

GM-CSF (mouse), (recombinant)

Granulocyte-macrophage colony-stimulating factor (GM-CSF) is hematopoietic growth factor produced by endothelial cells, monocytes, fibroblasts, and T cells. GM-CSF stimulates the production of neutrophilic granulocytes, macrophages, and mixed granulocyte-macrophage colonies from bone marrow cells. GM-CSF promotes immune system development and regulates neutrophil function during infection.

Ordering Information

[Order Online »](#)

ENZ-PRT253-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

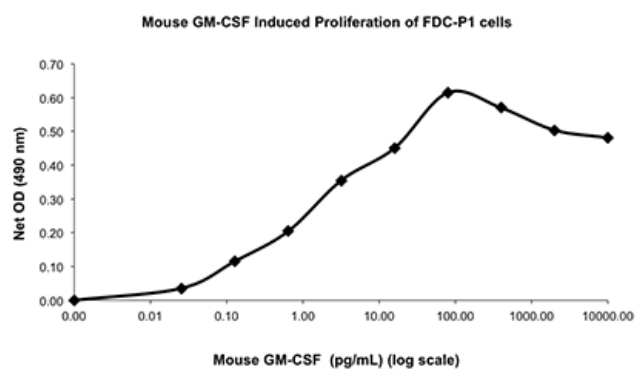


Figure 1: Serial dilutions of Mouse GM-CSF, starting at 1000 ng/mL, were added to FDC-P1 cells. After 93 hours, cell proliferation was measured and the linear portion of the curve can be used to calculate the ED50.

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	Granulocyte-macrophage colony stimulating factor, CSF-2, Pluripoietin- α , MGI1GM
Endotoxin Content	≤ 1 EU/ μ g protein (LAL test)
Formulation	Lyophilized from a sterile (0.2 micron) filtered aqueous solution containing 10 mM acetic acid.
MW	~ 14.3 kDa
Purity	$\geq 95\%$ (SDS-PAGE)
Reconstitution	Reconstitute in sterile water at 0.1 mg/mL.
Source	Produced in <i>E. coli</i> . Contains 125 amino acids.
UniProt ID	P01587



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com