

Product Number:	101599
Lot Number:	501276
Expiration:	1-Mar-2019
Platform(s):	HD-1 & SR-X



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

Component	Item Number	Lot Number
Bead Reagent	101601	81520
Detector Reagent	102012	81551
SBG Reagent	101602	81550
Sample Diluent	101603	81521
Calibrator Diluent	102014	81003
Calibrator Concentrate	100778	81562
RGP Reagent	103159	N/A ¹

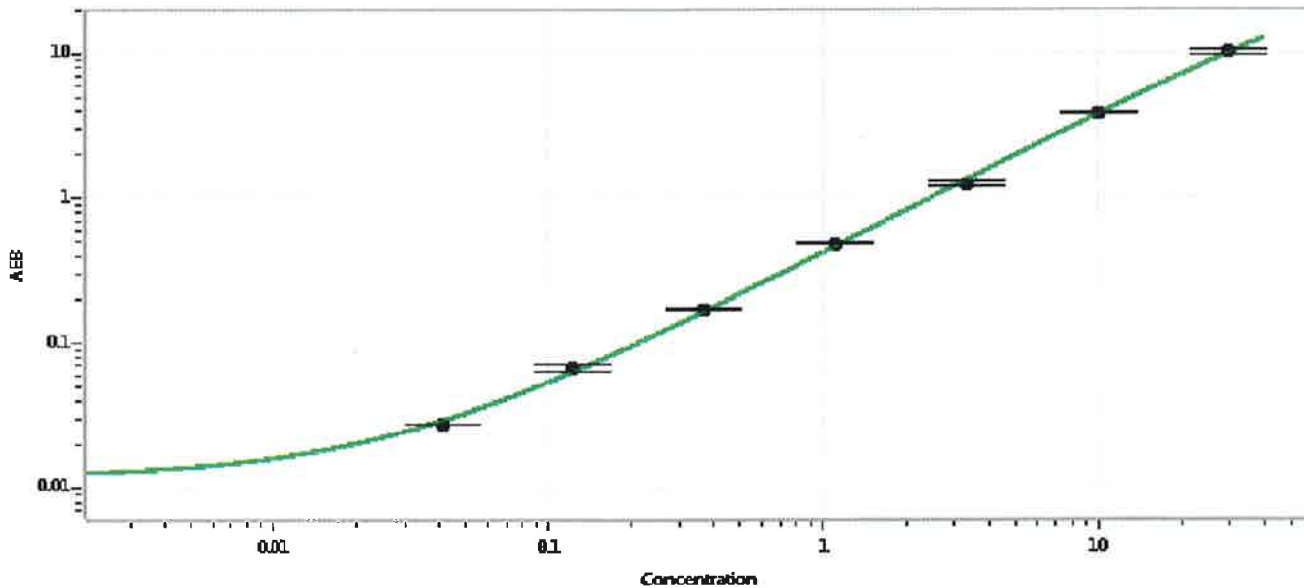
¹ RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	40582	0.535	0.364-0.646
Control 2	40582	27.8	19.2-33.0
Panel 1	MF-071615-01	0.336	0.325-0.513
Panel 2	MF-071615-02	6.63	5.58-8.37
Panel 3	MF-071615-03	2.03	1.68-2.52

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

Calibrator Concentrate Value Assignment (pg/mL)	800
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.041
C	0.123
D	0.370
E	1.11
F	3.33
G	10.0
H	30.0

Review/Approval: Sivasankari Murugan Name QC Supervisor Title [Signature] Signature/Date 16/02/2018 Date