

Product Number:	101160
Lot Number:	502649
Expiration:	1-Dec-2021
Platform(s):	HD-1/HD-X, SR-X

Component	Part Number	Lot Number
Bead Reagent †	101263	033607
Detector Reagent	101264	112712
SBG Reagent	101296	033903
Sample Diluent	103329	033902
Calibrators	103228	115504
Control 1	103229	115505
Control 2	103230	115505
SBG Diluent (SR-X)††	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> Reagents are not Kit Lot Specific

† Follow the Simoa™ Cytokine 3-Plex A Advantage Kit Instructions (KI-0011 - HD1/KI-0101 - SR-X) to prepare the Bead Reagent.

†† Specific for use on the SR-X platform - refer to the Simoa™ Cytokine 3-Plex A SR-X Kit Instructions (KI-0101).



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

**Release Materials**

**IL-6**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	115505	0.832	0.636 - 0.954
Control 2	115505	21.0	16.8 - 25.4
Panel 1	016401	20.1	16.1 - 28.7
Panel 2	016401	1.53	1.13 - 1.98
Panel 3	016401	0.785	0.553 - 1.05
Panel 4	016401	1.07	0.848 - 1.41

**IL-10**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	115505	0.606	0.451 - 0.676
Control 2	115505	11.8	8.75 - 15.3
Panel 1	016401	0.350	0.270 - 0.581
Panel 2	016401	7.64	5.85 - 10.8
Panel 3	016401	0.354	0.303 - 0.535
Panel 4	016401	0.526	0.414 - 0.814

**TNFα**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	115505	2.16	1.61 - 2.41
Control 2	115505	51.4	40.3 - 62.3
Panel 1	016401	1.34	1.04 - 1.83
Panel 2	016401	2.98	1.94 - 3.66
Panel 3	016401	24.6	19.3 - 34.5
Panel 4	016401	5.80	4.64 - 9.23

Ranges shown are generated internally for new lot release only.

Customer should generate their own control ranges.

Review/Approval

Junpil Park

QC Supervisor

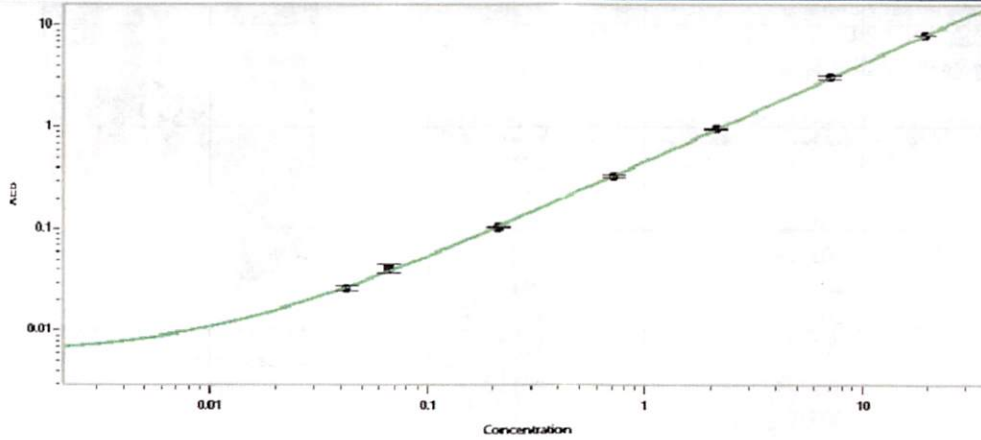
 14 Jun 2021

Name

Title

Signature/Date

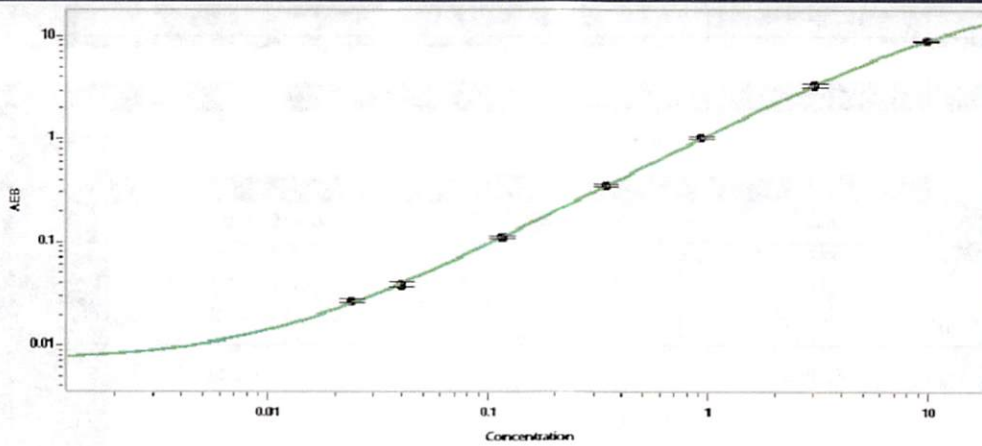
IL-6 Calibration Curve



Calibrator Levels  
(pg/mL)

A	0.000
B	0.042
C	0.066
D	0.212
E	0.719
F	2.11
G	7.03
H	19.4

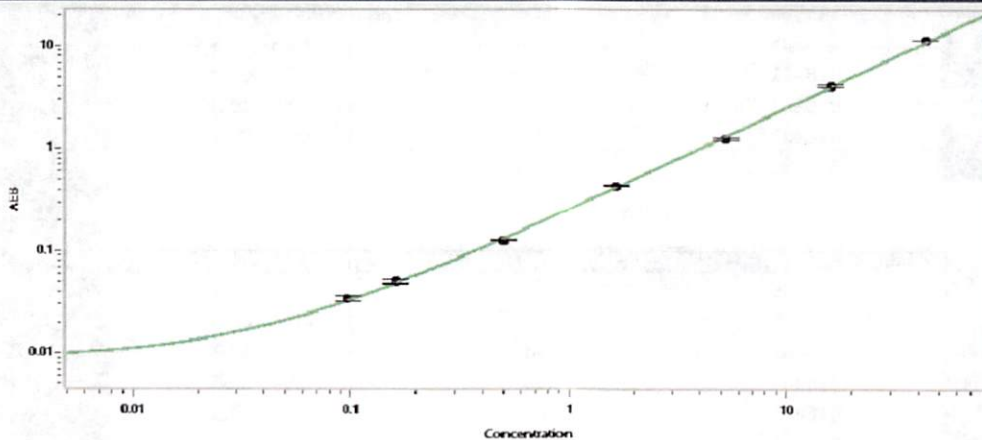
IL-10 Calibration Curve



Calibrator Levels  
(pg/mL)

A	0.000
B	0.024
C	0.040
D	0.116
E	0.341
F	0.922
G	3.02
H	9.79

TNFi Calibration Curve



Calibrator Levels  
(pg/mL)

A	0.000
B	0.097
C	0.162
D	0.504
E	1.64
F	5.24
G	15.9
H	43.5