

# IL-19 (mouse), (recombinant)

Interleukin-19 (IL-19) is produced by B cells and monocytes. IL-19 binds the interleukin 20 receptor complex to activate STAT3 signaling. IL-19 induces IL-6 and tumor necrosis factor alpha (TNF- $\alpha$ ) expression in monocytes, and promotes T helper 2 cell-mediated immune responses. IL-19 production is up-regulated in resting monocytes following granulocyte-macrophage colony-stimulating factor (GM-CSF) or lipopolysaccharide (LPS) stimulation.

## Ordering Information

[Order Online »](#)

ENZ-PRT246-0010	10 $\mu$ g
ENZ-PRT246-0100	100 $\mu$ g

## 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-19
Endotoxin Content	≤1 EU/μg protein (LAL test)
Formulation	Lyophilized from a sterile (0.2 micron) filtered solution containing 10 mM sodium phosphate and 50 mM sodium chloride, pH 7.5.
MW	~17.7 kDa
Purity	≥95% (SDS-PAGE)
Reconstitution	Reconstitute in sterile water at 0.1 mg/mL.
Source	Produced in <i>E. coli</i> . Contains 153 amino acids.
UniProt ID	Q8CJ70



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)