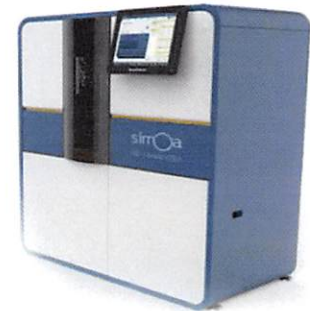


Product Number:	101160
Lot Number:	501198
Expiration:	16-Aug-2018
Platform(s):	HD-1



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

Component	Part Number	Lot Number
Bead Reagent†	101263	80931
Detector Reagent	101264	80990
SBG Reagent	101296	80761
Sample Diluent	101308	80472
Calibrators	101299	80880
Controls	101302 & 101316	80881
RGP Reagent	103159	N/A <sup>1</sup>

† Follow the Simoa™ Cytokine 3-Plex A Advantage Kit Instructions (KI-0011) to prepare the Bead Reagent.

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

**Release Materials**

**IL-6**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	80881	0.954	0.717-1.21
Control 2	80881	25.9	17.8-30.8
Panel 1	MF-072115-01	46.5	29.3-57.3
Panel 2	MF-072115-02	3.04	0.804-3.44
Panel 3	MF-072115-03	1.14	0.879-1.50
Panel 4	MF-072115-04	0.880	0.669-1.13

**IL-10**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	80881	0.656	0.418-0.692
Control 2	80881	13.2	9.44-14.5
Panel 1	MF-072115-01	1.19	0.828-1.28
Panel 2	MF-072115-02	11.6	6.21-14.2
Panel 3	MF-072115-03	1.97	1.47-2.21
Panel 4	MF-072115-04	2.13	1.54-2.31

**TNFα**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	80881	1.93	1.37-2.30
Control 2	80881	47.6	34.5-56.3
Panel 1	MF-072115-01	1.60	0.658-2.39
Panel 2	MF-072115-02	1.62	1.10-2.16
Panel 3	MF-072115-03	18.3	7.87-34.9
Panel 4	MF-072115-04	6.42	5.73-10.1

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

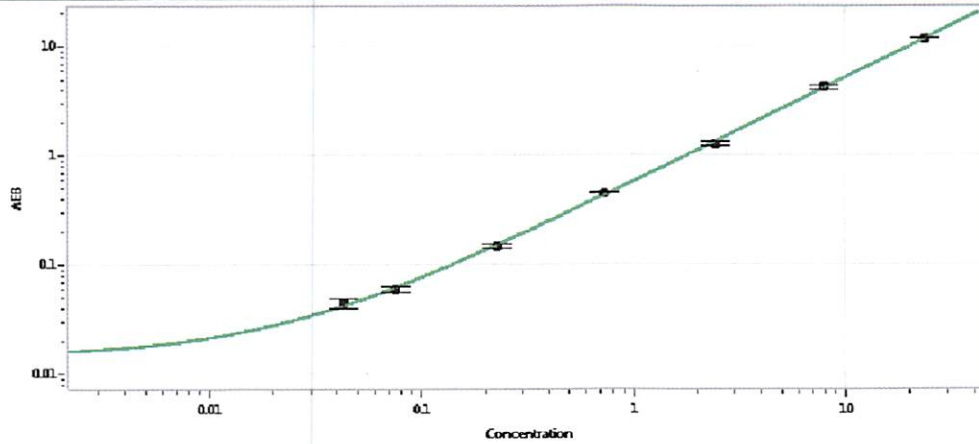
Review/Approval

Sivasankari Murugan  
Name

QC Supervisor  
Title

*[Signature]* 16 APR 2018  
Signature/Date

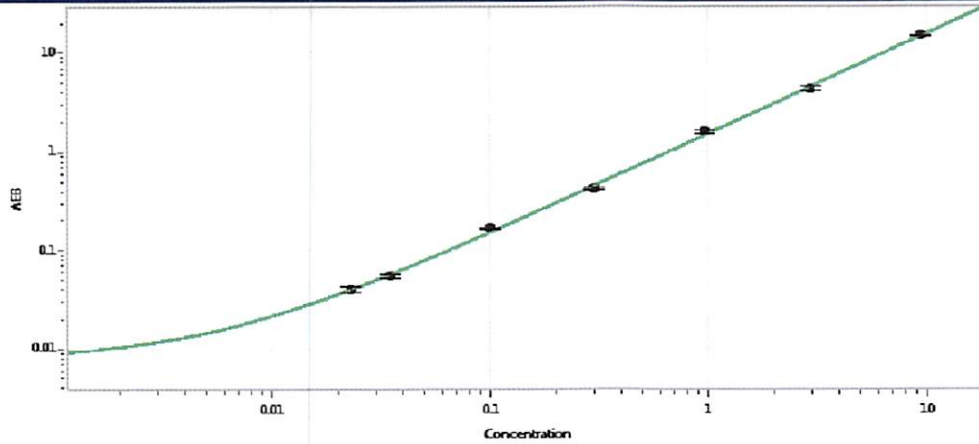
**IL-6 Calibration Curve**



**Calibrator Levels (pg/mL)**

A	0.000
B	0.043
C	0.076
D	0.229
E	0.729
F	2.44
G	7.99
H	23.6

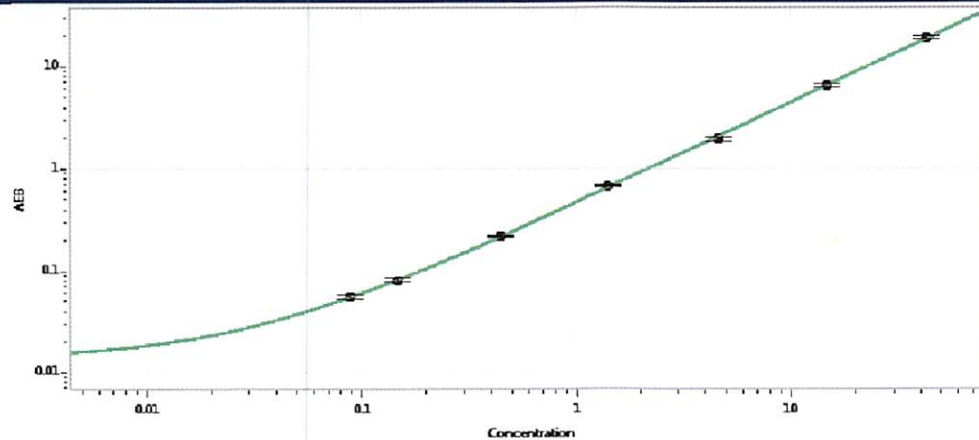
**IL-10 Calibration Curve**



**Calibrator Levels (pg/mL)**

A	0.000
B	0.023
C	0.035
D	0.100
E	0.301
F	0.977
G	2.96
H	9.40

**TNFα Calibration Curve**



**Calibrator Levels (pg/mL)**

A	0.000
B	0.088
C	0.148
D	0.446
E	1.41
F	4.66
G	14.8
H	42.5