

Product Number:	101599
Lot Number:	501136
Expiration:	8-Oct-2018
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	80450
Detector Reagent	102012	80440
SBG Reagent	101602	80471
Sample Diluent	101603	80434
Calibrator Diluent	102014	80088
Calibrator Concentrate	100778	73457
RGP Reagent	103159	N/A <sup>1</sup>

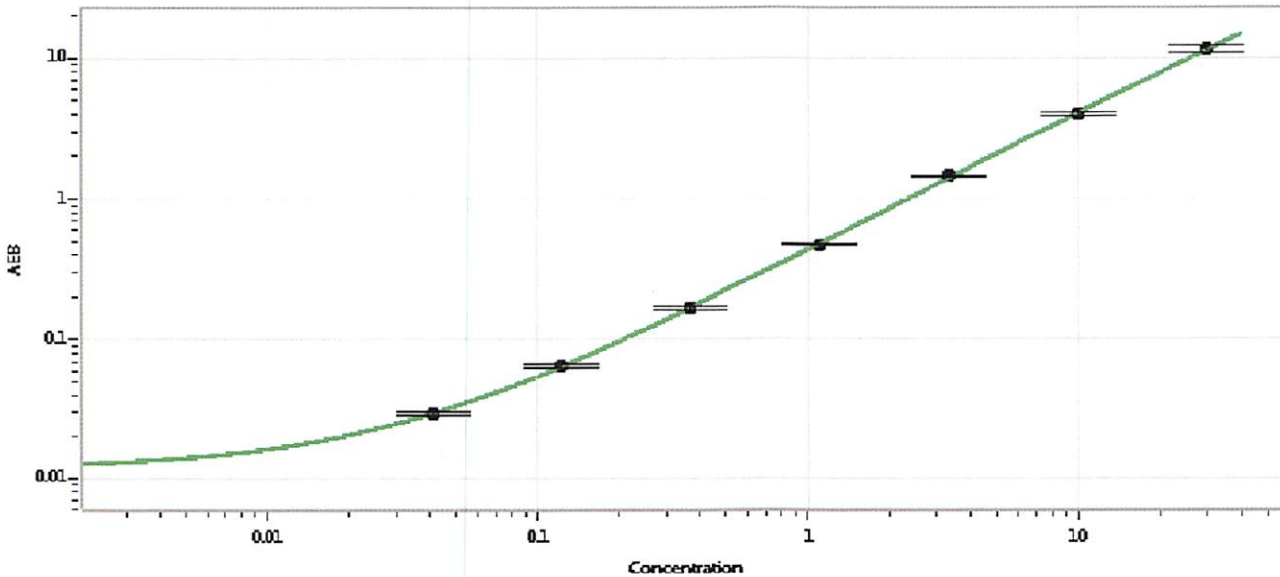
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	52442	0.617	0.485-0.727
Control 2	52442	29.2	19.5-34.7
Panel 1	MF-071615-01	0.341	0.325-0.513
Panel 2	MF-071615-02	7.71	5.58-8.37
Panel 3	MF-071615-03	2.32	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)	852
---	-----

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), 22 FEB 2018 (Date)