

Product Number:	101160
Lot Number:	502355
Expiration:	21-Feb-2021
Platform (s) :	HD-1/HD-X, SR-X



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

Component	Part Number	Lot Number
Bead Reagent †	101263	012101
Detector Reagent	101264	015604
SBG Reagent	101296	013405
Sample Diluent	103329	011601
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

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	0.613	0.477-0.715
Control 2	011502	16.4	13.5-20.3
Panel 2	MF-062818-02	0.908	0.777-1.17
Panel 3	MF-062818-03	1.54	1.37-2.06
Panel 4	MF-062818-04	12.3	10.4-15.7

IL-10

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	011502	0.486	0.313-0.577
Control 2	011502	10.9	7.88-12.1
Panel 2	MF-062818-02	12.7	9.23-16.5
Panel 3	MF-062818-03	0.788	0.302-1.18
Panel 4	MF-062818-04	1.32	0.908-1.69

TNFα

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	011502	1.79	1.44-2.16
Control 2	011502	49.7	40.2-60.2
Panel 2	MF-062818-02	1.86	1.54-2.41
Panel 3	MF-062818-03	43.4	31.9-47.8
Panel 4	MF-062818-04	6.97	5.51-8.26

Ranges shown are generated internally for new lot release only.
Customer should generate their own control ranges.

Review/Approval

Patrizia Stadler Assoc. Director, QC

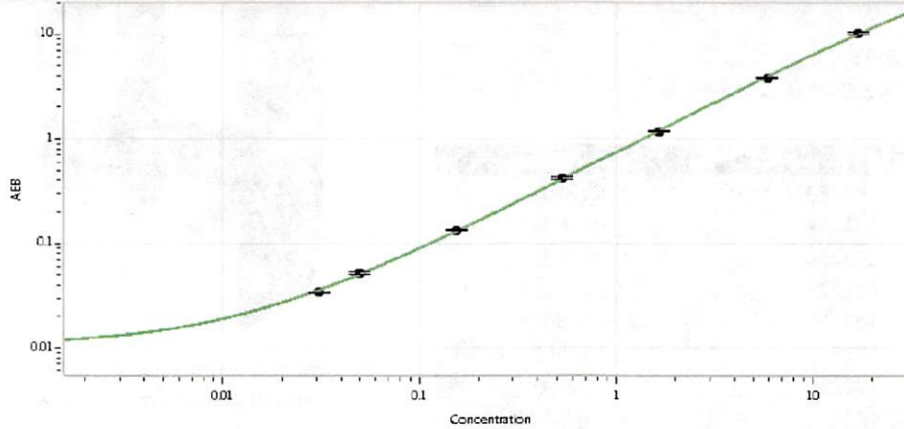
Name

Title

P. Stadler 17 Jul 2020

Signature/Date

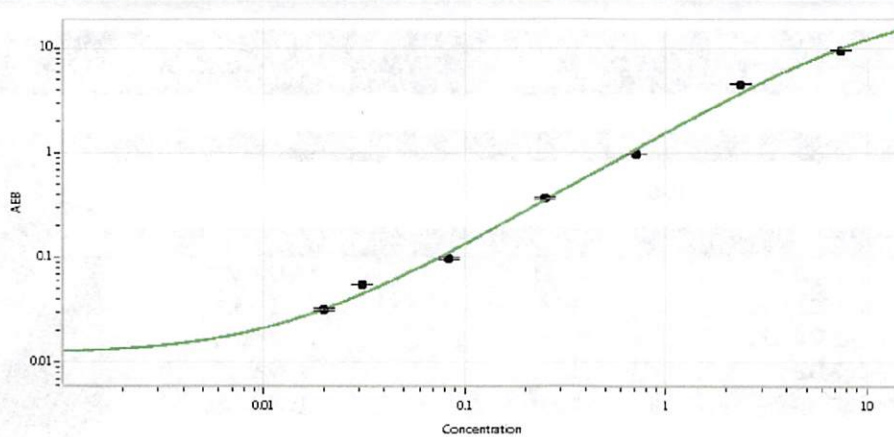
IL-6 Calibration Curve



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

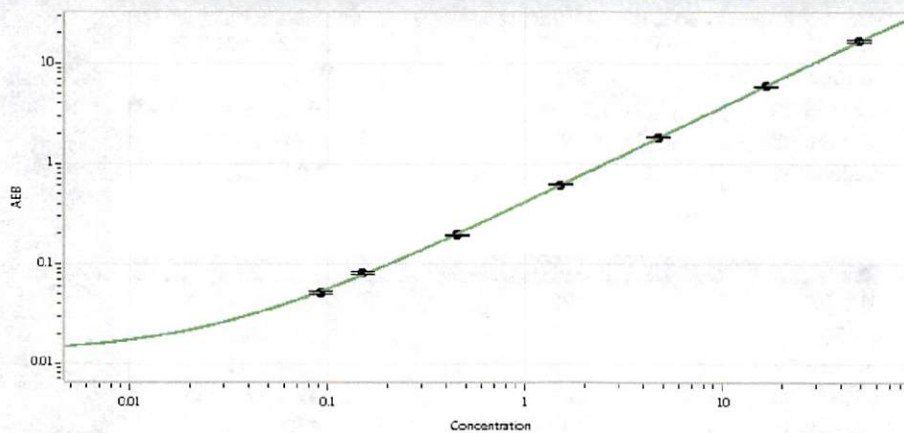
IL-10 Calibration Curve



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