

Product Number:	100988
Lot Number:	501216
Expiration:	19-Sep-2018
Platform(s):	HD-1



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

Component	Part Number	Lot Number
Bead Reagent	100974	80790
Detector Reagent	100976	81030
SBG Reagent	100987	80792
Sample Diluent	100985	80655
Calibrator Diluent	100980	80783
Calibrator Concentrate	100978	71812
RGP Reagent	103159	N/A ¹

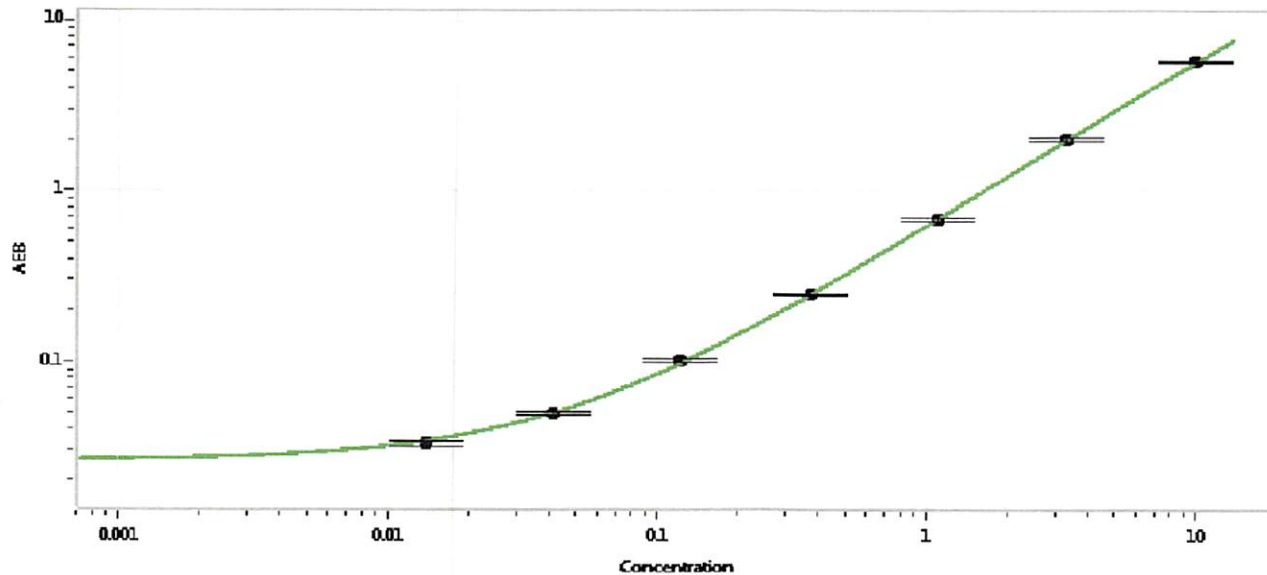
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	MF-101816-02	0.153	0.053-0.167
Control 2	MF-101816-02	8.64	6.14-10.2
Panel 1	MF-101716-04	0.292	0.211-0.343
Panel 2	MF-101716-05	0.073	0.030-0.095
Panel 3	MF-101816-03	17.6	15.0-22.5

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

Calibrator Concentrate Value Assignment (pg/mL)	285
---	-----

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.014
C	0.041
D	0.123
E	0.370
F	1.11
G	3.33
H	10.0

Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

S. Murugan
Signature/Date
24 APR 2018