

VeriKine Human Interferon Gamma Receptor 1 ELISA Kit

The VeriKine Human IFN Gamma Receptor 1 ELISA quantitates Human IFN Gamma Receptor 1 (IFNGR1) in serum, plasma, tissue culture media (TCM) and cell lysates by sandwich enzyme linked immunosorbent assay (ELISA). IFNGR1 binds to plates coated with antibody and detection is accomplished using a detection antibody followed by streptavidin conjugated to horseradish peroxidase (HRP). This ELISA kit utilizes Tetramethyl-benzidine (TMB) as the substrate. The standard provided is recombinant Human IFNGR1 expressed in mammalian cells.

Ordering Information

[Order Online »](#)

PBL-41580-1	96 wells
-------------	----------

Manuals, SDS & CofA

[View Online »](#)

- Able to quantify soluble and membrane IFNGR1
- Compatible with serum, plasma, cell lysates and tissue culture media (TCM)
- Provides high sensitivity detection of Human IFNGR1 to < 10 pg/ml in Sample Diluent and Cell Lysates
- < 4 hour assay time
- Conserve high value specimens with low sample requirements

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name IFN- γ receptor, Interferon- γ receptor

Application ELISA

Application Notes Quantitative determination of Hu-IFN Gamma Receptor 1 (IFNGR1) in serum, plasma, tissue culture media and cell lysates

Assay Time Incubation time: 3 hours 15 minutes

Contents Pre-coated microtiter plate, Plate Sealers, Wash Solution Concentrate, Human IFN Gamma Receptor 1 (IFNGR1) Standard (500,000 pg/ml), Sample Diluent, Antibody Concentrate, Antibody Diluent, HRP Conjugate Concentrate, HRP Diluent, TMB Substrate, Stop Solution

Crossreactivity No measurable IFNGR1 was observed when using this kit with sera of cynomolgus, rhesus, pig, bovine, rabbit, mouse, sheep or donkey. No cross-reactivity was detected with mouse IFNGR1-Fc.

Sensitivity 9.38-300 pg/ml (sample diluent and cell lysates), 234.38-7500 pg/ml (serum and plasma), 37.5-1200 pg/ml (TCM)

Species Reactivity Human



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