

Product Number:	101605
Lot Number:	501700
Expiration:	27-Feb-2020
Platform(s):	HD-1



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

Component	Part Number	Lot Number
Bead Reagent	101607	905810
Detector Reagent	101606	906403
SBG Reagent	101608	906304
Sample Diluent	103200	906309
Calibrator Diluent	101610	906014
Calibrator Concentrate	100895	901403
RGP Reagent	103159	N/A ¹

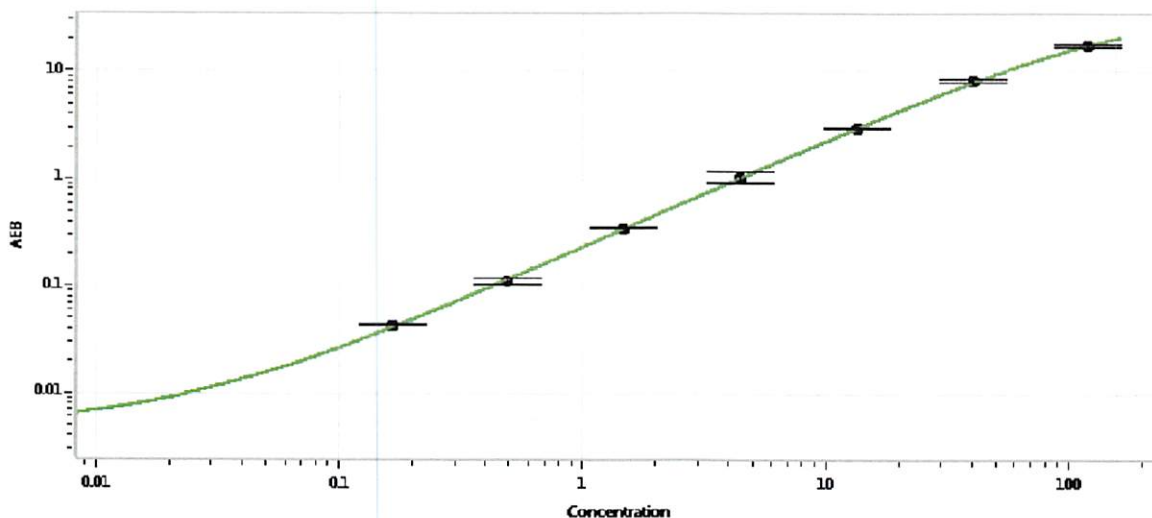
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-112515-02	5.80	3.13-7.30
Control 2	MF-112515-02	200	93.3-218
Panel 1	829001	2.87	1.85-4.32
Panel 2	829002	2.37	2.17-5.07
Panel 3	829003	9.74	6.71-15.7
Panel 4	829004	55.6	40.7-60.1

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

Calibrator Concentrate Value Assignment (pg/mL)	4,722
---	-------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.165
C	0.494
D	1.48
E	4.44
F	13.3
G	40.0
H	120

Review/Approval	Patrizia Stadler	Associate Director, QC	<i>P. Stadler</i>	08 Nov 2019
	Name	Title	Signature/Date	