

# IL-9 (mouse), (recombinant)

Interleukin 9 (IL-9) is a cytokine produced by type 2 T helper cells and regulates hematopoietic cells. IL-9 functions to induce cell proliferation and prevent cell apoptosis. IL-9 signals through the interleukin 9 receptor to activate STAT signaling.

**Manuals, SDS & CofA**

[View Online »](#)

## Handling & Storage

<b>Handling</b>	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.
<b>Long Term Storage</b>	-80°C
<b>Shipping</b>	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Interleukin-9, T-cell growth factor p40, P40 cytokine
<b>Biological Activity</b>	M07e cell proliferation; $\leq 2$ ng/mL ( $\geq 5.0 \times 10^5$ units/mg).
<b>Endotoxin Content</b>	$\leq 1$ EU/ $\mu$ g protein (LAL test)
<b>Formulation</b>	Lyophilized from a sterile (0.2 micron) filtered solution containing 0.1% Trifluoroacetic Acid (TFA).
<b>MW</b>	~7.9 kDa
<b>Purity</b>	$\geq 90\%$ (SDS-PAGE)
<b>Reconstitution</b>	Reconstitute in sterile water at 0.1 mg/mL.
<b>Source</b>	Produced in <i>E. coli</i> . Contains 127 amino acids.
<b>UniProt ID</b>	P15247

Last modified: May 29, 2024



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)