

# Macrophage inflammatory protein 2 (mouse), (recombinant)

Macrophage inflammatory protein 2 (MIP-2 or CXCL2), is a small cytokine secreted by monocytes and neutrophils at sites of inflammation. MIP-2 functions through CXCR2, a chemokine receptor, to act as a chemotactic agent for leukocytes and hematopoietic cells.

## Ordering Information

[Order Online »](#)

ENZ-PRT199-0020	
-----------------	--

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	Reconstituted solution is stable for 3 months when stored at -20°C to -80°C and for 1 month when stored at 4°C after reconstituting as directed.
<b>Handling</b>	Centrifuge vial before opening. Suspend the product by gently pipetting the above recommended solution and allow several minutes for complete reconstitution. Do not vortex. For prolonged storage, dilute to working aliquots in a 0.1% BSA solution, store at -80°C. Avoid freeze/thaw cycles.
<b>Short Term Storage</b>	-20°C
<b>Long Term Storage</b>	-80°C
<b>Shipping</b>	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	MIP-2, CXCL2
<b>Endotoxin Content</b>	≤1 EU/μg (LAL)
<b>Formulation</b>	Lyophilized from a sterile (0.2 micron) filtered aqueous solution containing 10 mM sodium phosphate, pH 7.5.
<b>MW</b>	Monomer, 7.9 kDa (73 amino acids)
<b>Purity</b>	≥95% (SDS-PAGE)
<b>Reconstitution</b>	Reconstitute with sterile water at 0.1 mg/mL.
<b>Sequence</b>	AVV ASELRCQCLK TLPRVDFKNI QSLSVTPPGP HCAQTEVIAT LKGGQKVCLD PEAPLVQKII QKILNKGKAN
<b>Source</b>	Produced in <i>E. coli</i> .
<b>UniProt ID</b>	P10889



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)