

Product Number:	101160
Lot Number:	501837
Expiration:	21-Feb-2020
Platform(s):	HD-1 & SR-X



Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent †	101263	908701
Detector Reagent	101264	908702
SBG Reagent	101296	907212
Sample Diluent	103329	907211
Calibrators	101299	805202
Controls	101302/101316	805203
SBG Diluent (SR-X)††	103158	N/A ¹
RGP Reagent	103159	N/A ¹

† 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).

¹ Reagents are not Kit Lot Specific

Release Materials

IL-6

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	805203	1.10	0.873-1.31
Control 2	805203	31.1	23.2-36.5
Panel 2	MF-062818-02	0.976	0.775-1.16
Panel 3	MF-062818-03	1.61	1.38-2.08
Panel 4	MF-062818-04	15.0	10.3-15.5

IL-10

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	805203	0.667	0.382-0.749
Control 2	805203	15.1	10.4-16.7
Panel 2	MF-062818-02	13.4	9.37-15.8
Panel 3	MF-062818-03	0.631	0.393-0.769
Panel 4	MF-062818-04	1.35	0.995-1.49

TNFα

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	805203	2.06	1.57-2.35
Control 2	805203	54.4	41.3-61.9
Panel 2	MF-062818-02	2.36	1.52-2.43
Panel 3	MF-062818-03	38.6	31.9-47.8
Panel 4	MF-062818-04	7.07	5.48-8.22

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

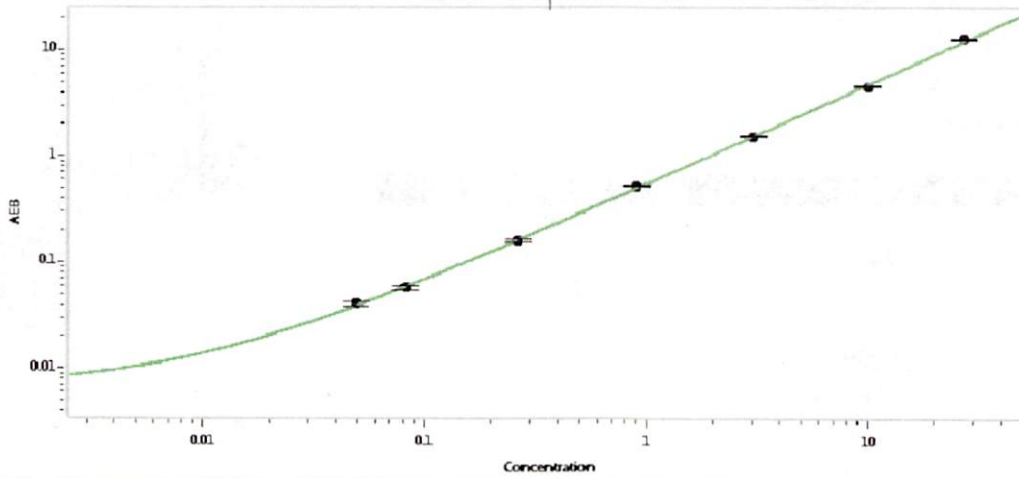
Review/Approval

Patrizia Stadler
Name

Assoc. Director, QC
Title

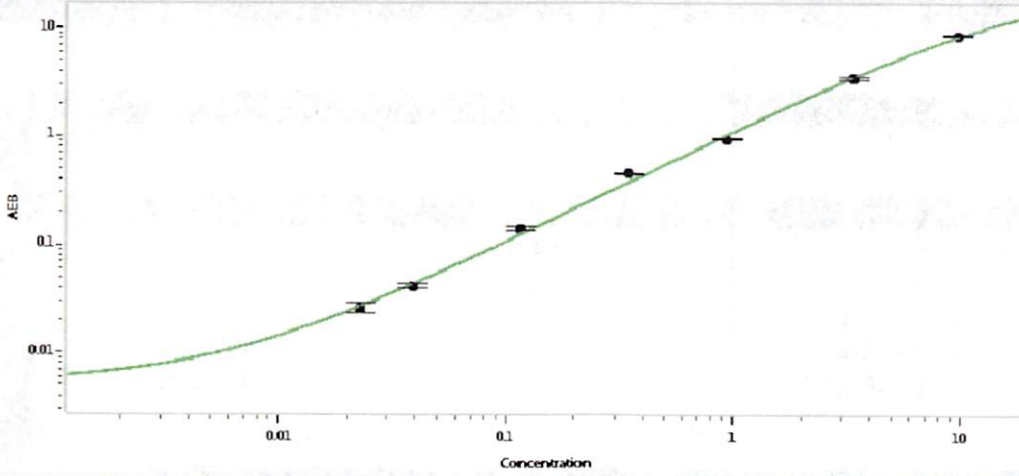
P. Stadler *20 May 2019*
Signature/Date

IL-6 Calibration Curve



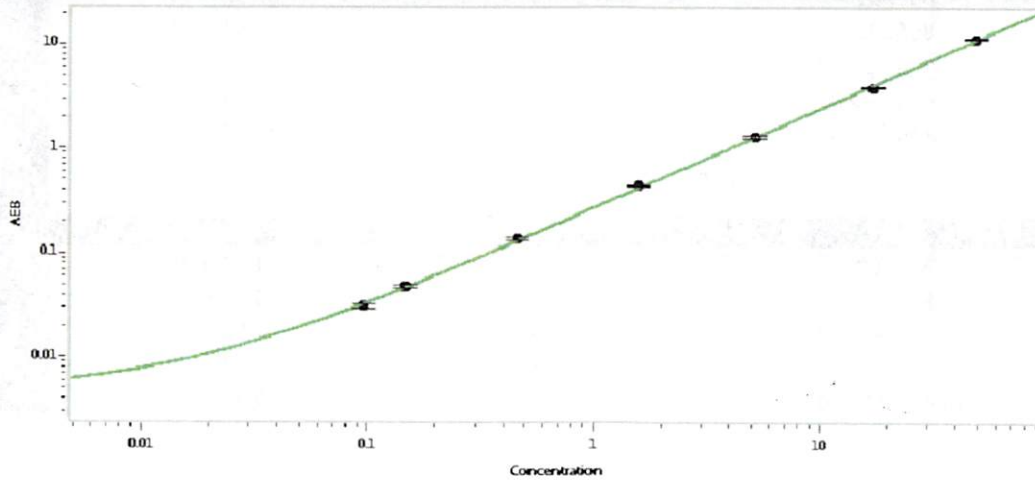
Calibrator Levels (pg/mL)	
A	0.000
B	0.049
C	0.083
D	0.263
E	0.899
F	3.01
G	10.0
H	27.1

IL-10 Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.023
C	0.039
D	0.116
E	0.345
F	0.941
G	3.36
H	9.75

TNFα Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.096
C	0.148
D	0.463
E	1.57
F	5.15
G	17.0
H	48.5