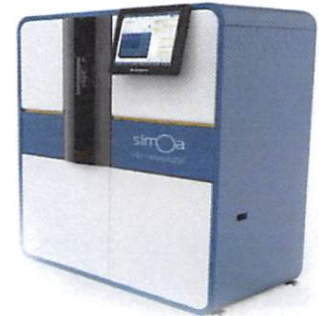


Product Number:	101160
Lot Number:	501281
Expiration:	27-Oct-2018
Platform(s):	HD-1 & SR-X



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

Component	Part Number	Lot Number
Bead Reagent †	101263	81631
Detector Reagent	101264	820604
SBG Reagent	101296	81212
Sample Diluent	101308	81175
Calibrators	101299	80880
Controls	101302 & 101316	80881
SBG Diluent (SR-X)††	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

† Follow the Simoa™ Cytokine 3-Plex A Advantage Kit Instructions (KI-0011 - HD1/KI-0101 - SR-X) to prepare the Bead Reagent.

†† Specific for use on the SR-X platform - refer to the Simoa™ Cytokine 3-Plex A SR-X Kit Instructions (KI-0101).

<sup>1</sup> Reagents are not Kit Lot Specific

**Release Materials**

**IL-6**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	80881	1.01	0.728-1.21
Control 2	80881	27.0	18.0-31.1
Panel 1	MF-072115-01	47.9	29.8-57.6
Panel 3	MF-072115-03	1.21	0.893-1.49
Panel 4	MF-072115-04	0.840	0.670-1.12

**IL-10**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	80881	0.644	0.418-0.707
Control 2	80881	12.3	9.54-14.4
Panel 1	MF-072115-01	0.957	0.818-1.28
Panel 3	MF-072115-03	1.86	1.47-2.21
Panel 4	MF-072115-04	1.78	1.53-2.29

**TNFα**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	80881	1.96	1.37-2.30
Control 2	80881	50.5	34.5-56.3
Panel 1	MF-072115-01	1.40	0.658-2.39
Panel 3	MF-072115-03	22.3	7.87-34.9
Panel 4	MF-072115-04	7.30	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

QC Supervisor

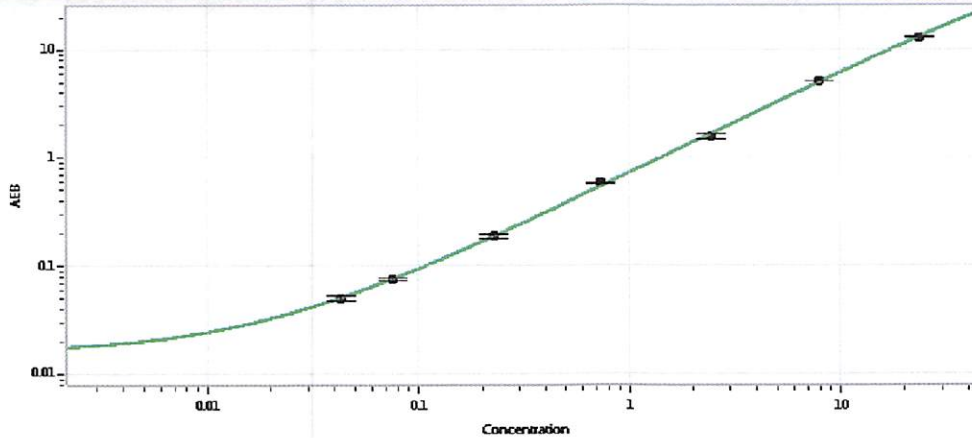
Name

Title

Signature/Date

*[Signature]* 13 AUG 2018

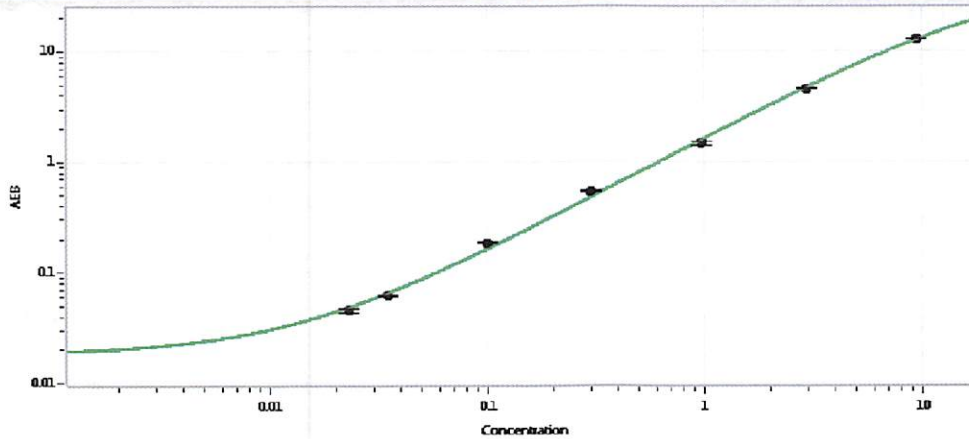
**IL-6 Calibration Curve**



**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

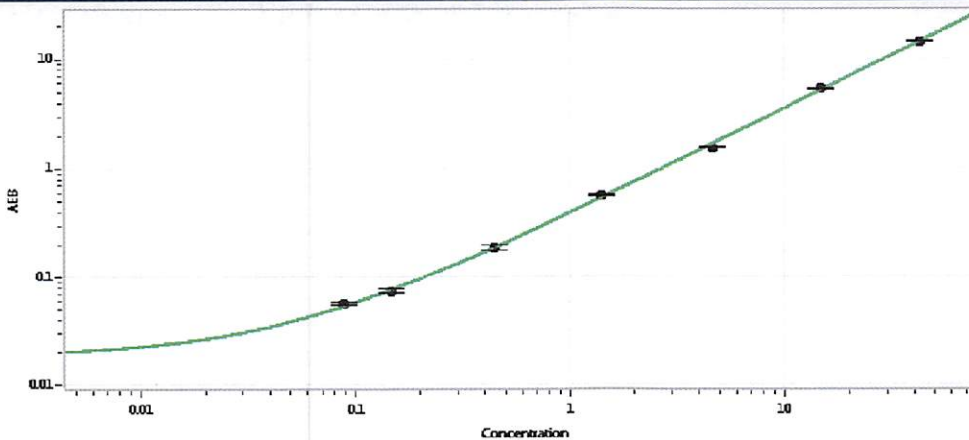
**IL-10 Calibration Curve**



**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