

Product Number:	101319
Lot Number:	502154
Expiration:	25-Nov-2020
Platform(s):	HD-1/HD-x



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

Component	Part Number	Lot Number
Bead Reagent †	101265	935405
Detector Reagent	101266	935212
SBG Reagent	101297	932912
Sample Diluent	103330	932909
Calibrators	103228	931101
Controls	103229-103230	931504
RGP Reagent	103159	N/A ¹

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

¹ Reagents are not Kit Lot Specific

Release Materials

IL-6

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	931504	0.726	0.560-0.841
Control 2	931504	21.0	17.0-25.6
Panel 2	MF-062818-02	0.912	0.753-1.13
Panel 3	MF-062818-03	1.28	1.22-1.91
Panel 4	MF-062818-04	12.8	10.9-16.4

IL-17A

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	931504	0.524	0.423-0.634
Control 2	931504	12.4	9.89-14.8
Panel 2	MF-062818-02	19.7	14.0-20.9
Panel 3	MF-062818-03	1.48	1.17-1.75
Panel 4	MF-062818-04	2.17	1.60-2.40

Ranges shown are generated internally for new lot release only.

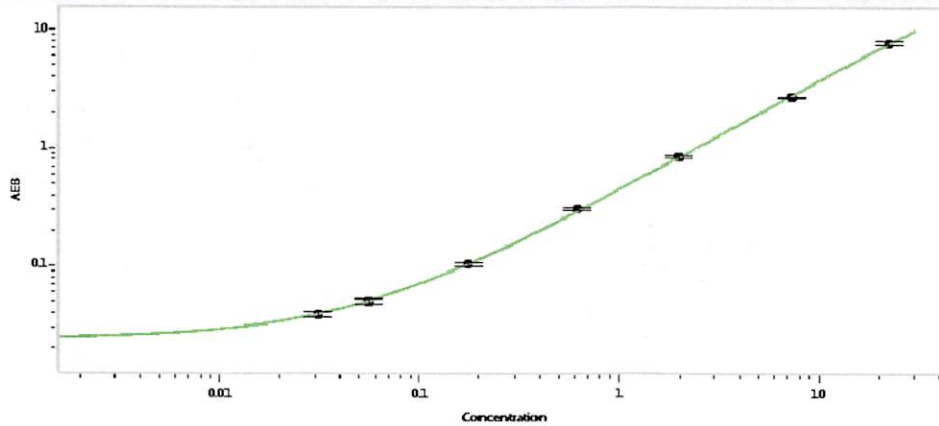
Customer should generate their own control ranges.

TNFα

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	931504	2.00	1.63-2.45
Control 2	931504	55.0	43.6-65.4
Panel 2	MF-062818-02	2.33	1.34-2.53
Panel 3	MF-062818-03	41.1	30.6-51.8
Panel 4	MF-062818-04	7.44	5.44-8.15

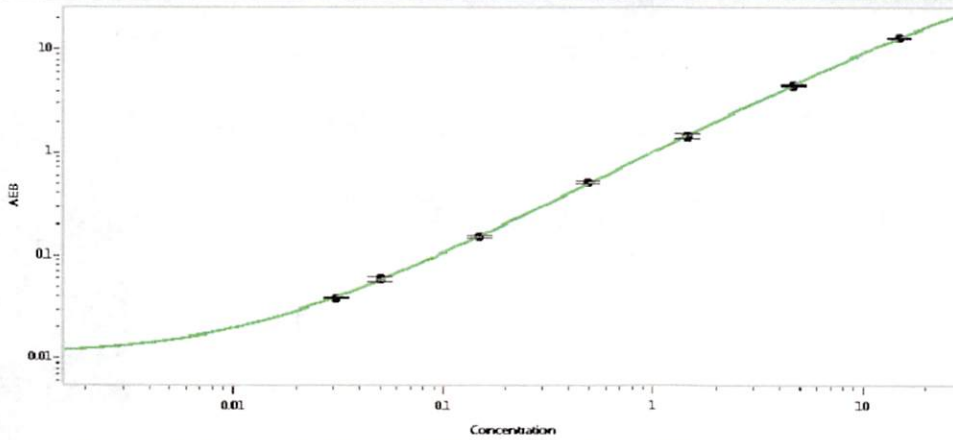
Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i>	10 Jan 2020
	Name	Title	Signature/Date	

IL-6 Calibration Curve



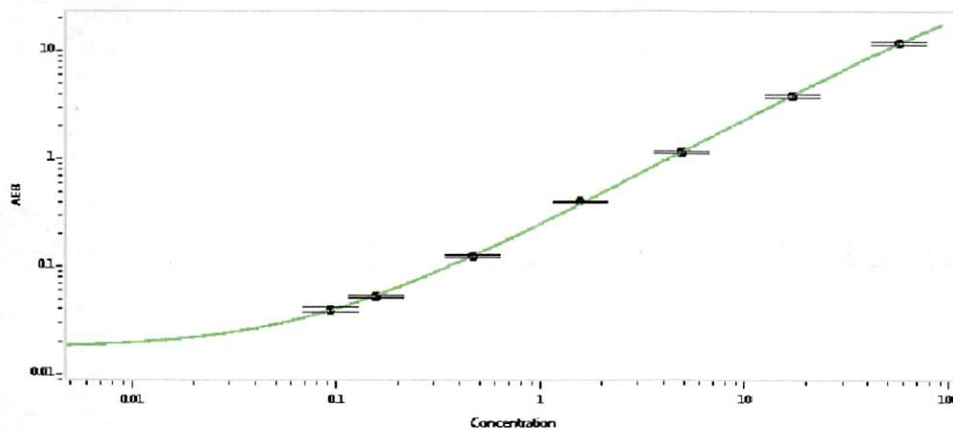
Calibrator Levels (pg/mL)	
A	0.000
B	0.031
C	0.056
D	0.176
E	0.617
F	1.99
G	7.33
H	22.5

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