

Product Number:	102329
Lot Number:	501200
Expiration:	30-Sep-2018
Platform(s):	HD-1



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

Component	Part Number	Lot Number
Bead Reagent	102322	80930
Detector Reagent	102324	80891
SBG Reagent	102325	80892
Sample Diluent	102326	80863
Calibrator Diluent	102327	80475
Calibrator Concentrate	100849	80923
RGP Reagent	103159	N/A ¹

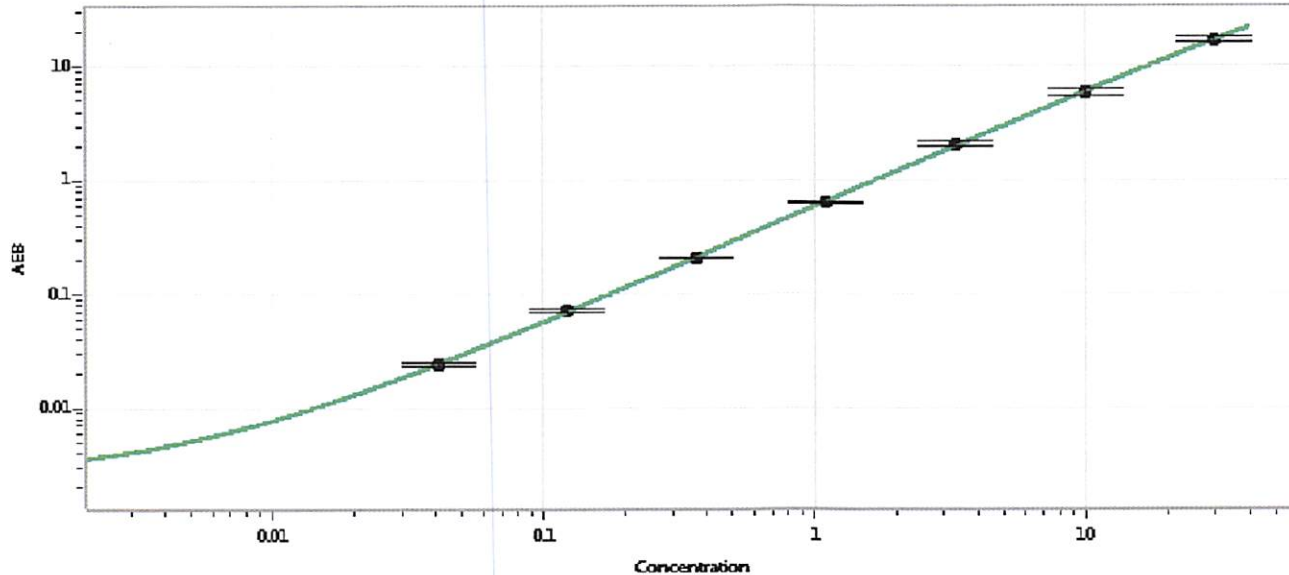
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-092716-01	0.214	0.150-0.287
Control 2	MF-092716-01	25.7	19.1-37.9
Panel 1	MF-092116-03	0.099	0.065-0.137
Panel 2	MF-092116-04	0.099	0.069-0.135
Panel 3	MF-092116-05	13.8	10.3-25.8

* 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

Sivasankari Murugan
Name

QC Supervisor
Title

[Signature] 14 APR 2018
Signature/Date