

Product Number:	101160
Lot Number:	502039
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	917201
Controls	103229-103230	917202
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	921802	0.820	0.618-0.927
Control 2	921802	21.4	12.8-24.0
Panel 2	MF-062818-02	1.14	0.777-1.17
Panel 3	MF-062818-03	1.80	1.37-2.06
Panel 4	MF-062818-04	14.5	10.4-15.7

**IL-10**

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	921802	0.459	0.426-0.638
Control 2	921802	9.56	6.38-14.7
Panel 2	MF-062818-02	14.5	9.23-16.5
Panel 3	MF-062818-03	1.006	0.302-1.18
Panel 4	MF-062818-04	1.52	0.897-1.70

**TNFα**

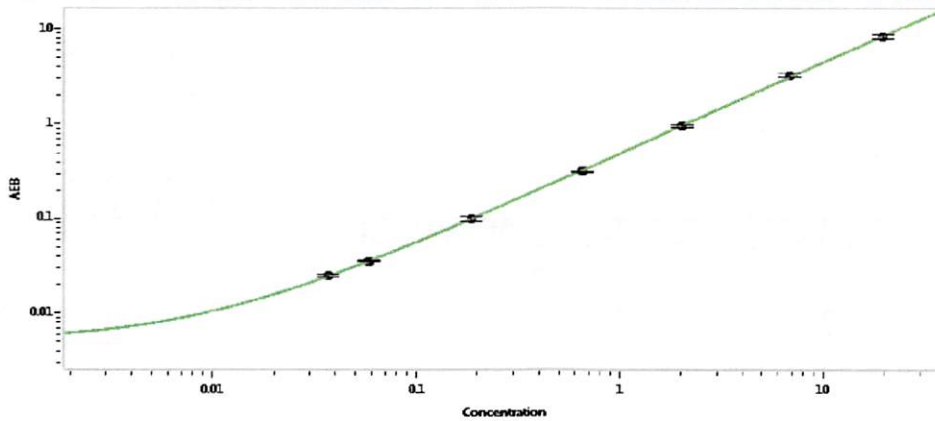
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	921802	2.78	2.06-3.09
Control 2	921802	63.8	41.3-74.9
Panel 2	MF-062818-02	1.99	1.54-2.41
Panel 3	MF-062818-03	43.4	31.9-47.8
Panel 4	MF-062818-04	8.01	5.51-8.26

Ranges shown are generated internally for new lot release only.

Customer should generate their own control ranges.

<b>Review/Approval</b>	Patrizia Stadler	Assoc. Director, QC	<i>P. Stadler</i> 26 Sep 2019
	<b>Name</b>	<b>Title</b>	<b>Signature/Date</b>

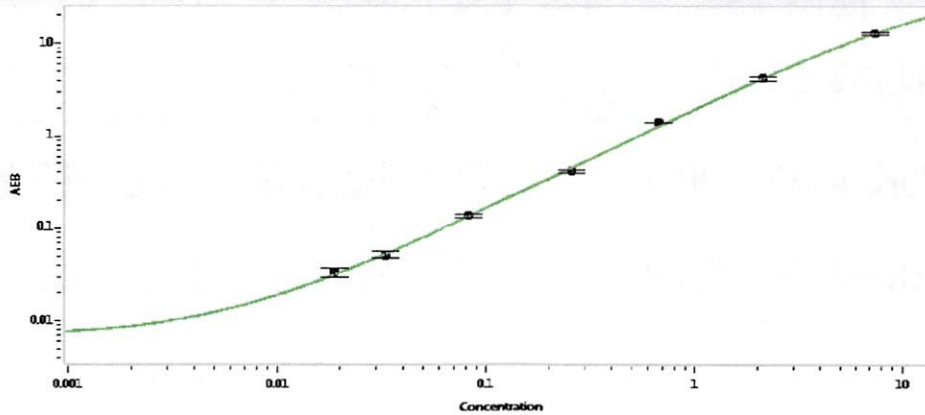
**IL-6 Calibration Curve**



**Calibrator Levels  
(pg/mL)**

A	0.000
B	0.041
C	0.071
D	0.199
E	0.658
F	2.13
G	6.82
H	19.6

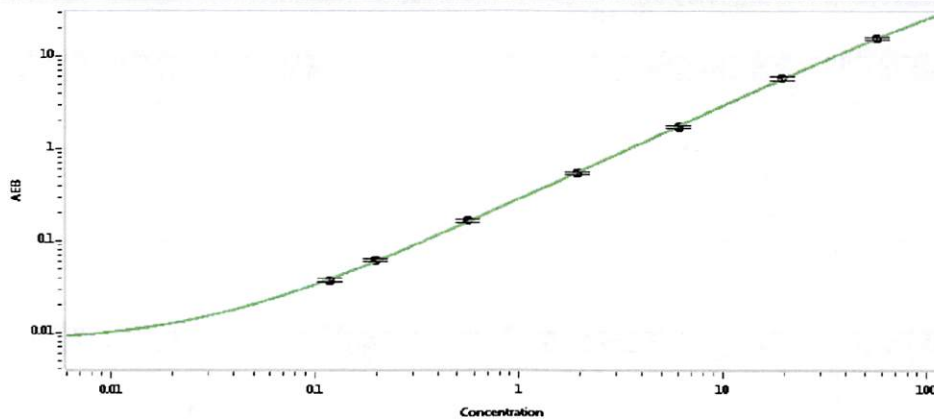
**IL-10 Calibration Curve**



**Calibrator Levels  
(pg/mL)**

A	0.000
B	0.023
C	0.036
D	0.106
E	0.292
F	0.837
G	2.71
H	9.41

**TNFα Calibration Curve**



**Calibrator Levels  
(pg/mL)**

A	0.000
B	0.130
C	0.207
D	0.606
E	2.00
F	6.42
G	20.3
H	54.7