



**Release Materials\***

**IFNgamma®**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	412221	0.743	0.538 - 0.999
Control 2 Concentrate	412222	34.4	27.3 - 50.7

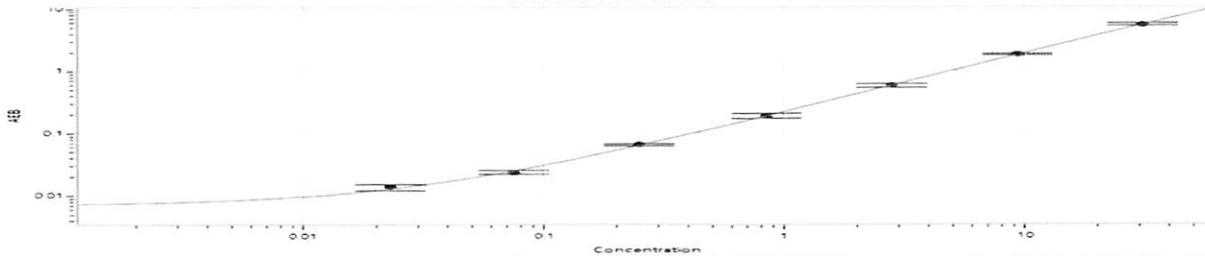
\* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibrator Concentrate Reconstitution Volume (uL)	1000
Control 1 Concentrate Reconstitution Volume (uL)	1000
Control 2 Concentrate Reconstitution Volume (uL)	1000

**Calibrator Levels (pg/mL)**

A	0.000
B	0.023
C	0.075
D	0.251
E	0.837
F	2.79
G	9.28
H	30.9

**Calibration Curve**



**Release Materials\***

**IL-6®**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	412221	1.01	0.744 - 1.38
Control 2 Concentrate	412222	48.2	37.0 - 68.7

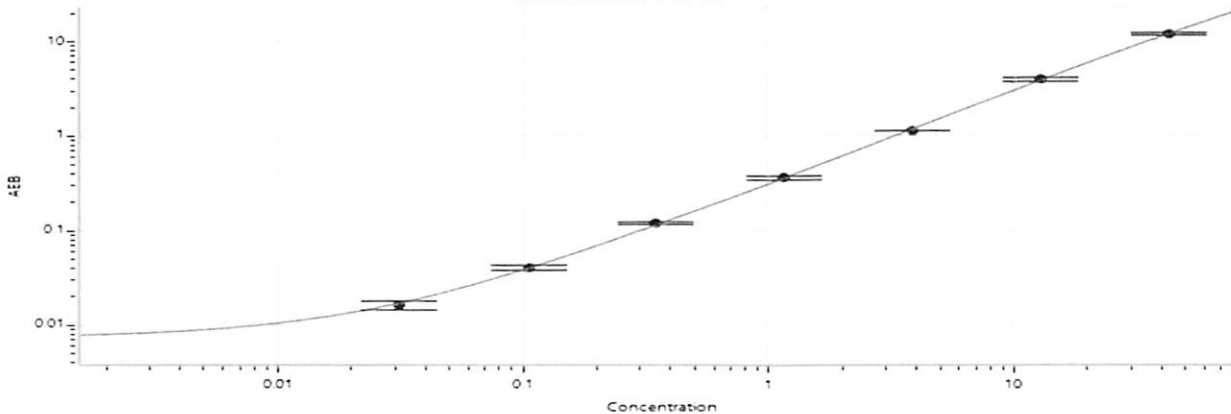
\* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibrator Concentrate Reconstitution Volume (uL)	1000
Control 1 Concentrate Reconstitution Volume (uL)	1000
Control 2 Concentrate Reconstitution Volume (uL)	1000

**Calibrator Levels (pg/mL)**

A	0.000
B	0.031
C	0.105
D	0.349
E	1.16
F	3.87
G	12.9
H	42.9

**Calibration Curve**



**Release Materials\***

**IL-8®**

	Lot	Result (pg/mL)	Spec. Range (pg/mL)
Control 1 Concentrate	412221	1.96	1.48 - 2.75
Control 2 Concentrate	412222	89.1	71.9 - 134

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

Calibrator Concentrate Reconstitution Volume (uL)	1000
Control 1 Concentrate Reconstitution Volume (uL)	1000
Control 2 Concentrate Reconstitution Volume (uL)	1000

**Calibrator Levels (pg/mL)**

A	0.000
B	0.063
C	0.210
D	0.699
E	2.33
F	7.75
G	25.8
H	85.9

**Calibration Curve**

