

Product Number:	101160
Lot Number:	502571
Expiration:	13-May-2021
Platform (s):	HD-1/HD-X, SR-X

Component	Part Number	Lot Number
Bead Reagent †	101263	028208
Detector Reagent	101264	015604
SBG Reagent	101296	013405
Sample Diluent	103329	014906
Calibrators	103228	011501
Control 1	103229	011502
Control 2	103230	011502
SBG Diluent (SR-X)††	103158	N/A ¹
RGP Reagent	103159	N/A ¹



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

¹ 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).

Release Materials

IL-6

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	011502	0.636	0.475 - 0.712
Control 2	011502	16.8	13.4 - 20.1
Panel 1	016401	20.7	16.3 - 29.0
Panel 2	016401	1.48	1.11 - 2.01
Panel 3	016401	0.785	0.541 - 1.07
Panel 4	016401	1.09	0.844 - 1.43

IL-10

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	011502	0.458	0.310 - 0.609
Control 2	011502	8.58	5.86 - 12.9
Panel 1	016401	0.356	0.287 - 0.582
Panel 2	016401	7.39	5.86 - 11.0
Panel 3	016401	0.377	0.322 - 0.533
Panel 4	016401	0.517	0.431 - 0.819

TNFα

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	011502	1.93	1.44 - 2.16
Control 2	011502	50.3	39.9 - 59.8
Panel 1	016401	1.36	1.04 - 1.86
Panel 2	016401	2.58	1.96 - 3.60
Panel 3	016401	25.6	19.5 - 34.9
Panel 4	016401	6.22	4.90 - 9.25

Ranges shown are generated internally for new lot release only.

Customer should generate their own control ranges.

Review/Approval

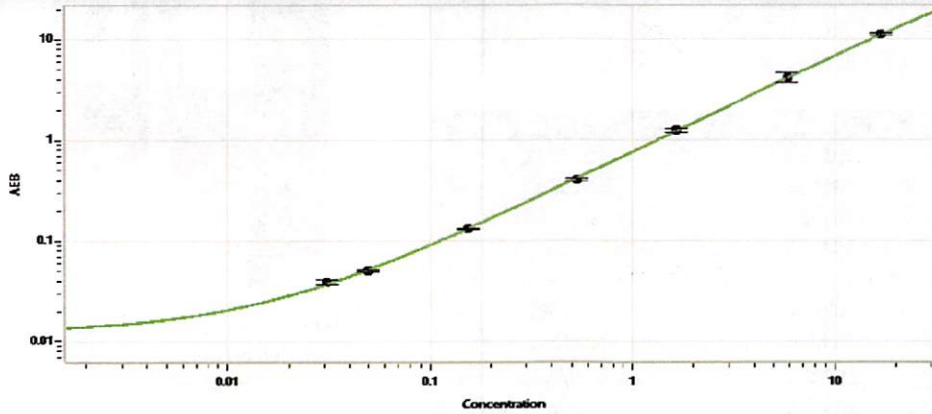
Junpil Park
Name

QC Supervisor
Title

Signature/Date

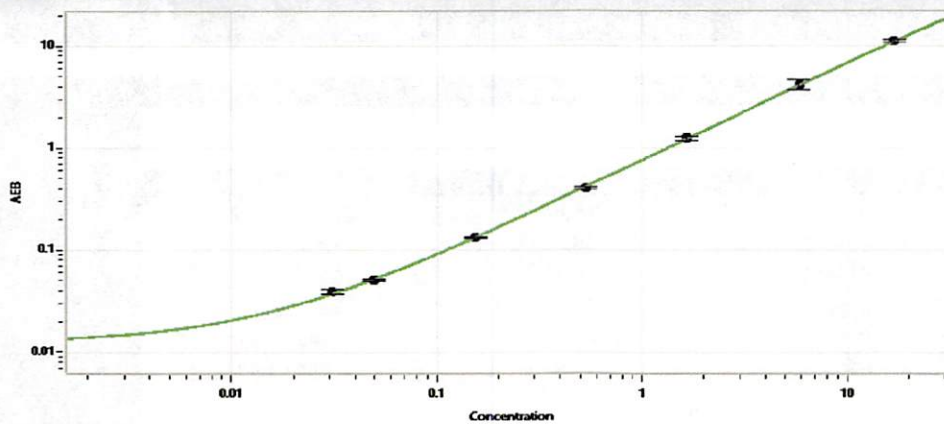
[Handwritten Signature] 03 Dec 2020

IL-6 Calibration Curve



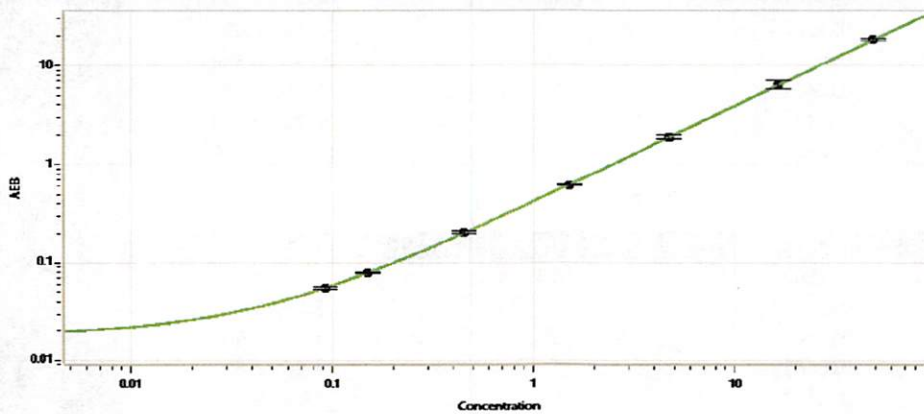
Calibrator Levels (pg/mL)	
A	0.000
B	0.031
C	0.050
D	0.154
E	0.531
F	1.65
G	5.85
H	17.0

IL-10 Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.020
C	0.031
D	0.083
E	0.249
F	0.710
G	2.35
H	7.42

TNFα Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.092
C	0.150
D	0.449
E	1.50
F	4.73
G	16.5
H	48.8