

Product Number:	101319
Lot Number:	502071
Expiration:	17-Jul-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	930208
Detector Reagent	101266	930215
SBG Reagent	101297	930811
Sample Diluent	103330	919806
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	0.672	0.554-0.830
Control 2	931504	20.0	17.1-25.7
Panel 2	MF-062818-02	0.846	0.753-1.13
Panel 3	MF-062818-03	1.33	1.22-1.91
Panel 4	MF-062818-04	13.1	10.9-16.4

IL-17A

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	931504	0.549	0.424-0.636
Control 2	931504	12.1	9.88-14.8
Panel 2	MF-062818-02	18.5	14.0-20.9
Panel 3	MF-062818-03	1.40	1.17-1.75
Panel 4	MF-062818-04	2.10	1.60-2.40

TNFα

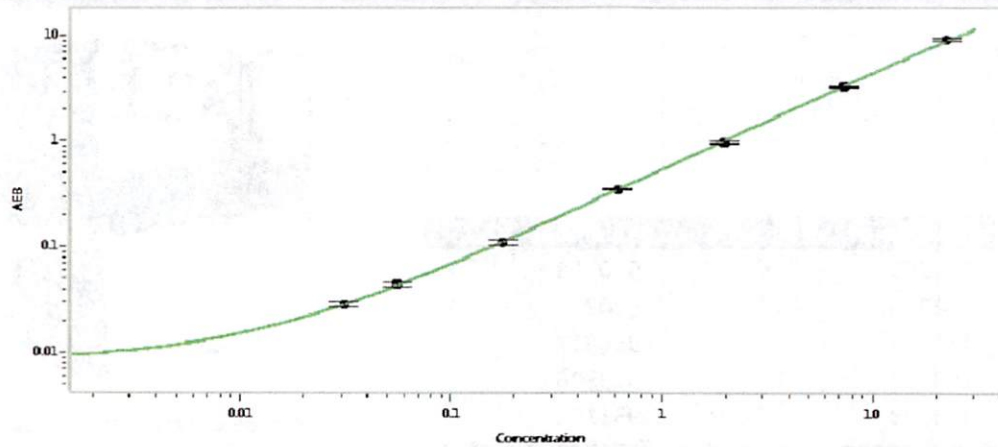
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	931504	1.96	1.63-2.44
Control 2	931504	51.6	43.9-65.8
Panel 2	MF-062818-02	1.68	1.34-2.53
Panel 3	MF-062818-03	35.4	30.6-51.8
Panel 4	MF-062818-04	6.32	5.44-8.15

Ranges shown are generated internally for new lot release only.

Customer should generate their own control ranges.

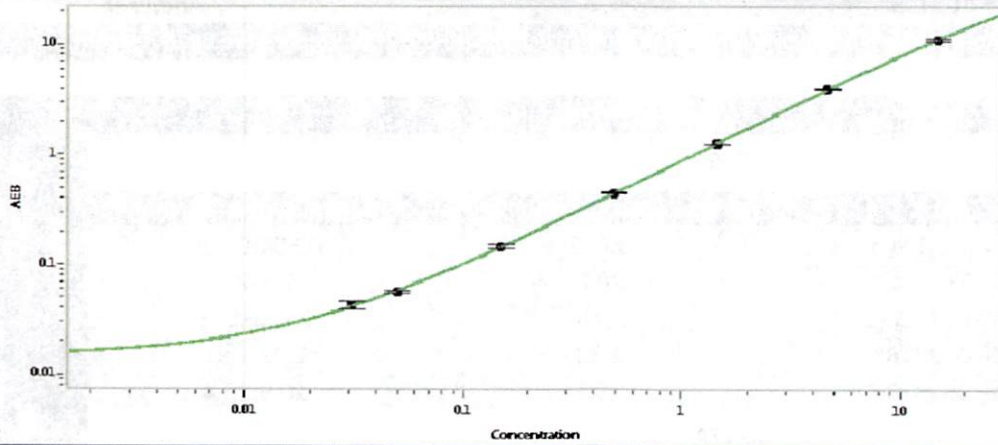
Review/Approval	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i> 26 NOV 2019
	Name	Title	Signature/Date

IL-6 Calibration Curve



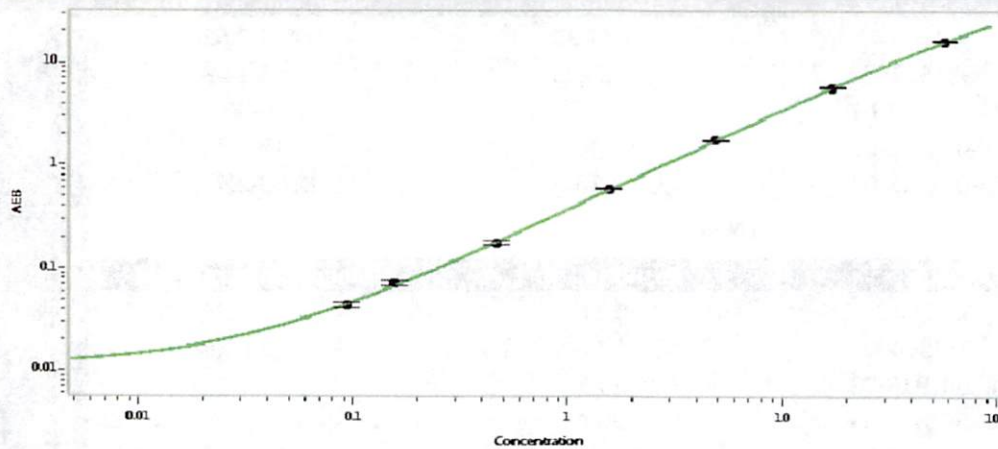
Calibrator Levels (pg/mL)	
A	0.000
B	0.031
C	0.056
D	0.176
E	0.617
F	1.99
G	7.33
H	22.5

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

TNFi 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