

Release Materials*

IL-12p70

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	148	108 - 185
Panel 2	319839	47.1	33.5 - 51.7
Panel 3	319839	19.3	12.4 - 21.1
Panel 4	319839	4.68	2.73 - 6.45
Panel 5	319839	1.60	0.781 - 2.23

IL-1β

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	13.2	9.83 - 20.2
Panel 2	319839	4.13	2.87 - 5.85
Panel 3	319839	1.75	1.07 - 2.32
Panel 4	319839	0.458	0.250 - 0.738
Panel 5	319839	0.176	0.088 - 0.272

IL-4

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	78.8	51.1 - 98.8
Panel 2	319839	26.4	15.2 - 29.5
Panel 3	319839	10.4	6.18 - 11.4
Panel 4	319839	3.10	1.58 - 3.73
Panel 5	319839	1.16	0.531 - 1.50

IL-5

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	40.2	31.1 - 80.5
Panel 2	319839	12.9	8.64 - 23.8
Panel 3	319839	6.01	3.24 - 9.83
Panel 4	319839	1.81	0.78 - 3.01
Panel 5	319839	0.762	0.273 - 1.23

IFNγ

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	12.4	5.46 - 15.5
Panel 2	319839	3.91	1.56 - 4.89
Panel 3	319839	1.69	0.631 - 2.10
Panel 4	319839	0.497	0.200 - 0.670
Panel 5	319839	0.234	0.095 - 0.334

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

Release Materials*
IL-6

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	157	148 - 294
Panel 2	319839	46.5	41.0 - 69.5
Panel 3	319839	19.9	15.0 - 28.9
Panel 4	319839	6.97	4.50 - 11.1
Panel 5	319839	2.80	1.50 - 4.46

IL-8

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	290	189 - 393
Panel 2	319839	82.3	51.3 - 106
Panel 3	319839	32.8	20.3 - 40.9
Panel 4	319839	9.43	6.02 - 13.0
Panel 5	319839	3.90	2.18 - 5.17

IL-22

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	39.0	32.0 - 101
Panel 2	319839	7.80	5.67 - 14.9
Panel 3	319839	3.83	2.42 - 7.55
Panel 4	319839	0.905	0.54 - 1.50
Panel 5	319839	0.259	0.165 - 0.439

TNFα

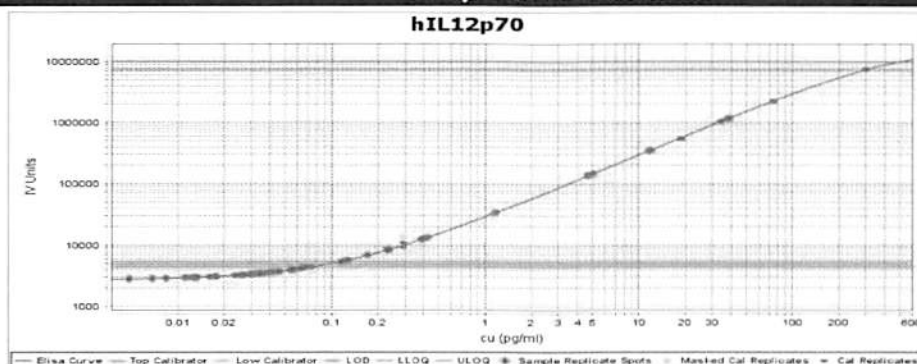
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	61.4	35.8 - 101
Panel 2	319839	19.7	11.6 - 28.3
Panel 3	319839	7.67	4.28 - 11.3
Panel 4	319839	1.90	1.16 - 3.19
Panel 5	319839	0.964	0.413 - 1.34

IL-10

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	49.9	30.3 - 61.7
Panel 2	319839	15.6	9.48 - 17.7
Panel 3	319839	6.20	3.41 - 7.65
Panel 4	319839	1.65	0.944 - 2.59
Panel 5	319839	0.688	0.437 - 1.03

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

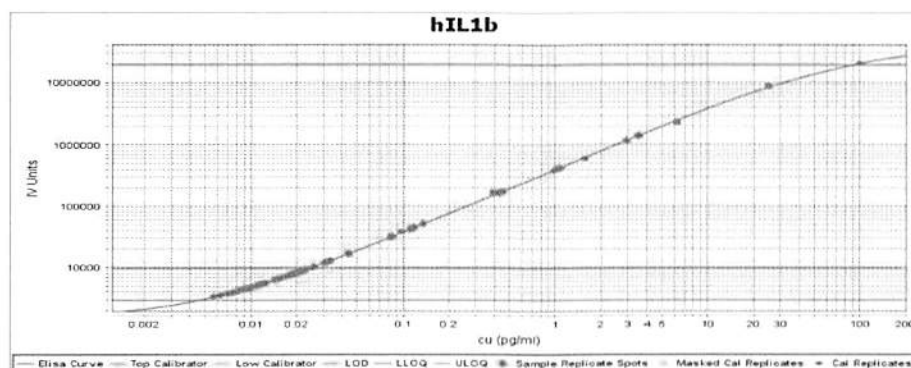
IL-12p70 Calibrator Curve



Calibrator Levels (pg/mL)

A	300
B	75.0
C	18.8
D	4.69
E	1.17
F	0.293
G	0.073
H	0.000

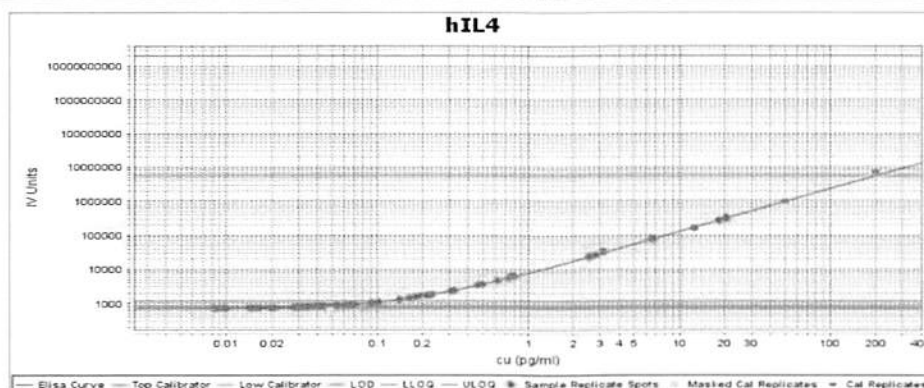
IL-1β Calibrator Curve



Calibrator Levels (pg/mL)

A	100
B	25.0
C	6.25
D	1.56
E	0.391
F	0.098
G	0.024
H	0.000

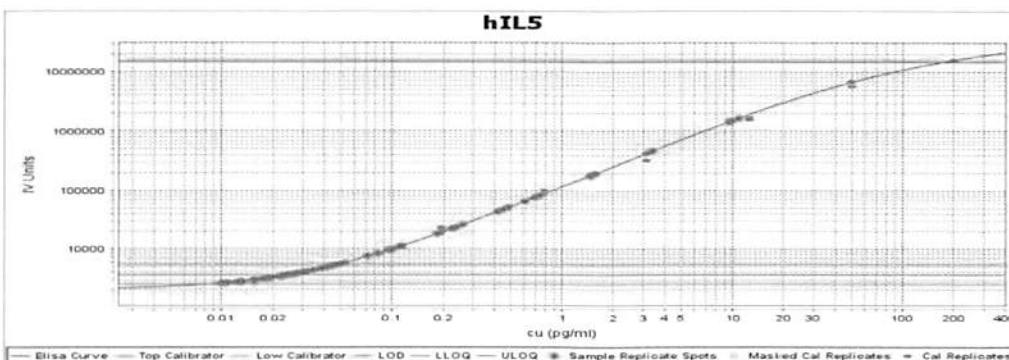
IL-4 Calibrator Curve



Calibrator Levels (pg/mL)

A	200
B	50.0
C	12.5
D	3.13
E	0.781
F	0.195
G	0.049
H	0.000

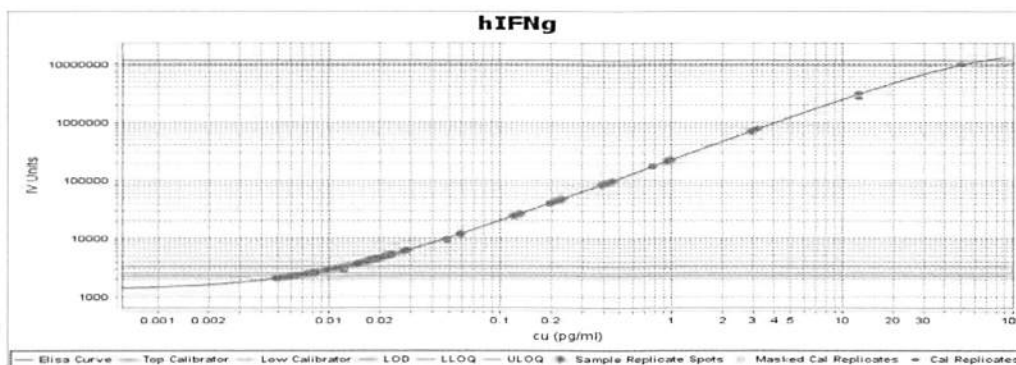
IL-5 Calibrator Curve



**Calibrator Levels
(pg/mL)**

A	200
B	50.0
C	12.5
D	3.13
E	0.781
F	0.195
G	0.049
H	0.000

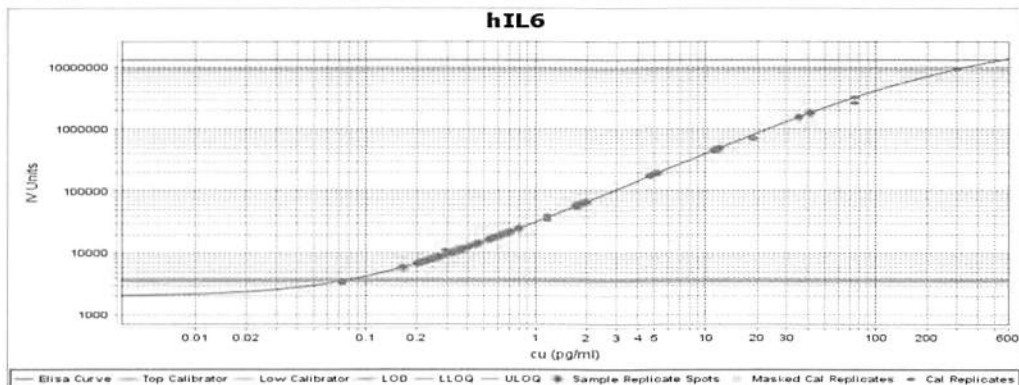
IFN γ Calibrator Curve



**Calibrator Levels
(pg/mL)**

A	50.0
B	12.5
C	3.13
D	0.781
E	0.195
F	0.049
G	0.012
H	0.000

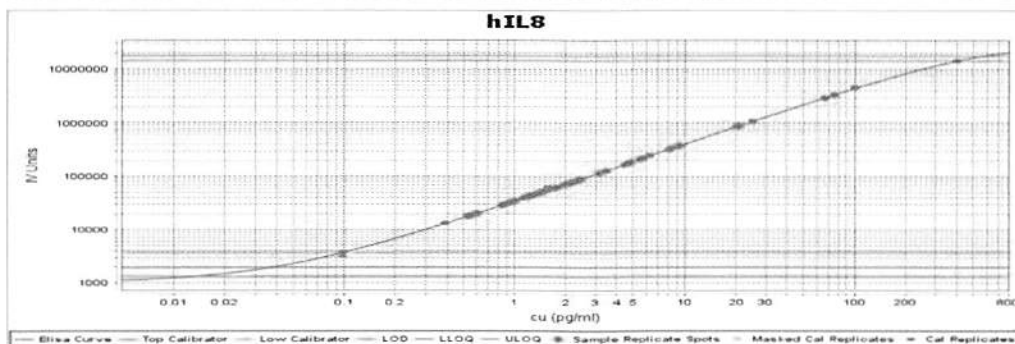
IL-6 Calibrator Curve



**Calibrator Levels
(pg/mL)**

A	300
B	75.0
C	18.8
D	4.69
E	1.17
F	0.293
G	0.073
H	0.000

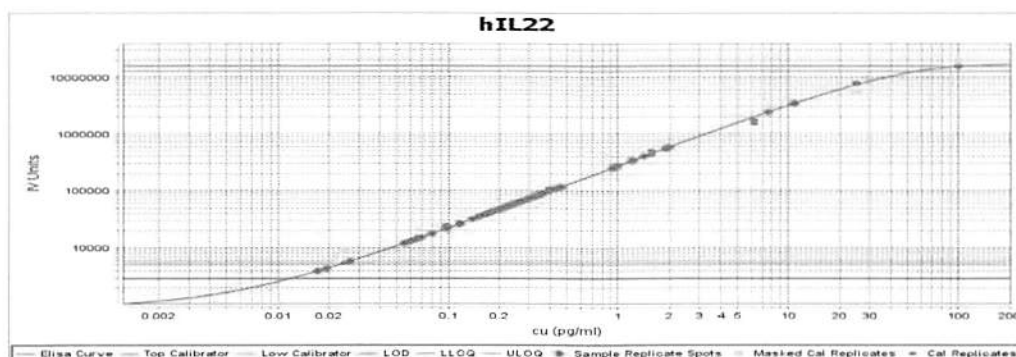
IL-8 Calibrator Curve



Calibrator Levels (pg/mL)

A	400
B	100
C	25.0
D	6.25
E	1.56
F	0.391
G	0.098
H	0.000

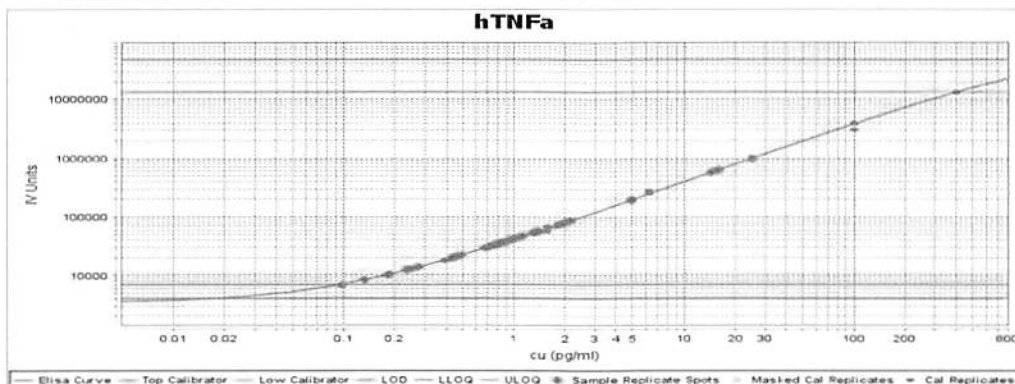
IL-22 Calibrator Curve



Calibrator Levels (pg/mL)

A	100
B	25.0
C	6.25
D	1.56
E	0.391
F	0.098
G	0.024
H	0.000

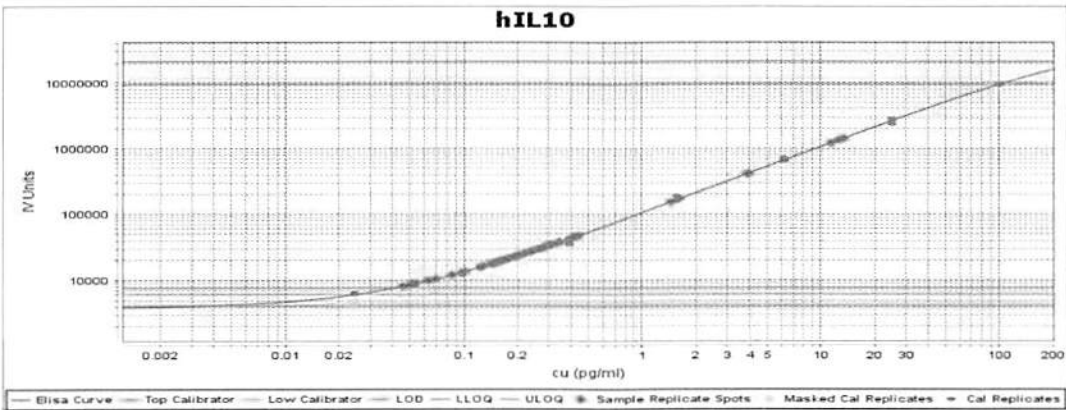
TNFα Calibrator Curve



Calibrator Levels (pg/mL)

A	400
B	100
C	25.0
D	6.25
E	1.56
F	0.391
G	0.098
H	0.000

IL-10 Calibrator Curve



Calibrator Levels (pg/mL)	
A	100
B	25.0
C	6.25
D	1.56
E	0.391
F	0.098
G	0.024
H	0.000