

Product Number:	101160
Lot Number:	501650
Expiration:	25-Oct-2019
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	830413
Detector Reagent	101264	830414
SBG Reagent	101296	829811
Sample Diluent	103329	829809
Calibrators	101299	81720   826802
Controls	101302 & 101316	81721
SBG Diluent (SR-X)††	103158	N/A <sup>1</sup>
RGP Reagent	103159	N/A <sup>1</sup>

† 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).

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

**Release Materials**

**IL-6**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	81721	0.949	0.711-1.07
Control 2	81721	25.4	19.5-29.2
Panel 2	MF-062818-02	1.02	0.774-1.16
Panel 3	MF-062818-03	1.68	1.41-2.12
Panel 4	MF-062818-04	14.5	10.1-15.1

**IL-10**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	81721	0.657	0.534-0.801
Control 2	81721	15.1	10.8-18.2
Panel 2	MF-062818-02	13.6	9.22-15.5
Panel 3	MF-062818-03	0.613	0.406-0.780
Panel 4	MF-062818-04	1.34	0.989-1.48

**TNFα**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	81721	1.96	1.51-2.26
Control 2	81721	50.3	38.6-57.9
Panel 2	MF-062818-02	1.92	1.52-2.29
Panel 3	MF-062818-03	44.9	31.8-47.7
Panel 4	MF-062818-04	7.81	5.39-8.09

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

Review/Approval

Patrizia Stadler

Assoc. Director, QC

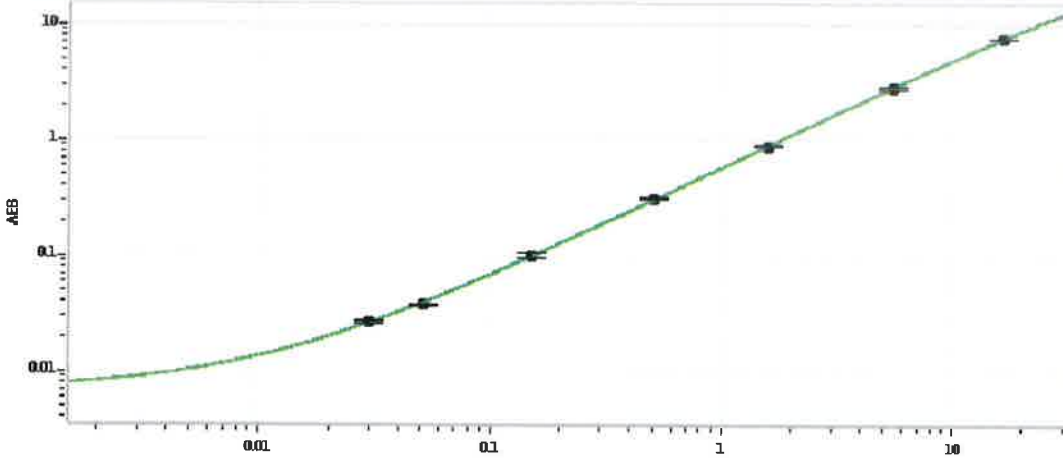
*P. Stadler 08 Feb 2019*

Name

Title

Signature/Date

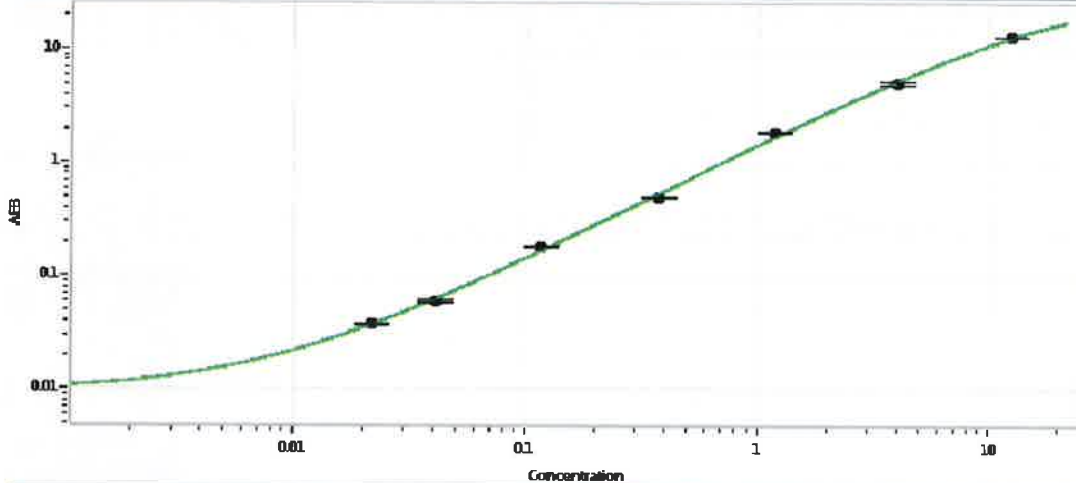
IL-6 Calibration Curve



Calibrator Levels  
(pg/mL)

A	0.000
B	0.030
C	0.051
D	0.150
E	0.511
F	1.58
G	5.50
H	16.6

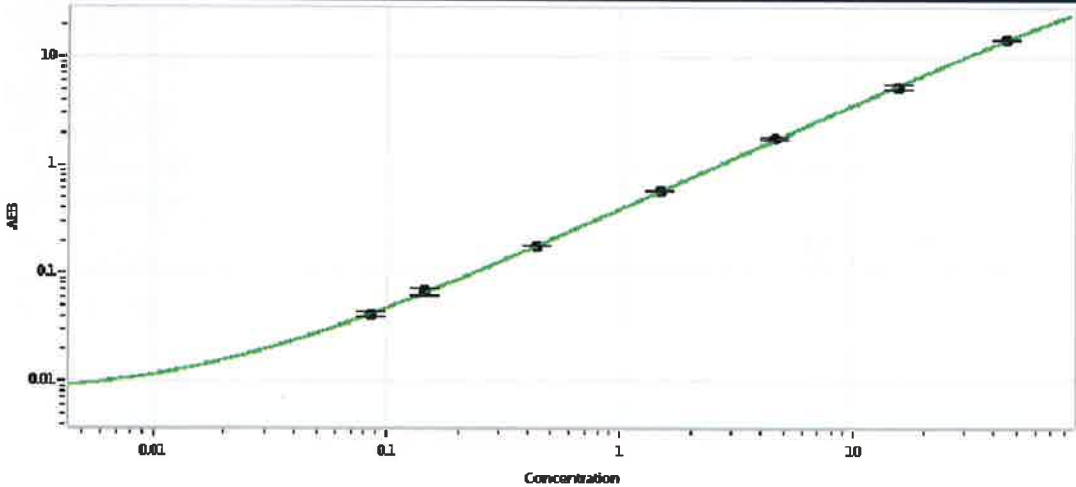
IL-10 Calibration Curve



Calibrator Levels  
(pg/mL)

A	0.000
B	0.022
C	0.041
D	0.117
E	0.375
F	1.20
G	4.02
H	12.6

TNFα Calibration Curve



Calibrator Levels  
(pg/mL)

A	0.000
B	0.086
C	0.145
D	0.440
E	1.47
F	4.56
G	15.4
H	44.9