

Product Number:	101319
Lot Number:	501894
Expiration:	21-Feb-2020
Platform(s):	HD-1/ HD-X



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

Component	Part Number	Lot Number
Bead Reagent	101265	907202
Detector Reagent	101266	907005
SBG Reagent	101297	906306
Sample Diluent	103330	906305
Calibrators	101300	80520
Controls	101303 & 101317	80521
RGP Reagent	103159	N/A ¹

¹ RGP is not Kit Lot Specific

Release Materials

IL-6

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	80521	1.00	0.791-1.19
Control 2	80521	28.4	22.3-33.4
Panel 1	MF-062818-02	0.913	0.762-1.14
Panel 2	MF-062818-03	1.48	1.29-1.93
Panel 3	MF-062818-04	12.9	11.1-16.6

IL-17A

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	80521	0.783	0.618-0.926
Control 2	80521	17.0	13.1-19.7
Panel 1	MF-062818-02	16.5	13.7-20.5
Panel 2	MF-062818-03	1.27	1.15-1.73
Panel 3	MF-062818-04	1.92	1.61-2.42

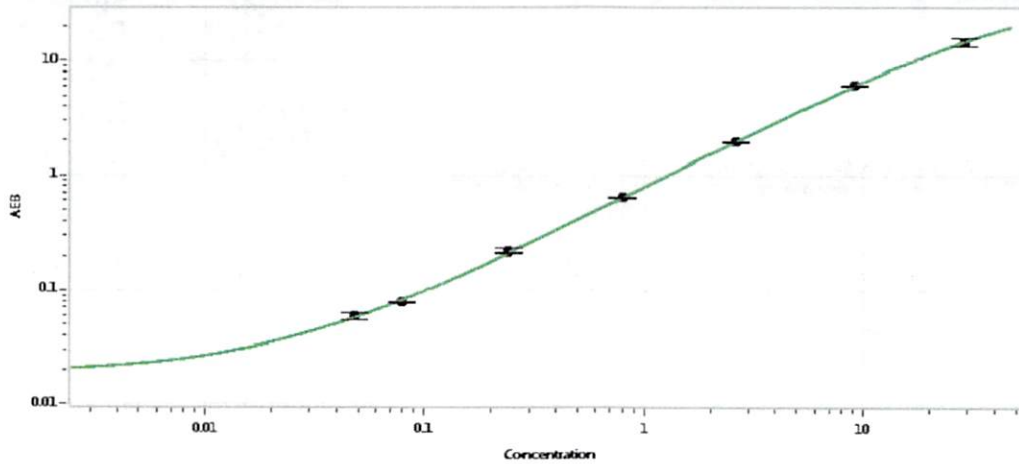
TNFα

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	80521	1.92	1.51-2.26
Control 2	80521	48.3	37.9-56.8
Panel 1	MF-062818-02	2.04	1.59-2.38
Panel 2	MF-062818-03	32.8	32.8-49.2
Panel 3	MF-062818-04	6.57	5.54-8.31

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

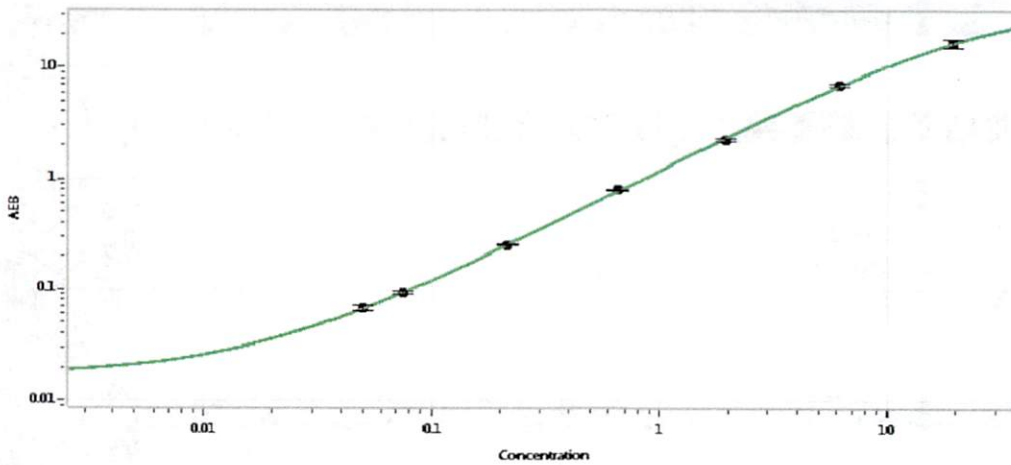
Review/Approval	Patrizia Stadler Assoc. Director, QC	<i>P. Stadler</i>	16 Jul 2019
	Name	Title	Signature/Date

IL-6 Calibration Curve



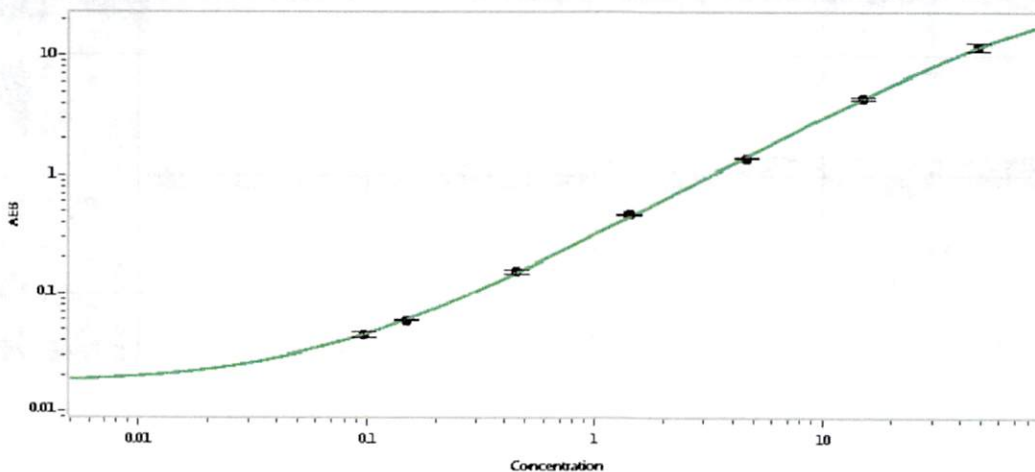
Calibrator Levels (pg/mL)	
A	0.000
B	0.048
C	0.079
D	0.241
E	0.799
F	2.63
G	9.16
H	29.0

IL-17A Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.050
C	0.075
D	0.214
E	0.648
F	1.94
G	6.13
H	19.4

TNFα Calibration Curve



Calibrator Levels (pg/mL)	
A	0.000
B	0.097
C	0.150
D	0.452
E	1.42
F	4.59
G	15.1
H	48.1