

Product Number: 101580
 Lot Number: 501639
 Expiration: 29-Jan-2020
 Platform(s): HD-1 & SR-X



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

Component	Part Number	Lot Number
Bead Reagent	101582	903604
Detector Reagent	101581	903605
SBG Reagent	101583	903606
Sample Diluent	101584	902925
Calibrator Diluent	101585	902909
Calibrator Concentrate	100655	834101
RGP Reagent	103159	N/A ¹

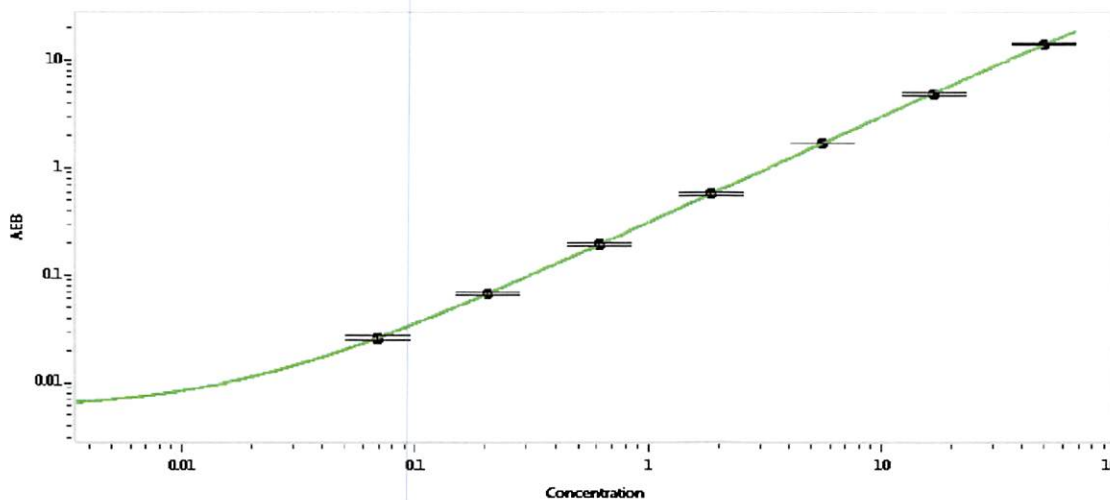
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-042817-01	2.88	2.05-3.08
Control 2	MF-051617-02	52.2	35.8-53.7
Panel 1	MF-051017-03	29.9	20.9-32.4
Panel 2	MF-032217-11	2.05	1.33-2.15
Panel 3	MF-033018-06	3.32	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) 1155 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: Patrizia Stadler Title: QC Signature/Date: P. Stadler 19 Feb 2019