

Product Number:	102860
Lot Number:	501329
Expiration:	14-Nov-2018
Platform(s):	HD-1 & SR-X



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

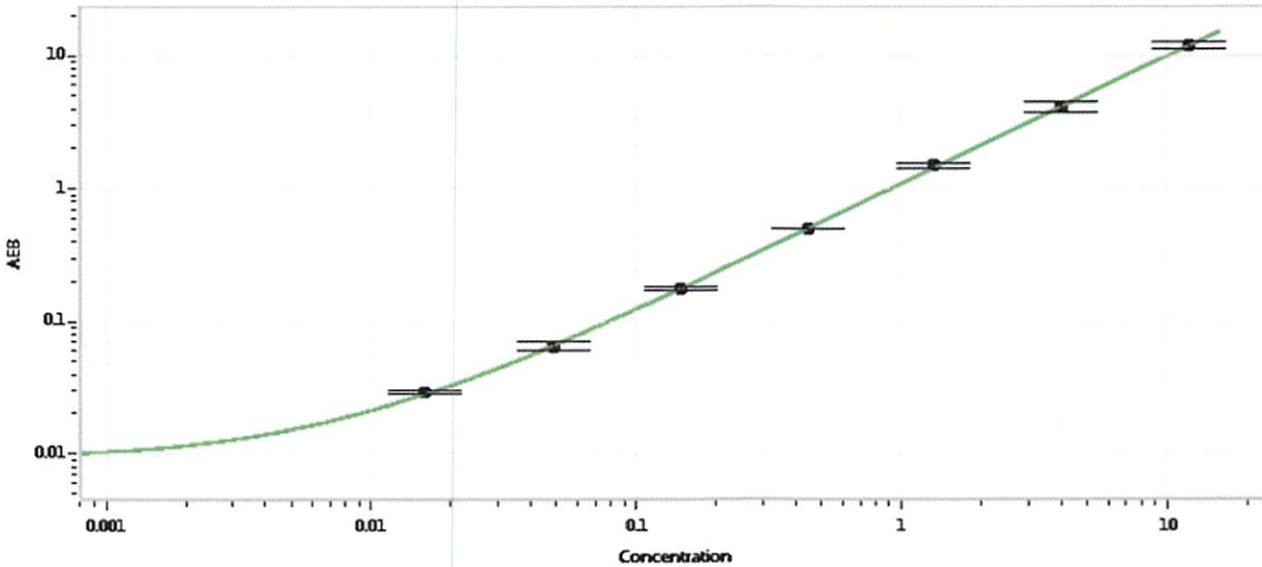
Component	Part Number	Lot Number
Bead Reagent	102852	81590
Detector Reagent	102854	81410
SBG Reagent	102855	81412
Sample Diluent	102856	81341
Calibrator Diluent	102859	81342
Calibrator Concentrate	102857	81423
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials*				* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.
	Lot	Result (pg/mL)	Mean Range (pg/mL)	
Control 1	MF-110817-02	0.289	0.251-0.377	
Control 2	MF-110817-02	4.76	4.19-6.47	
Panel 1	MF-100517-01	0.582	0.480-0.721	
Panel 2	MF-100517-02	1.30	1.12-1.69	
Panel 3	MF-100517-03	4.49	4.01-6.29	

Calibrator Concentrate Value Assignment (pg/mL)	217
---	-----

Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.016
C	0.049
D	0.148
E	0.444
F	1.33
G	4.00
H	12.0

Review/Approval: Sivasankari Murugan, QC Supervisor. Signature: [Handwritten Signature], Date: 07/09/2018.