

# 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.

## Ordering Information

[Order Online »](#)

ENZ-PRT261-0100	100µg
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)

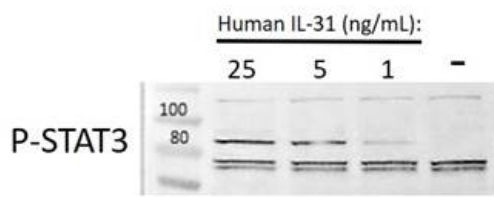


Figure 1: Human IL-31 induced activation of STAT in U-87MG cells.

## 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



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-be@enzolifesciences.com](mailto:info-be@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)