

Product Number:	101580
Lot Number:	501935
Expiration:	14-Mar-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	101582	908803
Detector Reagent	101581	908510
SBG Reagent	101583	908407
Sample Diluent	103201	907309
Calibrator Diluent	101585	909205
Calibrator Concentrate	100655	909204
RGP Reagent	103159	N/A ¹

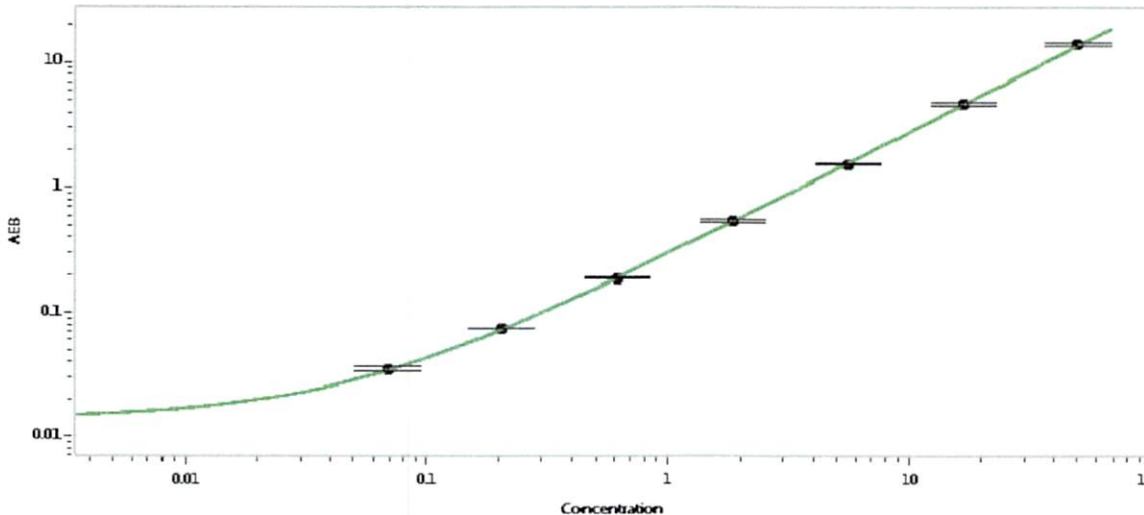
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-042817-01	2.90	2.05-3.08
Control 2	MF-051617-02	44.9	35.8-53.7
Panel 1	MF-051017-03	30.6	20.9-32.4
Panel 2	MF-032217-11	2.03	1.33-2.15
Panel 3	MF-033018-06	3.21	1.68-3.50

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

Calibrator Concentrate Value Assignment (pg/mL)	1065 pg/mL
---	------------

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.069
C	0.206
D	0.617
E	1.85
F	5.56
G	16.7
H	50.0

Review/Approval

Patrizia Stadler Assoc. Director, QC

p. stadler 25 Jul 2019

Name

Title

Signature/Date