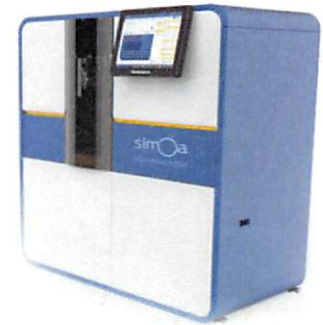


Product Number:	101319
Lot Number:	501105
Expiration:	3-Jun-2018
Platform(s):	HD-1



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

Component	Part Number	Lot Number
Bead Reagent	101265	80470
Detector Reagent	101266	80602
SBG Reagent (HD-1)†	101297	80152
Sample Diluent	101309	73373
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	1.06	0.765-1.15
Control 2	80521	27.3	21.6-32.3
Panel 1	MF-072115-01	39.5	23.4-46.9
Panel 3	MF-072115-03	1.17	0.814-1.37
Panel 4	MF-072115-04	0.719	0.562-0.993

IL-17A

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	80521	0.807	0.601-0.901
Control 2	80521	16.5	12.6-19.0
Panel 1	MF-072115-01	0.603	0.422-1.45
Panel 3	MF-072115-03	0.703	0.506-0.768
Panel 4	MF-072115-04	1.78	1.38-2.07

TNFα

	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	80521	2.01	1.46-2.20
Control 2	80521	46.9	36.9-55.4
Panel 1	MF-072115-01	0.760	0.618-2.05
Panel 3	MF-072115-03	20.7	15.9-28.2
Panel 4	MF-072115-04	6.75	5.76-8.64

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

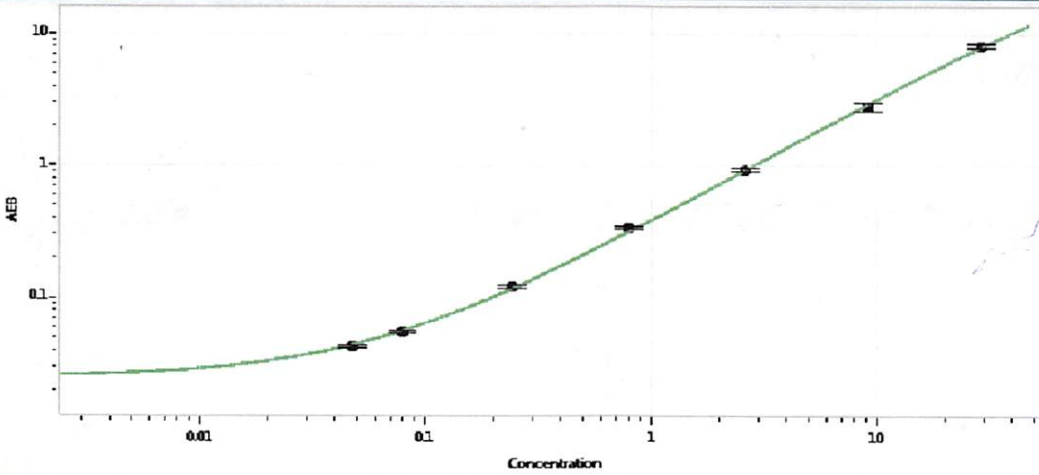
Review/Approval

Sivasankari Murugan
Name

QC Supervisor
Title

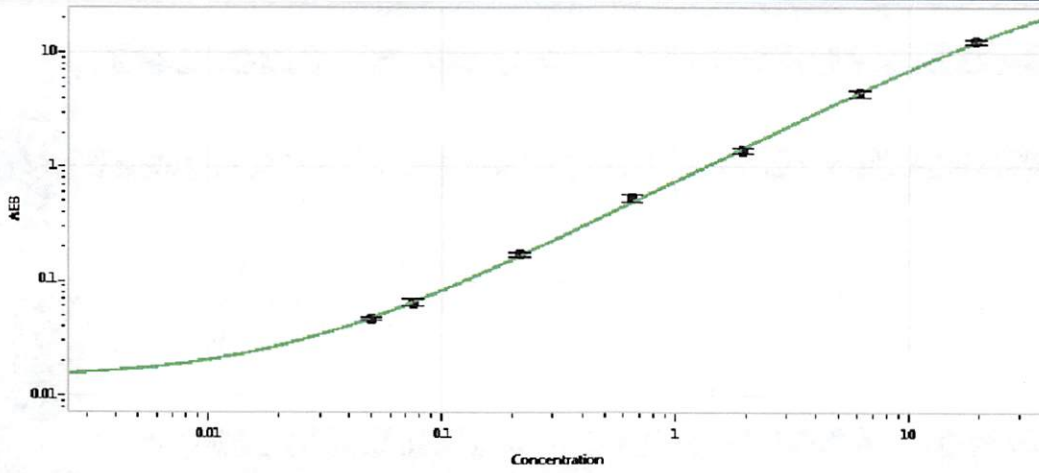
[Signature] 07 MAR 2018
Signature/Date

IL-6 Calibration Curve



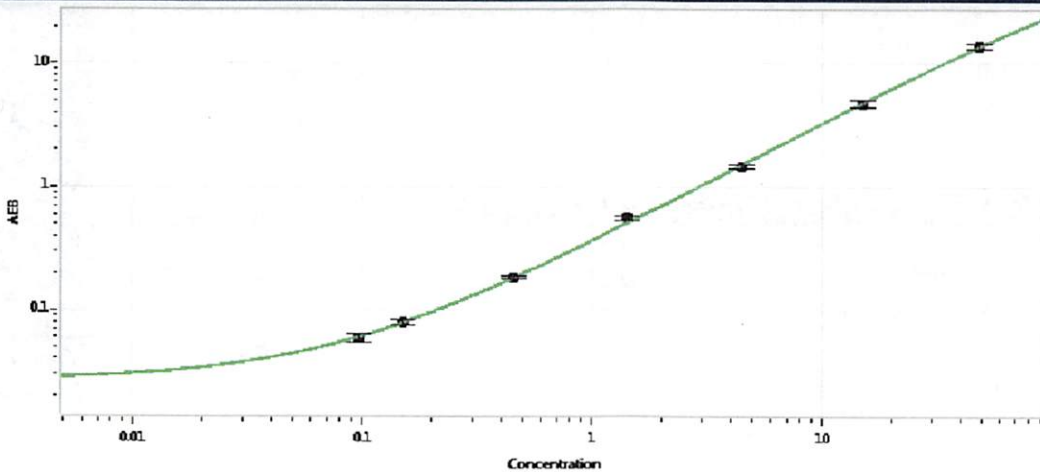
Calibrator Levels (pg/mL)	
A	0.000
B	0.048
C	0.079
D	0.241
E	0.799
F	2.63
G	9.16
H	29.0

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