## IFN-γ (human), (recombinant)

Human interferon- $\gamma$  (IFN- $\gamma$ ) is an inflammatory cytokine secreted by T lymphocytes and natural killer (NK) cells. IFN- $\gamma$  is primarily characterized as an activator of macrophages, and is the principle cytokine produced by CD4+ inflammatory TH1 cells. IFN- $\gamma$  signals are transmitted via a heterodimeric IFNGR1/IFNGR2 receptor which activates transcription of immunoregulatory genes via the Jak/Stat pathway.

Citations: 4

View Online »

**Ordering Information** 

Order Online »

ENZ-PRT141-0100

100µg

Manuals, SDS & CofA

**View Online »** 

## **Handling & Storage**

**Handling** Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -80°C. For

long term storage, it is recommended to dilute to working aliquots in a 0.1% BSA

solution.

Short Term Storage -20°C

Long Term Storage -20°C

**Shipping** Ambient Temperature

## Regulatory Status RUO - Research Use Only

## **Product Details**

Alternative Name Interferon-γ

Endotoxin Content ≤1 EU/μg

**Formulation** Lyophilized from 10mM sodium phosphate, pH 7.4.

**MW** ~16.8 kDa

Purity ≥98% (SDS-PAGE gel & HPLC)

**Reconstitution** Reconstitute in sterile water at 1 mg/mL. Do not vortex.

**Source** Produced in *E.coli*.

UniProt ID P01579

Last modified: May 29, 2024

