

Product Number:	101160
Lot Number:	501951
Expiration:	17-Jul-2020
Platform(s):	HD-1/HD-X, SR-X



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

Component	Part Number	Lot Number
Bead Reagent †	101263	923409
Detector Reagent	101264	923405
SBG Reagent	101296	920307
Sample Diluent	103329	919810
Calibrators	103228	921801
Controls	103229-103230	921802
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	921802	0.735	0.600-0.900
Control 2	921802	20.8	16.8-25.2
Panel 2	MF-062818-02	0.992	0.777-1.17
Panel 3	MF-062818-03	1.62	1.37-2.06
Panel 4	MF-062818-04	14.4	10.4-15.7

IL-10

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	921802	0.466	0.379-0.568
Control 2	921802	10.6	6.85-12.5
Panel 2	MF-062818-02	15.1	10.0-15.8
Panel 3	MF-062818-03	1.08	0.302-1.18
Panel 4	MF-062818-04	1.65	0.897-1.70

TNFα

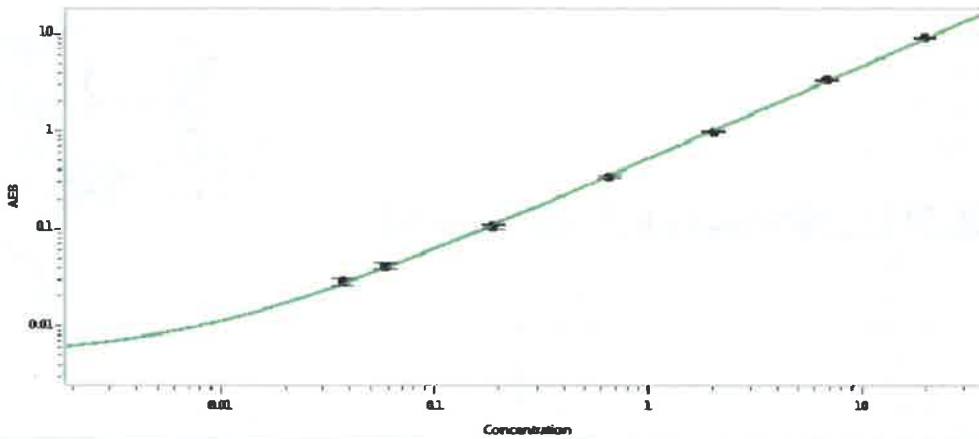
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	921802	2.59	2.03-3.04
Control 2	921802	63.5	51.0-76.5
Panel 2	MF-062818-02	1.98	1.54-2.41
Panel 3	MF-062818-03	39.8	31.9-47.8
Panel 4	MF-062818-04	7.17	5.51-8.26

Ranges shown are generated internally for new lot release only.

Customer should generate their own control ranges.

Review/Approval Patrizia Stadler Assoc. Director, QC P. Stadler 05 Sept 2019
Name Title Signature/Date

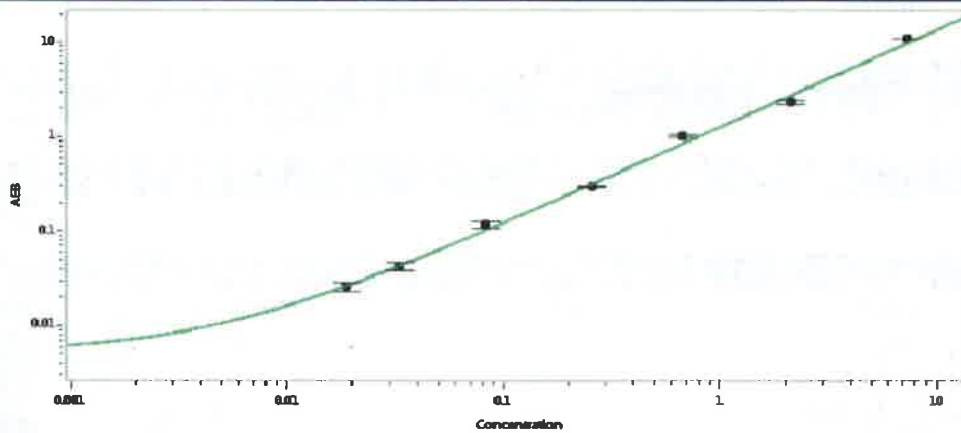
IL-6 Calibration Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	0.037
C	0.059
D	0.186
E	0.653
F	2.01
G	6.80
H	19.6

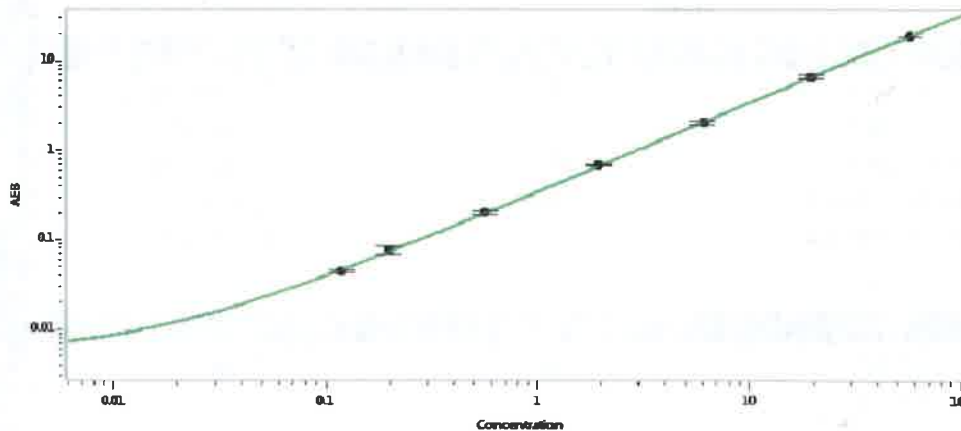
IL-10 Calibration Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	0.019
C	0.033
D	0.083
E	0.258
F	0.677
G	2.12
H	7.33

TNFα Calibration Curve



**Calibrator Levels
(pg/mL)**

A	0.000
B	0.118
C	0.197
D	0.563
E	1.93
F	6.04
G	19.6
H	56.5