

Product Number:	101319
Lot Number:	502737
Expiration:	08Feb2022
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	103901
Detector Reagent	101266	111908
SBG Reagent	101297	105403
Sample Diluent	103330	103914
Calibrators	103228	115504
Control 1	103229	115505
Control 2	103230	115505
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	115505	0.811	0.583-0.875
Control 2	115505	23.4	15.6-25.5
Panel 2	016401	1.35	1.13-1.70
Panel 3	016401	0.673	0.562-0.843
Panel 4	016401	1.05	0.858-1.29

IL-17A

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	115505	0.580	0.423-0.635
Control 2	115505	11.0	8.33-14.2
Panel 2	016401	11.7	9.72-19.7
Panel 3	016401	0.944	0.786-1.18
Panel 4	016401	1.18	1.05-1.57

TNFα

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	115505	2.14	1.54-2.31
Control 2	115505	53.0	37.9-56.9
Panel 2	016401	1.78	1.08-3.11
Panel 3	016401	19.6	12.5-38.9
Panel 4	016401	5.00	4.03-7.40

Ranges shown are generated internally for new lot release only.

Customer should generate their own control ranges.

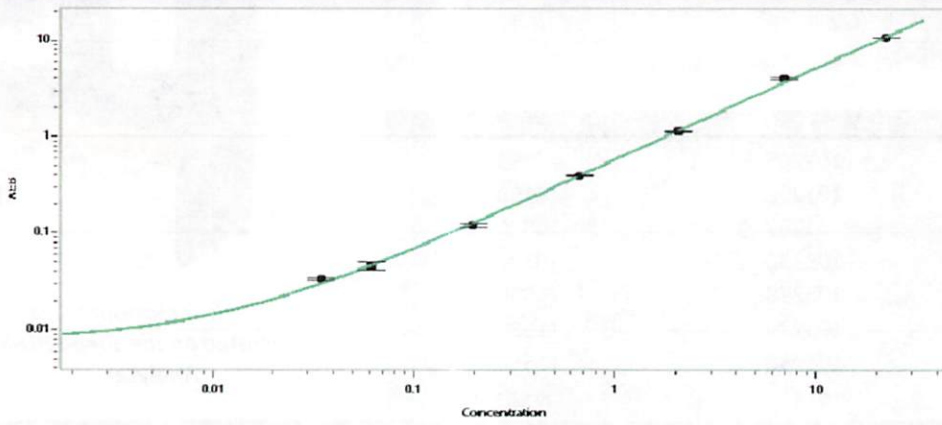
Review/Approval

Junpil Park
Name

QC Supervisor
Title

 15 Jun 2021
Signature/Date

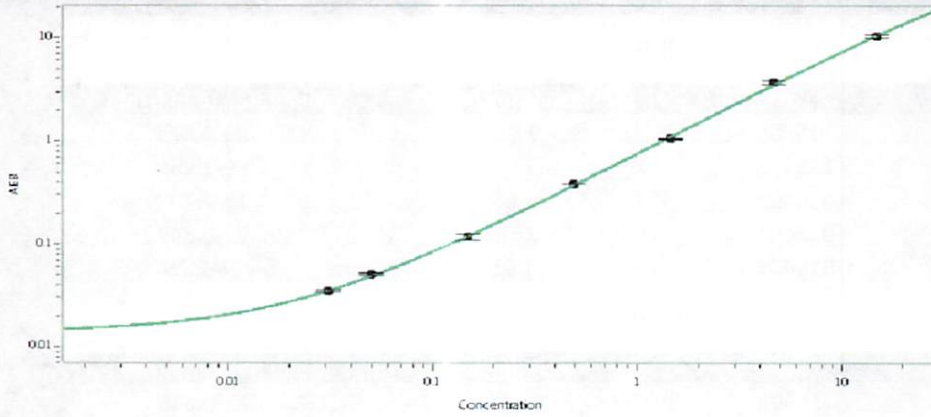
IL-6 Calibration Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	0.035
C	0.062
D	0.199
E	0.667
F	2.07
G	6.94
H	22.2

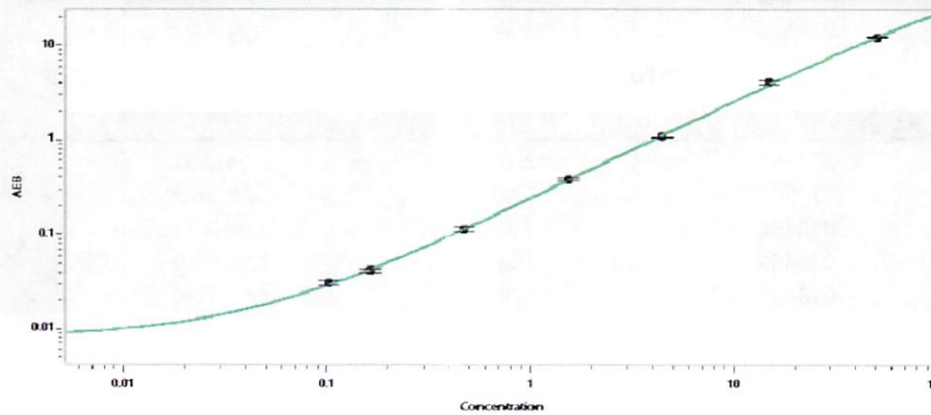
IL-17A Calibration Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	0.035
C	0.055
D	0.158
E	0.491
F	1.37
G	4.22
H	13.0

TNFα Calibration Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	0.103
C	0.164
D	0.471
E	1.54
F	4.46
G	14.9
H	50.4