

Product Number:	101580
Lot Number:	502141
Expiration:	4-Oct-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	932604
Detector Reagent	101581	932606
SBG Reagent	101583	932618
Sample Diluent	103201	927723
Calibrator Diluent	101585	932616
Calibrator Concentrate	100655	932607
RGP Reagent	103159	N/A ¹

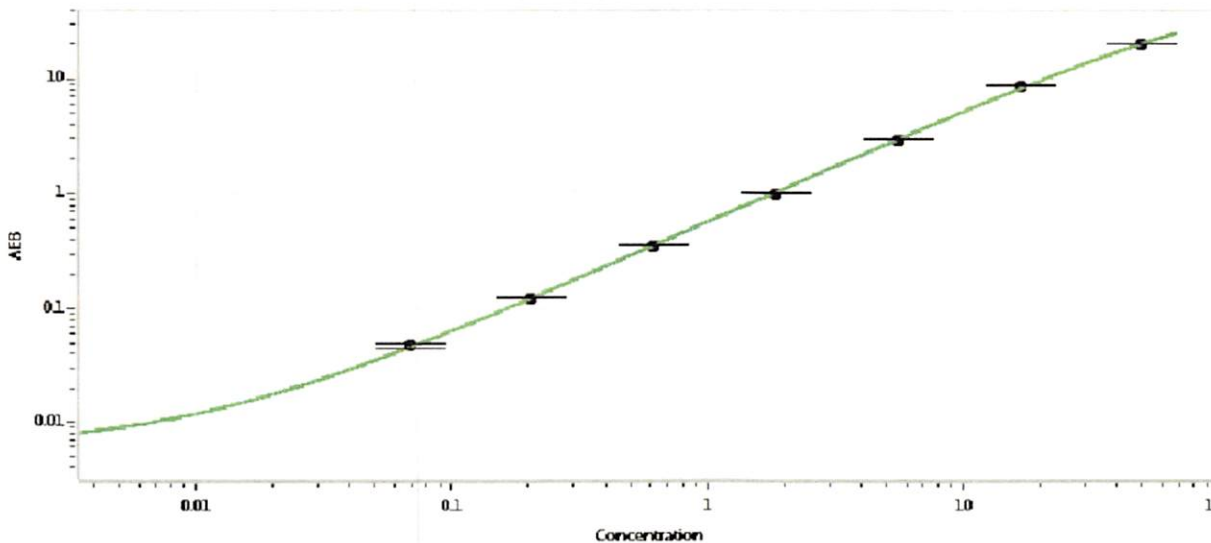
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-042817-01	2.60	2.05-3.08
Control 2	MF-051617-02	47.0	35.8-53.7
Panel 1	MF-051017-03	24.8	20.9-32.4
Panel 2	MF-032217-11	1.61	1.33-2.15
Panel 3	MF-033018-06	2.64	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)	1311 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
Name Title

P. Stadler 05 Dec 2019
Signature/Date