

<b>Product Number:</b>	101319
<b>Lot Number:</b>	502326
<b>Expiration:</b>	25-Nov-2020
<b>Platform(s):</b>	HD-1/HD-x



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

Component	Part Number	Lot Number
Bead Reagent †	101265	010203
Detector Reagent	101266	935212
SBG Reagent	101297	932912
Sample Diluent	103330	932909
Calibrators	103228	931101
Controls	103229-103230	931504
RGP Reagent	103159	N/A <sup>1</sup>

† Follow the Simoa™ Cytokine 3-Plex B Advantage Kit Instructions (KI-0012) to prepare the Bead Reagent.

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

**Release Materials**

**IL-6**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	931504	0.613	0.527-0.839
Control 2	931504	19.6	16.8-25.2
Panel 2	MF-062818-02	0.853	0.753-1.13
Panel 3	MF-062818-03	1.34	1.22-1.91
Panel 4	MF-062818-04	14.6	10.9-16.4

**IL-17A**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	931504	0.482	0.415-0.623
Control 2	931504	12.1	9.85-14.8
Panel 2	MF-062818-02	18.3	14.0-20.9
Panel 3	MF-062818-03	1.43	1.17-1.75
Panel 4	MF-062818-04	2.23	1.60-2.40

Ranges shown are generated internally for new lot release only.

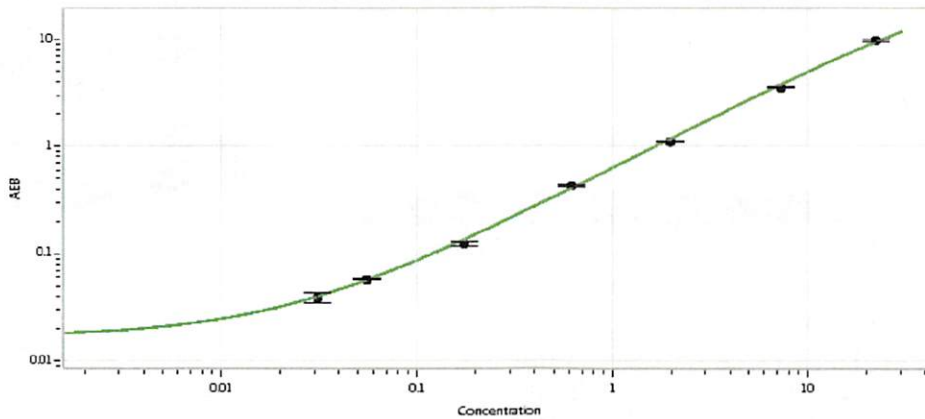
Customer should generate their own control ranges.

**TNFα**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	931504	1.77	1.59-2.38
Control 2	931504	51.3	43.1-64.7
Panel 2	MF-062818-02	1.87	1.34-2.53
Panel 3	MF-062818-03	47.2	30.6-51.8
Panel 4	MF-062818-04	7.95	5.44-8.15

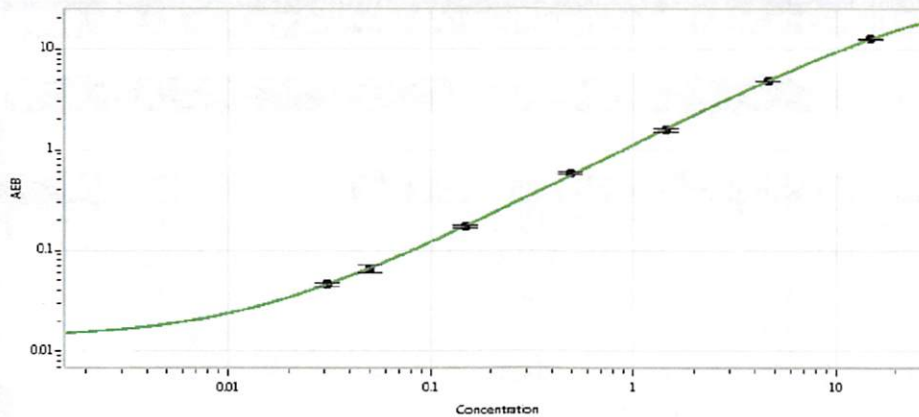
Review/Approval Shivani Goel QC Manager Shivani Goel 23 Apr 2020  
Name Title Signature/Date

**IL-6 Calibration Curve**



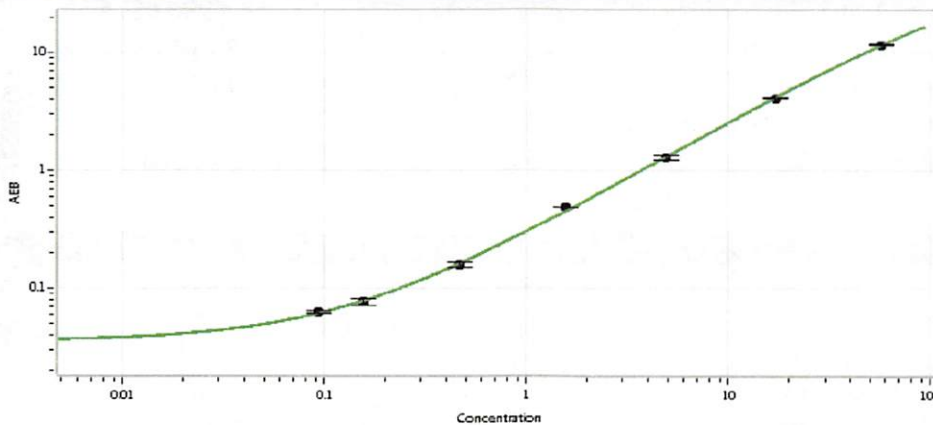
Calibrator Levels (pg/mL)	
A	0.000
B	0.031
C	0.056
D	0.176
E	0.617
F	1.99
G	7.33
H	22.5

**IL-17A Calibration Curve**



Calibrator Levels (pg/mL)	
A	0.000
B	0.031
C	0.050
D	0.150
E	0.488
F	1.45
G	4.69
H	14.8

**TNFα Calibration Curve**



Calibrator Levels (pg/mL)	
A	0.000
B	0.094
C	0.156
D	0.466
E	1.55
F	4.90
G	17.1
H	56.6