

IFN- γ (human), ELISA kit

**Ultrasensitive IFN-Gamma ELISA kit
measuring natural and recombinant human
IFN- γ , < 2 pg/ml**

This IFN-gamma ELISA kit is a colorimetric immunometric enzyme immunoassay kit with results in 3 hours. Absorbance is read at 450 nm.

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 »](#)

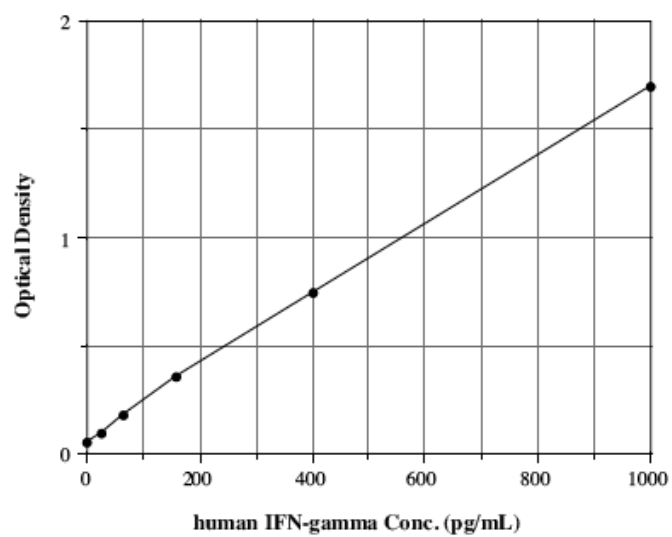
ADI-900-136

96 wells

Manuals, SDS & CofA

[View Online »](#)

- Ultrasensitive - quantify < 2 pg/ml
- Efficient - only 50 μ L diluted sample needed per well
- Specific - measure natural and recombinant human IFN- γ
- Results from up to 37 samples in duplicate in just 3 hours



IFN- γ (human), ELISA kit Kit graph

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Interferon-γ

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of human IFN-γ in culture supernatants, plasma, serum, and urine.

Assay Time 3 hours

Compatibility This product is compatible with the [Absorbance 96 Plate Reader](#).

Contents Microtiter plate, SA-HRP concentrate, SA-HRP dilution buffer, Biotinylated antibody, Standard, Standard diluent, Wash buffer concentrate, TMB Substrate, Stop solution

Sensitivity < 2 pg/ml (range 25.6 - 1,000 pg/ml)

Species Reactivity Human

Technical Info / Product Notes **Cited samples:**
[Cytokine ELISAs Cited Samples](#)

UniProt ID P01579

Wavelength 450 nm



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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