

# 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.

## Ordering Information

[Order Online »](#)

ENZ-PRT239-0010	10µg
ENZ-PRT239-0100	100µg

## 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 ≤ 2 ng/mL ( $\geq 5.0 \times 10^5$ units/mg).
Endotoxin Content	≤1 EU/μg 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	≥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



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