

Product Number:	103186
Lot Number:	501258
Expiration:	27-Dec-2018
Platform(s):	HD-1 & SR-X



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

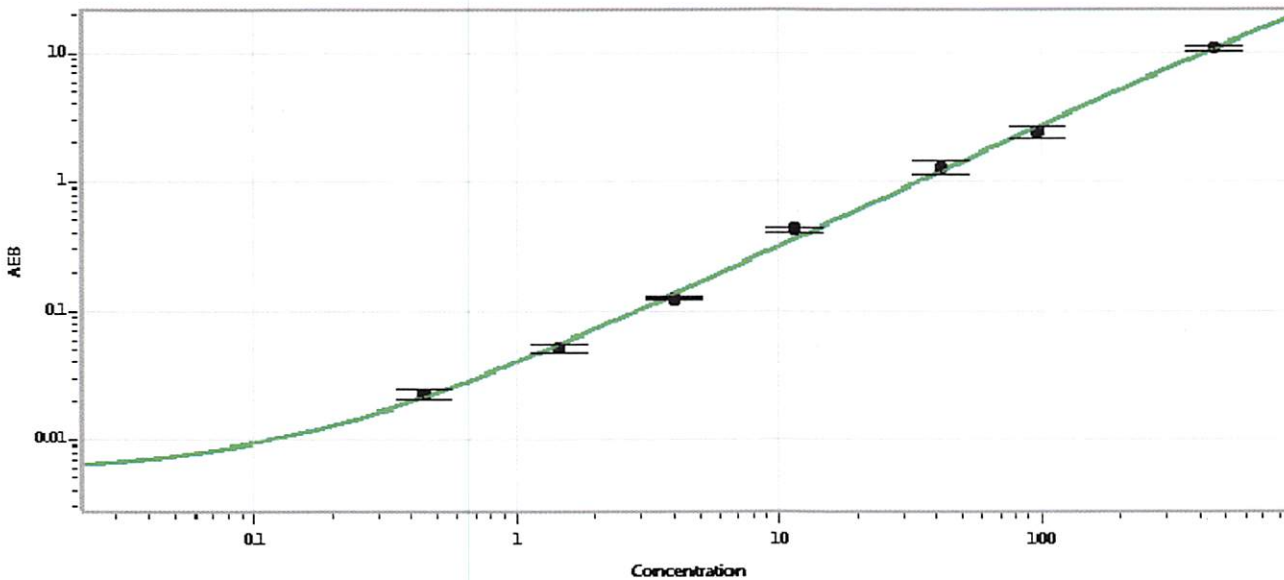
Component	Part Number	Lot Number
Bead Reagent	102246	81000
Detector Reagent	102248	80862
SBG Reagent	102250	80861
Sample Diluent	102252	80860
Reference Calibrators	102255	80942/81240
Controls	102256 - 102257	81090
RGP Reagent	103159	N/A <sup>1</sup>

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

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	81090	5.01	3.30-6.99
Control 2	81090	182	123-267
Panel 1	MF-092717-01	4.8	4.27-7.42
Panel 2	MF-092717-02	60.1	46.8-78.1
Panel 3	MF-092717-03	225	124-291

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

Calibration Curve



Calibrator Levels (pg/mL)

A	0.000
B	0.442
C	1.45
D	4.00
E	11.6
F	41.8
G	97.5
H	455

Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

*[Signature]*  
Signature/Date  
24 May 2018