

## IFN- $\gamma$

### Interferon $\gamma$

#### human, recombinant, *E. coli*

| Cat. No. | Amount     |
|----------|------------|
| PR-420   | 50 $\mu$ g |

For *in vitro* use only  
Quality guaranteed for 12 months  
Store at -80°C

#### Avoid freeze / thaw cycles

#### Form

Liquid. Supplied in 20mM Tris-HCl pH 8.0, 5mM MgCl<sub>2</sub>, 10% glycerol and BSA as carrier.

#### Specific activity

10<sup>7</sup> units/mg.

#### Purity

≥95% by SDS-PAGE

#### Description

The cytokine Interferon  $\gamma$  (IFN- $\gamma$ ) plays an essential role in cell mediated immune responses to viral and mycobacterial infections. IFN- $\gamma$  has no viral specificity and inhibits the spread of viruses containing either RNA or DNA. Upon binding to macrophages IFN- $\gamma$  enables them to kill intracellular pathogens. It was also found that IFN- $\gamma$  exhibits antitumor properties.

#### Selected References:

- Meager *et al.* (1987) Quantification of interferons by anti-viral assays and their standardization. *Lymphokines and Interferons-A Practical Approach*. IRL, Oxford:129.
- Roilides *et al.* (2003) Immunomodulation of invasive fungal infections. *Infectious Disease Clinics of North America* **16**:193.
- Rotondi *et al.* (2003) Role for interferon-gamma inducible chemokines in endocrine autoimmunity: An expanding field. *Journal of Endocrinological Investigation* **26**:177.
- Agnello *et al.* (2003) Cytokines and transcription factors that regulate T helper cell differentiation: New players and new insights. *Journal of Clinical Immunology* **23**:147.