

Product Number:	101319
Lot Number:	501822
Expiration:	6-May-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	912904
Detector Reagent	101266	913403
SBG Reagent	101297	912603
Sample Diluent	103330	912602
Calibrators	103228	917201
Controls	103229-102330	917202
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	917202	0.744	0.596-0.894
Control 2	917202	21.0	17.0-25.5
Panel 2	MF-062818-02	0.880	0.754-1.13
Panel 3	MF-062818-03	1.25	1.20-1.92
Panel 4	MF-062818-04	12.6	10.9-16.4

IL-17A

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	917202	0.515	0.426-0.638
Control 2	917202	11.4	9.26-13.9
Panel 2	MF-062818-02	18.2	13.8-20.7
Panel 3	MF-062818-03	1.52	1.16-1.75
Panel 4	MF-062818-04	2.04	1.62-2.43

TNFα

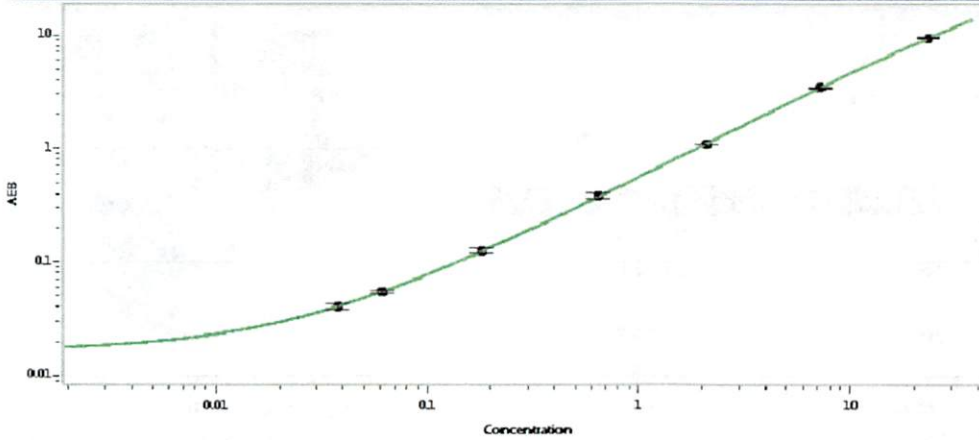
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	917202	2.63	2.05-3.08
Control 2	917202	64.0	52.5-78.7
Panel 2	MF-062818-02	2.62	1.37-2.80
Panel 3	MF-062818-03	45.0	33.3-49.9
Panel 4	MF-062818-04	7.63	5.63-8.44

Ranges shown are generated internally for new lot release only.

Customer should generate their own control ranges.

Review/Approval Patrizia Stadler Assoc. Director, QC P. Stadler 02 JUL 2019
Name Title Signature/Date

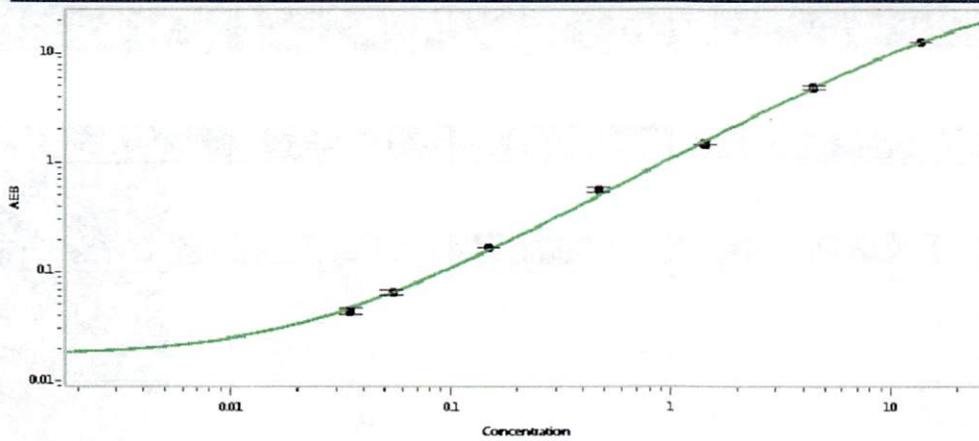
IL-6 Calibration Curve



Calibrator Levels
(pg/mL)

A	0.000
B	0.038
C	0.061
D	0.182
E	0.648
F	2.09
G	7.21
H	23.2

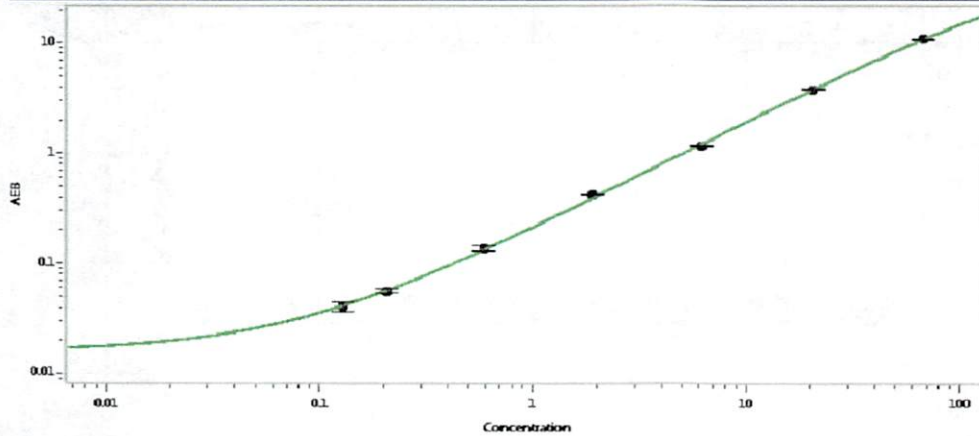
IL-17A Calibration Curve



Calibrator Levels
(pg/mL)

A	0.000
B	0.035
C	0.054
D	0.148
E	0.473
F	1.43
G	4.42
H	13.6

TNFα Calibration Curve



Calibrator Levels
(pg/mL)

A	0.000
B	0.130
C	0.207
D	0.591
E	1.92
F	6.16
G	20.6
H	67.2