

Product Number:	101319
Lot Number:	502372
Expiration:	27-Apr-2021
Platform(s):	HD-1/HD-x



Data below represents results generated on the Simoa HD-X® Analyzer.

Component	Part Number	Lot Number
Bead Reagent †	101265	012902
Detector Reagent	101266	016813
SBG Reagent	101297	018801
Sample Diluent	103330	011805
Calibrators	103228	931101
Control 1	103229	931504
Control 2	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	0.686	0.527-0.839
Control 2	931504	20.7	16.8-25.2
Panel 2	016401	1.45	1.15-1.73
Panel 3	016401	0.738	0.570-0.856
Panel 4	016401	1.07	0.842-1.32

IL-17A

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	931504	0.528	0.415-0.623
Control 2	931504	12.6	9.85-14.8
Panel 2	016401	15.5	12.5-18.8
Panel 3	016401	1.03	0.796-1.19
Panel 4	016401	1.34	1.08-1.63

TNFα

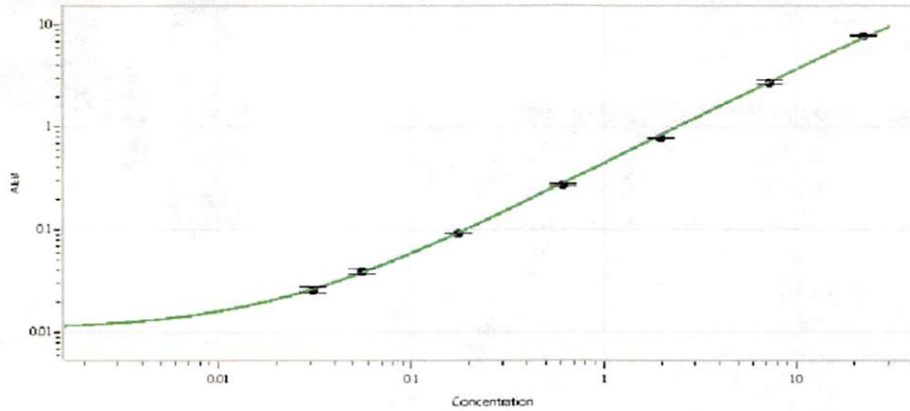
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	931504	1.97	1.59-2.38
Control 2	931504	54.4	43.1-64.7
Panel 2	016401	1.82	1.19-3.22
Panel 3	016401	25.6	16.6-38.8
Panel 4	016401	5.44	4.54-7.39

Ranges shown are generated internally for new lot release only.

Customer should generate their own control ranges.

Review/Approval	Shivani Goel	QC Manager	Shivani Goel	10/11/2020
	Name	Title	Signature/Date	

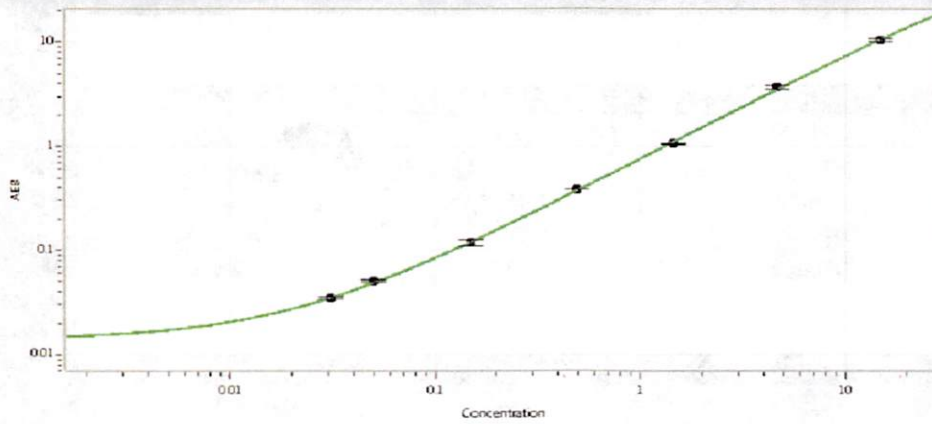
IL-6 Calibration Curve



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

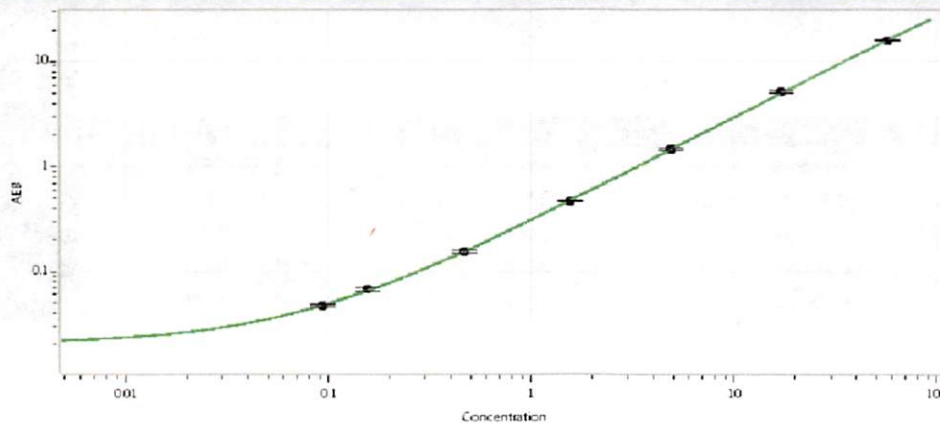
IL-17A Calibration Curve



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