

Product Number:	101160
Lot Number:	502040
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	926204
Detector Reagent	101264	926307
SBG Reagent	101296	924102
Sample Diluent	103329	919812
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	0.725	0.597-0.896
Control 2	921802	19.2	16.6-24.9
Panel 2	MF-062818-02	0.908	0.777-1.17
Panel 3	MF-062818-03	1.52	1.37-2.06
Panel 4	MF-062818-04	12.9	10.4-15.7

IL-10

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	921802	0.408	0.371-0.557
Control 2	921802	7.99	6.41-12.5
Panel 2	MF-062818-02	11.1	9.23-16.5
Panel 3	MF-062818-03	0.797	0.302-1.18
Panel 4	MF-062818-04	1.23	0.897-1.70

TNFα

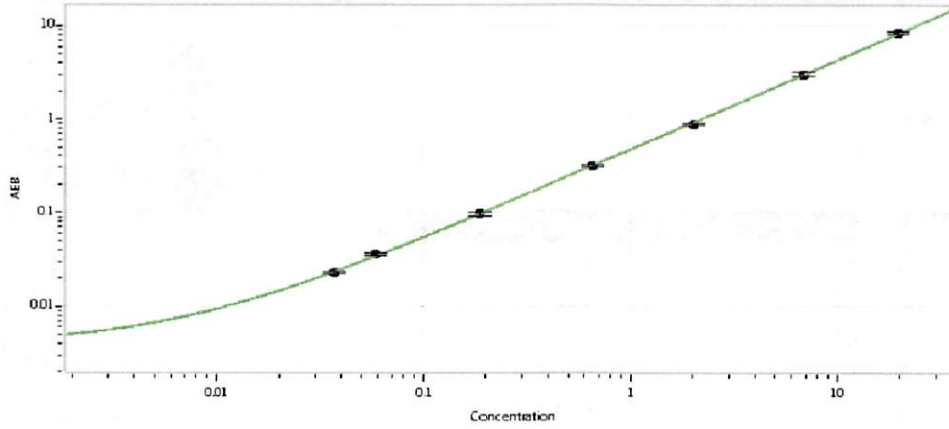
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	921802	2.38	2.01-3.02
Control 2	921802	59.3	50.5-75.7
Panel 2	MF-062818-02	1.73	1.54-2.41
Panel 3	MF-062818-03	40.7	31.9-47.8
Panel 4	MF-062818-04	6.61	5.51-8.26

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 25 Sep 2019
Name Title Signature/Date

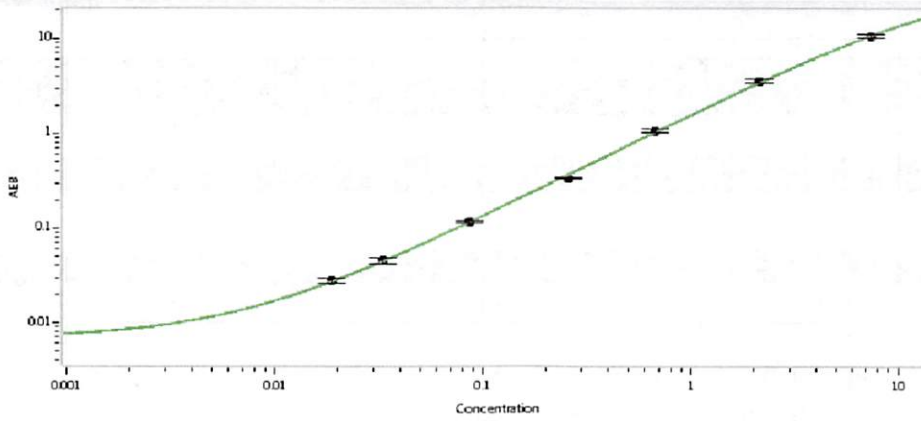
IL-6 Calibration Curve



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

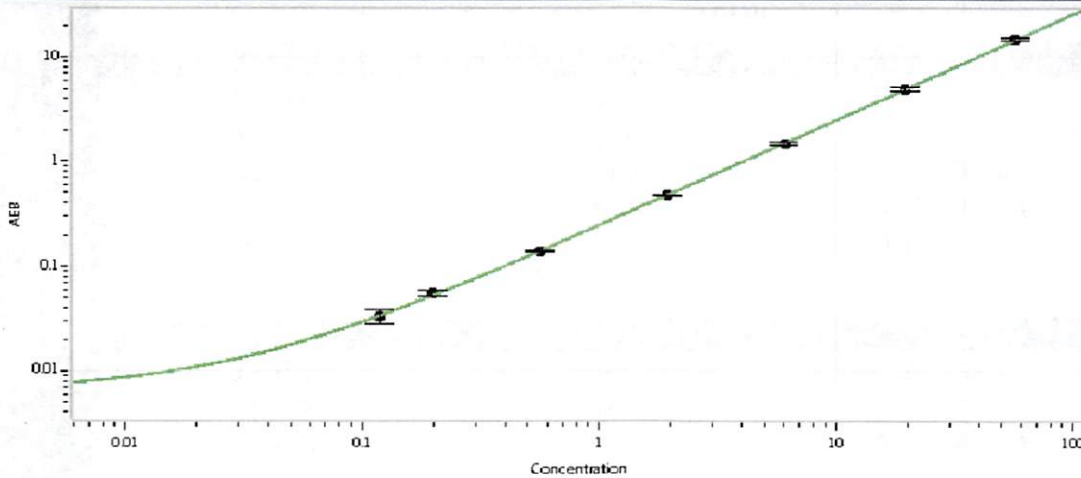
IL-10 Calibration Curve



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