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 »

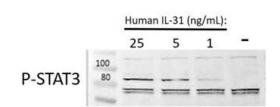


Figure 1: Human IL-31 induced activation of STAT in U-87MG cells.

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 *E.coli*.

UniProt ID Q6EBC2

Last modified: May 29, 2024

