

Product Number:	85-0329
Lot Number:	319977
Expiration:	25-Oct-2024
Platform(s):	SP-X



Data below represents results generated on the SP-X™ Imaging and Analysis System.

CorPlex Cytokine Plate	97-0329	319977P
Biotinylated Antibody Reagent	98-0329	319977B
Sample Diluent 21	1863329	319974
IL-12p70 Calibrator	18910034	319844
IL-1β Calibrator	18910040	319564
IL-4 Calibrator	1858452	319458
IL-5 Calibrator	1858462	319838
IFNγ Calibrator	1858260	319843
IL-6 Calibrator	18910044	319484
IL-8 Calibrator	1858454	319575
IL-22 Calibrator	18910036	319459
TNFα Calibrator	1880012	319460
IL-10 Calibrator	18910043	319842
Wash Buffer (12.5X) ¹	1863332	N/A
Streptavidin-HRP Reagent ¹	1858263	N/A
Super Signal® Luminol ^{1,2}	1858089	N/A
Super Signal® Peroxide ^{1,2}	1858090	N/A

¹ Reagents are not Kit Lot Specific

² SuperSignal™ technology is protected by US Patent #6,432,662 and is a registered trademark of ThermoFisher Scientific.

All reagents are specific for use on the SP-X™ platform - refer to the Corplex Cytokine Panel Kit Instructions (KI-0195).

Review/Approval

Robert Somogyi	Quality Control Manager	
Name	Title	Signature/Date

Release Materials*

IL12p70

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	147.9	101.2 - 193.3
Panel 2	319839	40.01	32.24 - 52.86
Panel 3	319839	15.11	12.45 - 20.78
Panel 4	319839	3.937	2.632 - 6.594
Panel 5	319839	1.230	0.764 - 2.272

IL-1β

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	15.61	10.37 - 21.04
Panel 2	319839	4.267	3.172 - 5.834
Panel 3	319839	1.619	1.201 - 2.331
Panel 4	319839	0.456	0.279 - 0.773
Panel 5	319839	0.167	0.096 - 0.301

IL-4

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	79.77	45.70 - 99.67
Panel 2	319839	20.12	13.41 - 27.61
Panel 3	319839	7.567	5.612 - 11.10
Panel 4	319839	2.274	1.515 - 3.657
Panel 5	319839	0.834	0.559 - 1.400

IL-5

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	47.28	27.65 - 80.96
Panel 2	319839	14.28	8.208 - 22.84
Panel 3	319839	4.933	2.848 - 9.877
Panel 4	319839	1.569	1.085 - 2.550
Panel 5	319839	0.587	0.320 - 1.170

IFNγ

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	9.029	5.622 - 14.47
Panel 2	319839	2.569	1.692 - 4.770
Panel 3	319839	1.067	0.731 - 2.000
Panel 4	319839	0.336	0.215 - 0.692
Panel 5	319839	0.150	0.071 - 0.348

* 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	241.9	141.9 - 312.1
Panel 2	319839	50.01	39.07 - 73.71
Panel 3	319839	18.59	14.61 - 30.61
Panel 4	319839	6.105	4.493 - 11.67
Panel 5	319839	2.337	1.407 - 4.765

IL-8

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	304.8	164.0 - 399.7
Panel 2	319839	75.74	45.67 - 107.4
Panel 3	319839	26.82	18.96 - 41.09
Panel 4	319839	8.022	6.105 - 13.08
Panel 5	319839	3.119	2.260 - 5.125

IL-22

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	47.04	36.22 - 101.9
Panel 2	319839	7.640	6.825 - 14.71
Panel 3	319839	3.718	3.227 - 7.790
Panel 4	319839	0.763	0.646 - 1.509
Panel 5	319839	0.211	0.180 - 0.453

TNFα

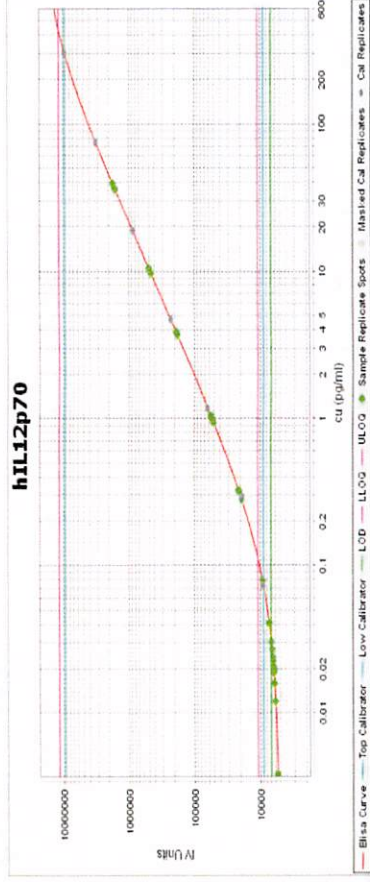
	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	53.47	37.50 - 105.1
Panel 2	319839	14.47	12.23 - 27.53
Panel 3	319839	5.732	4.430 - 11.09
Panel 4	319839	1.524	1.213 - 2.913
Panel 5	319839	0.596	0.543 - 1.056

IL-10

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Panel 1	319839	40.27	28.83 - 63.34
Panel 2	319839	11.22	9.000 - 18.28
Panel 3	319839	4.397	3.397 - 7.891
Panel 4	319839	1.385	1.141 - 2.386
Panel 5	319839	0.561	0.457 - 1.051

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

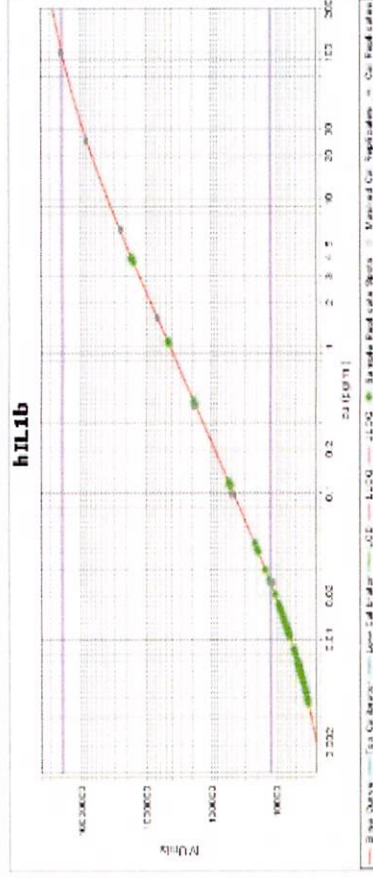
IL-12p70 Calibrator Curve



Calibrator Levels (pg/mL)

A	300.00
B	75.00
C	18.75
D	4.688
E	1.172
F	0.293
G	0.073
H	0.000

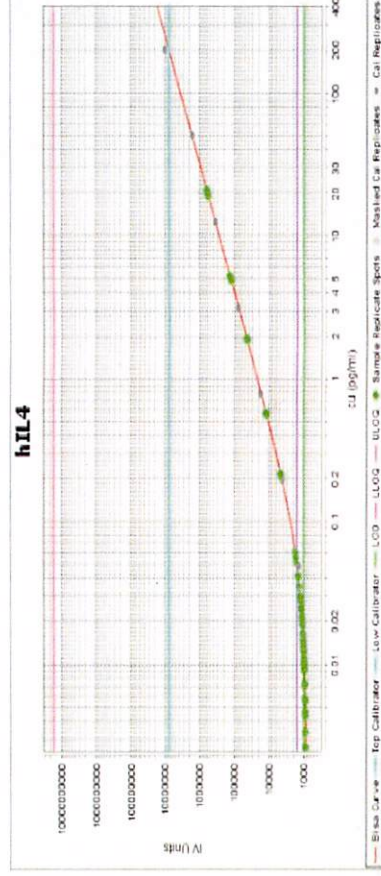
IL-1β Calibrator Curve



Calibrator Levels (pg/mL)

A	100.00
B	25.00
C	6.25
D	1.562
E	0.391
F	0.098
G	0.024
H	0.000

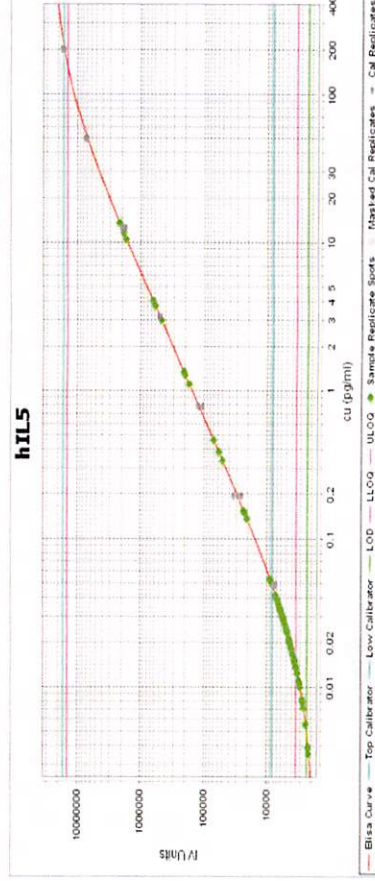
IL-4 Calibrator Curve



Calibrator Levels (pg/mL)

A	200.00
B	50.00
C	12.5
D	3.125
E	0.781
F	0.195
G	0.049
H	0.000

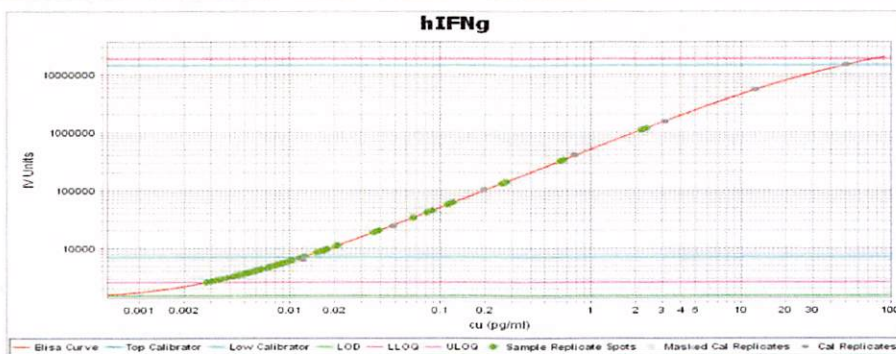
IL-5 Calibrator Curve



Calibrator Levels (pg/mL)

A	200.00
B	50.00
C	12.50
D	3.125
E	0.781
F	0.195
G	0.049
H	0.000

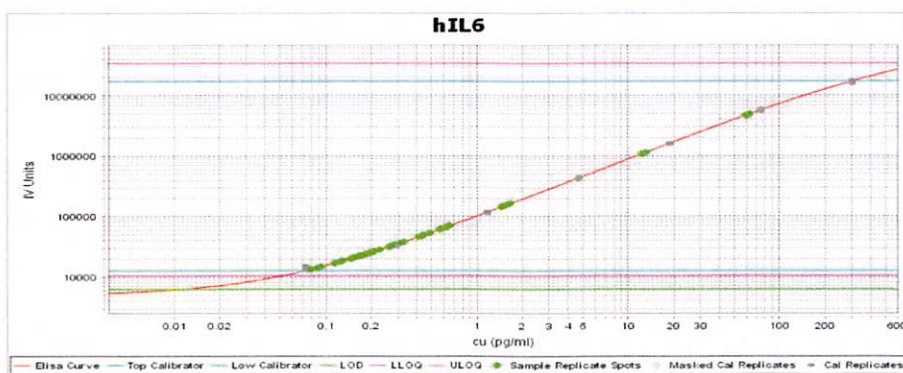
IFN γ Calibrator Curve



**Calibrator Levels
(pg/mL)**

A	50.00
B	12.50
C	3.125
D	0.781
E	0.195
F	0.049
G	0.012
H	0.000

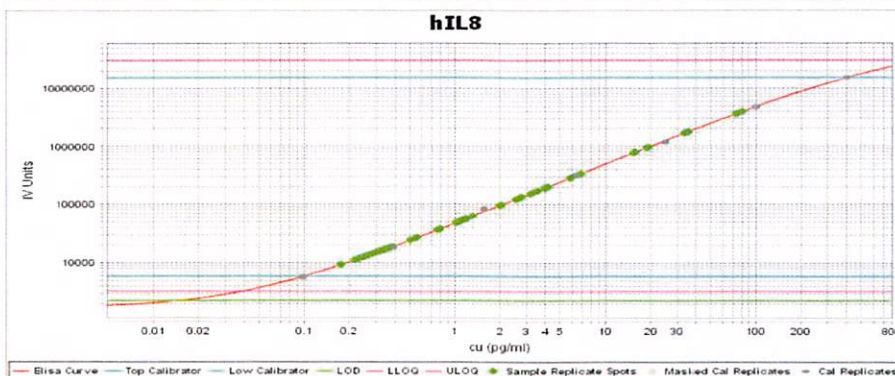
IL-6 Calibrator Curve



**Calibrator Levels
(pg/mL)**

A	300.00
B	75.00
C	18.75
D	4.688
E	1.172
F	0.293
G	0.073
H	0.000

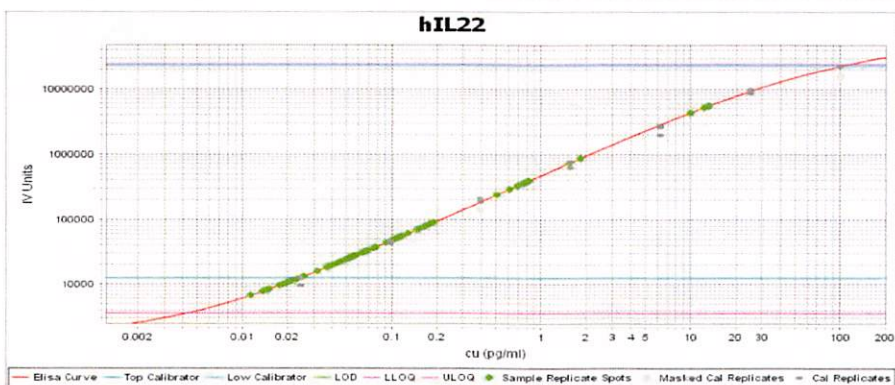
IL-8 Calibrator Curve



**Calibrator Levels
(pg/mL)**

A	400.00
B	100.00
C	25.00
D	6.25
E	1.562
F	0.391
G	0.098
H	0.000

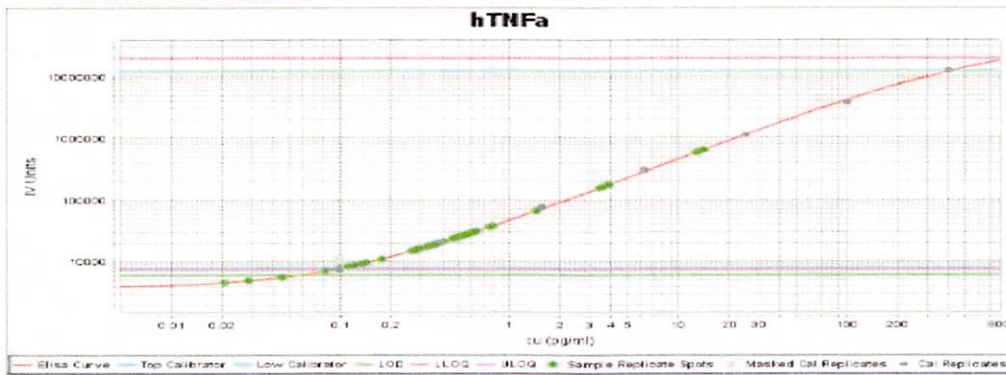
IL-22 Calibrator Curve



**Calibrator Levels
(pg/mL)**

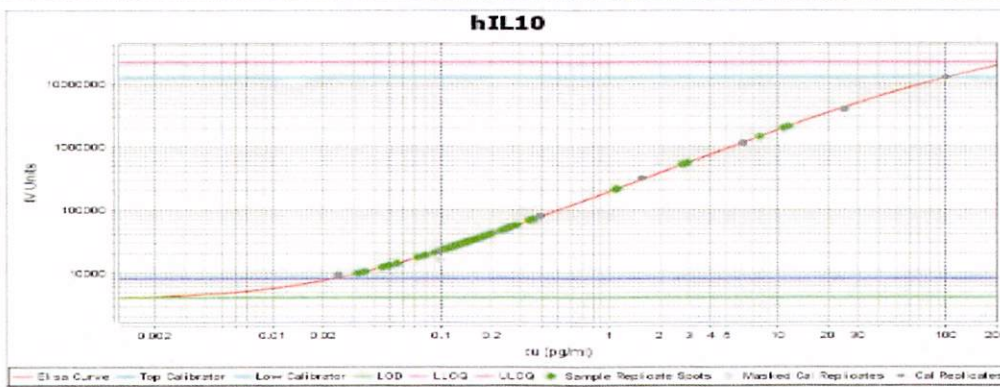
A	100.00
B	25.00
C	6.25
D	1.562
E	0.391
F	0.098
G	0.024
H	0.000

TNFα Calibrator Curve



Calibrator Levels (pg/mL)	
A	400.00
B	100.00
C	25.00
D	6.25
E	1.562
F	0.391
G	0.098
H	0.000

IL-10 Calibrator Curve



Calibrator Levels (pg/mL)	
A	100.00
B	25.00
C	6.25
D	1.562
E	0.391
F	0.098
G	0.024
H	0.000