

IL-5 (human), (recombinant)

Interleukin 5 (IL-5) is a hematopoietic growth factor that is expressed in type 2 T helper cells, mast cells, and eosinophils. IL-5 acts through the IL-5 receptor, stimulates B cell growth, and mediates eosinophil activation.

Manuals, SDS & CofA

[View Online »](#)

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-5, B-cell differentiation factor I, Eosinophil differentiaion factor, TFR
Biological Activity	TF-1 cell proliferation; ED50 \leq 2 ng/mL (\geq 5.0 x 10 ⁵ units/mg).
Endotoxin Content	\leq 1 EU/ μ g protein (LAL test)
Formulation	Lyophilized from a sterile (0.2 micron) filtered solution containing 10 mM sodium glycinate, pH 8.5.
MW	Dimer, ~13.3/26.6 kDa
Purity	\geq 95% (SDS-PAGE)
Reconstitution	Reconstitute in sterile water at 0.1 mg/mL.
Source	Produced in <i>E. coli</i> . Contains 116/232 amino acids.
UniProt ID	P05113

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