

Product Number:	102329
Lot Number:	501565
Expiration:	19-Sep-2019
Platform(s):	HD-1



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

Component	Part Number	Lot Number
Bead Reagent	102322	835108
Detector Reagent	102324	900906
SBG Reagent	102325	900907
Sample Diluent	102326	826226
Calibrator Diluent	102327	834712
Calibrator Concentrate	100849	80923
RGP Reagent	103159	N/A ¹

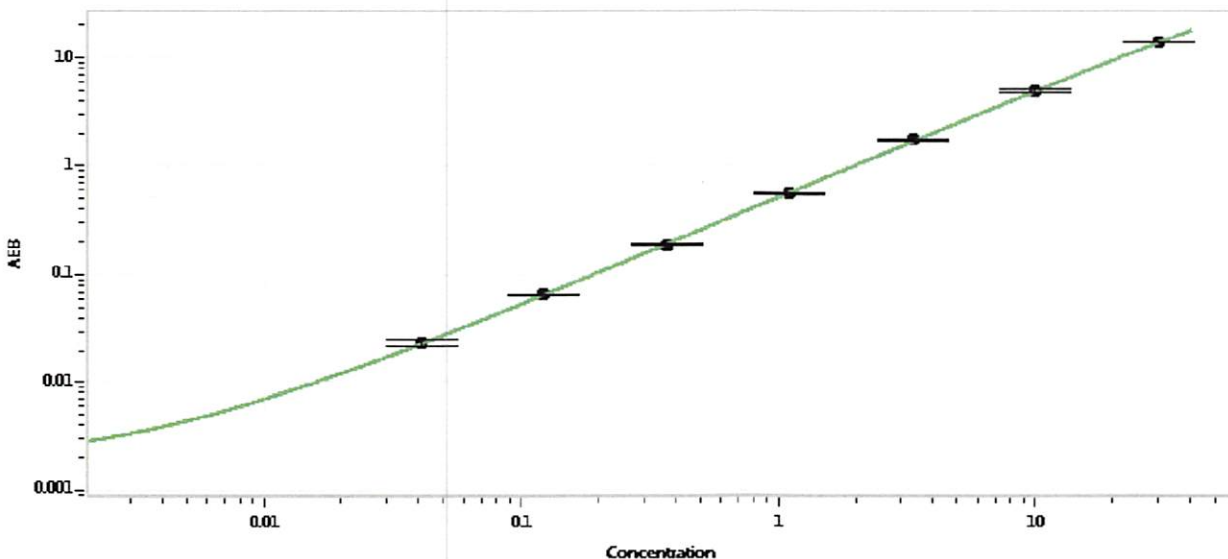
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-092716-01	0.142	0.108-0.300
Control 2	MF-092716-01	22.3	17.8-37.9
Panel 1	MF-092116-03	0.081	0.065-0.137
Panel 2	MF-092116-04	0.076	0.069-0.135
Panel 3	MF-092116-05	13.5	9.7-25.5

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

Calibrator Concentrate Value Assignment (pg/mL)	647
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.041
C	0.123
D	0.370
E	1.110
F	3.33
G	10.0
H	30.0

Review/Approval

Patrizia Stadler QC Assoc. Director

Name

Title

P. Stadler 25 Jan 2019

Signature/Date