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

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 *E. coli*.

UniProt ID P10889



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

Phone: +33 472 440 655 infofr@enzolifesciences.com

Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com