

IL-31 (human), (recombinant)

Interleukin 31 (IL-31) is an immunoregulatory cytokine that is expressed by activated type 2 T helper (Th2) cells. IL-31 signals through a heterodimer receptor consisting of the IL-31 Receptor A (IL-31RA) and the oncostatin M receptor (OSMR). IL-31 promotes allergic reactions and inflammatory skin diseases.

Manuals, SDS & CofA

[View Online »](#)

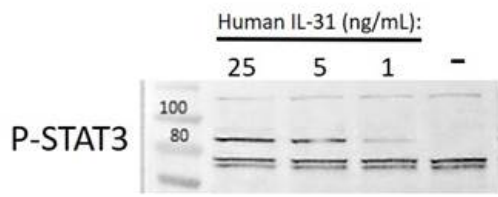


Figure 1: Human IL-31 induced activation of STAT in U-87MG cells.

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.
Long Term Storage	-80°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Interleukin-31
Endotoxin Content	≤1 EU/μg
Formulation	Lyophilized from a sterile (0.2 micron) filtered aqueous solution containing 10 mM sodium phosphate, 50 mM sodium chloride, pH 7.5.
MW	Monomer, 15.8 kDa (141 amino acids)
Purity	≥95% (SDS-PAGE)
Reconstitution	Reconstitute in sterile water at 0.1 mg/mL.
Source	Produced in <i>E.coli</i> .
UniProt ID	Q6EBC2

Last modified: May 29, 2024



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