

Recombinant Human M-CSF

Macrophage Colony Stimulating Factor, M-CSF, is a potent hematopoietic factor and regulator of cellular proliferation, differentiation, and survival for blood monocytes, tissue macrophages, and their respective progenitor cells. M-CSF has also been shown to play important roles in modulating dermal thickness and fertility, and is clinically used in the treatment of infection, malignancies and atherosclerosis.

Materials Provided

Catalog Number:	92-1295
Quantity Supplied:	1 vial(s), 10 µg per vial
Reconstitution Buffer:	1 mL

Description

DiscoverX control ligands have been functionally tested and validated for optimal use with all cell line targets and Assay Ready kits.

Product Information

Molecular Weight:	36800 g/mol
Source:	Recombinant
Purity:	≥ 98%
Endotoxin Level:	< 1EU/µg
Formulation:	N/A
Storage:	-20°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

To avoid condensation, equilibrate the vial to ambient temperature before opening.

Stock Concentration:	100 µg/mL
Reconstitution Volume:	0.100 mL
Reconstitution Solvent:	Reconstitution Buffer

For Research Use Only

Ordering: +1.510.979.1415 option 4 or e-mail CustomerServiceDRX@eurofins.com
Technical support: +1.510.979.1415 option 5 or e-mail DRX_SupportUS@eurofinsUS.com
General product information: www.discoverx.com