

IFN- γ (human), (recombinant)

Human interferon- γ (IFN- γ) is an inflammatory cytokine secreted by T lymphocytes and natural killer (NK) cells. IFN- γ is primarily characterized as an activator of macrophages, and is the principle cytokine produced by CD4+ inflammatory TH1 cells. IFN- γ 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	100 μ g
-----------------	-------------

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/ μ g
Formulation	Lyophilized from 10mM sodium phosphate, pH 7.4.
MW	~16.8 kDa
Purity	$\geq 98\%$ (SDS-PAGE gel & HPLC)
Reconstitution	Reconstitute in sterile water at 1 mg/mL. Do not vortex.
Source	Produced in <i>E.coli</i> .
UniProt ID	P01579



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)