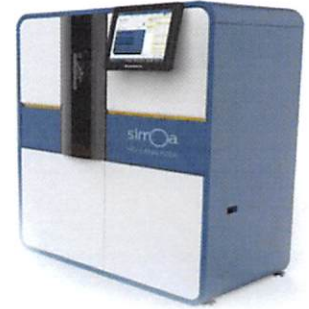


Product Number:	101319
Lot Number:	501425
Expiration:	21-Feb-2019
Platform(s):	HD-1



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

Component	Part Number	Lot Number
Bead Reagent	101265	826805
Detector Reagent	101266	826806
SBG Reagent (HD-1)†	101297	823307
Sample Diluent	103330	826206
Calibrators	101300	80520
Controls	101303 & 101317	80521
RGP Reagent	103159	N/A <sup>1</sup>

<sup>1</sup> RGP is not Kit Lot Specific

**Release Materials**

**IL-6**

	Lot	Result (pg/mL)	Historical Range (pg/mL)
Control 1	80521	1.05	0.838-1.26
Control 2	80521	29.4	22.9-34.4
Panel 2	MF-062818-02	0.927	0.782-1.17
Panel 3	MF-062818-03	1.55	1.27-1.91
Panel 4	MF-062818-04	13.1	10.8-16.2

**IL-17A**

	Lot	Result (pg/mL)	Historical Range (pg/mL)
Control 1	80521	0.803	0.651-0.976
Control 2	80521	17.4	13.6-20.4
Panel 1	MF-062818-02	17.1	14.0-21.1
Panel 3	MF-062818-03	1.52	1.18-1.78
Panel 4	MF-062818-04	2.07	1.67-2.51

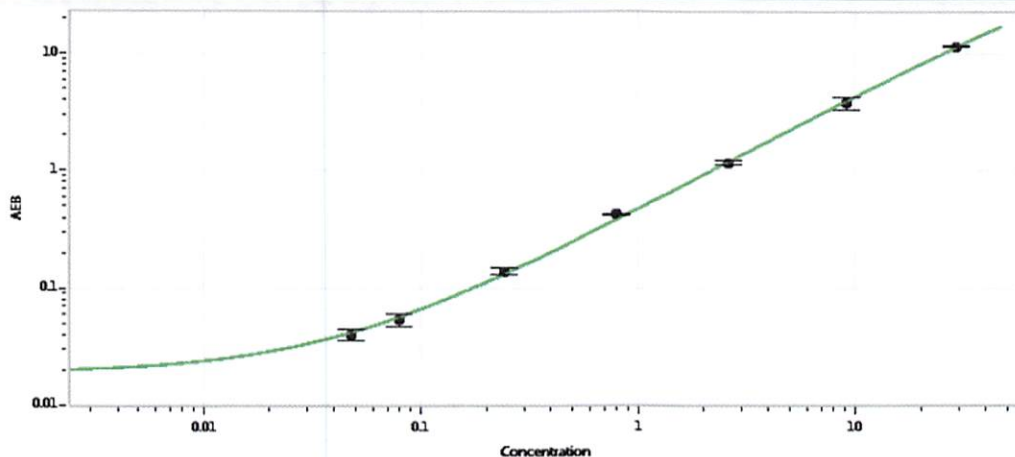
**TNFα**

	Lot	Result (pg/mL)	Historical Range (pg/mL)
Control 1	80521	2.00	1.57-2.36
Control 2	80521	49.8	39.2-58.7
Panel 1	MF-062818-02	2.02	1.51-2.27
Panel 3	MF-062818-03	40.1	33.7-50.6
Panel 4	MF-062818-04	7.11	5.63-8.44

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

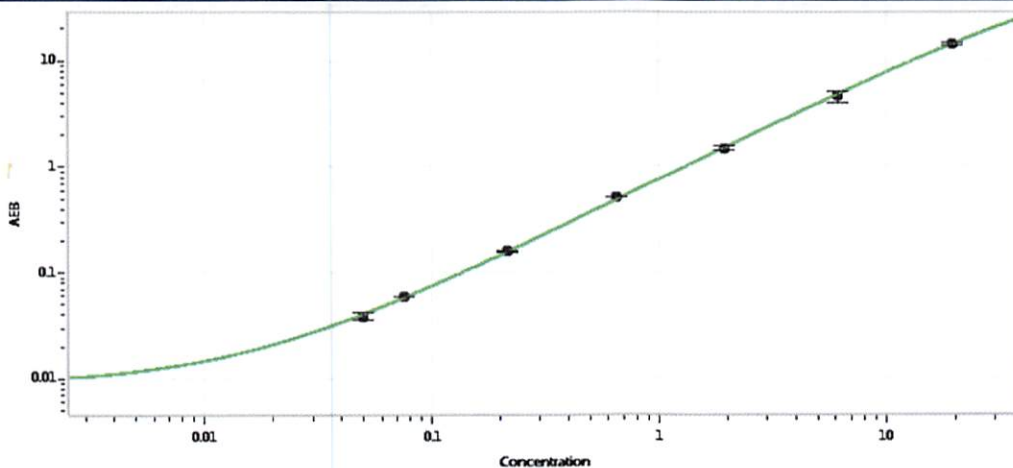
Review/Approval: Sivasankari Murugan, QC Supervisor. Signature: [Signature], Date: 03 OCT 2018.

**IL-6 Calibration Curve**



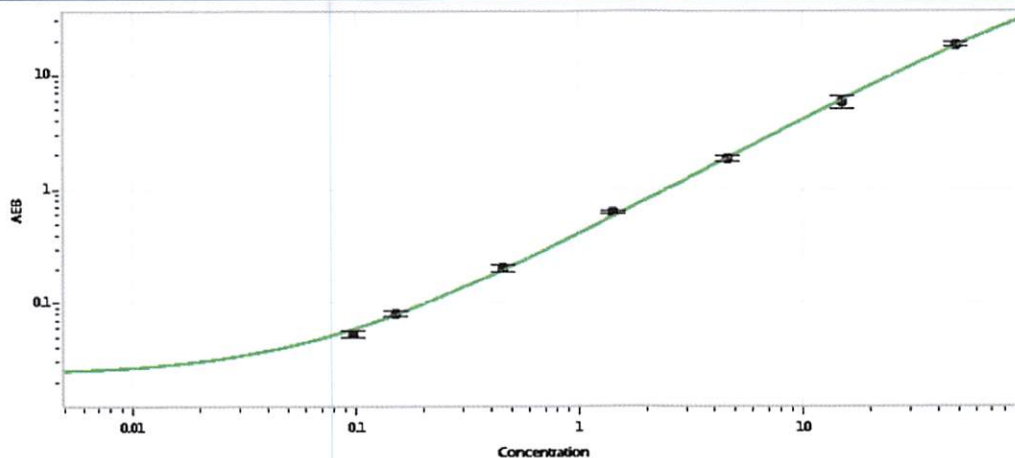
Calibrator Levels (pg/mL)	
A	0.000
B	0.048
C	0.079
D	0.241
E	0.799
F	2.63
G	9.16
H	29.0

**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