

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.

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	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.
Handling	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.
Short Term Storage	-20°C
Long Term Storage	-80°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

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

Last modified: May 29, 2024