GM-CSF (mouse), (recombinant)

Recombinant mouse GM-CSF that stimulates growth and differentiation of hematopoietic precursor cells.

The cytokine Granulocyte Macrophage-Colony Stimulating Factor (GM-CSF) controls the production, differentiation, and function of granulocytes and macrophages. The active form of GM-CSF functions extracellularly as a homodimer. Localized to a cluster of related genes at chromosome region 5q31, data links this gene to interstitial deletions in the 5q-syndrome and acute myelogenous leukemia. Other genes in the cluster include those encoding interleukins 4, 5, and 13. GM-CSF stimulates the growth and differentiation of hematopoietic precursor cells from various lineages, including granulocytes, macrophages, eosinophils and erythrocytes.

Ordering Information

Order Online »

ADI-908-064-005 5μg

Manuals, SDS & CofA

View Online »

Handling & Storage

Long Term Storage -20°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name CSF2, Colony stimulating factor 2 (granulocyte-

macrophage)

Formulation Lyophilized with no additives.

Purity ≥98% (RP-HPLC and SDS-PAGE)

Purity Detail Purified by multi-step chromatography.

Source Produced in E. coli.

UniProt ID P01587

Last modified: October 9, 2025

