reagent	101665	910510
SBG Reagent	101667	909816
Sample Diluent	101668	910107
Calibrator Diluent	101669	910106
Calibrator Concentrate	103217	912801
RGP Reagent	103159	N/A ¹

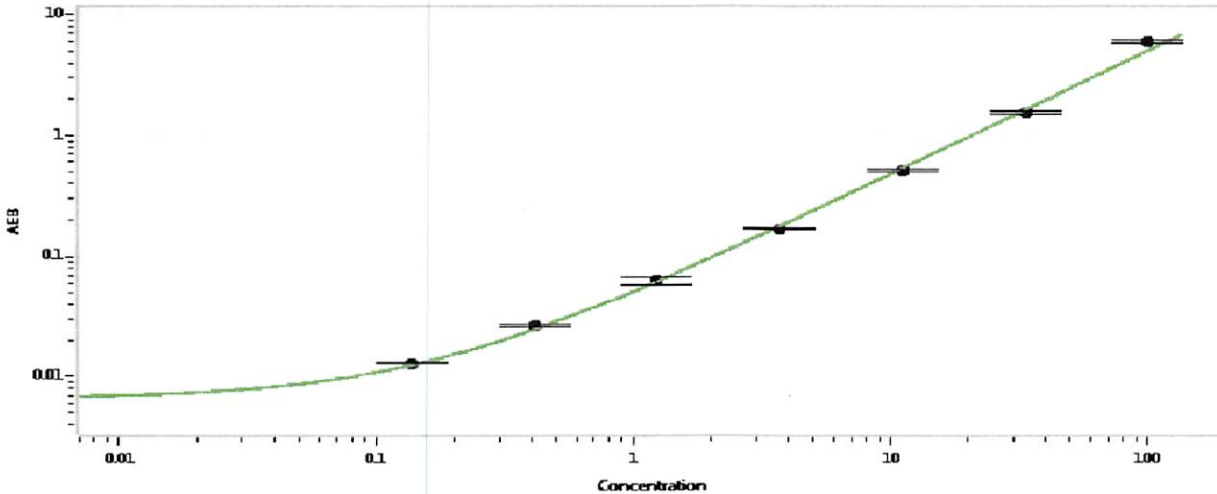
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	52520	0.982	0.740-1.33
Control 2	52520	34.2	27.1-40.6
Panel 1	MF-082015-01	1.02	0.942-1.53
Panel 2	MF-081415-02	8.55	7.46-11.2
Panel 3	MF-081415-03	10.1	9.20-15.0
Panel 4	MF-090315-09	42.5	27.6-66.8

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

Calibrator Concentrate Value Assignment (pg/mL)	2862
---	------

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. Stodley
Signature/Date

18 Sep 2019