

Recombinant Human Oncostatin M

Oncostatin M is a pleiotropic cytokine that belongs to the interleukin 6 group of cytokines. This human protein functions in growth regulator; inhibits the proliferation of a number of tumor cell lines; stimulates proliferation of AIDS-KS cells; and regulates cytokine production, including IL-6, G-CSF and GM-CSF from endothelial cells.

Materials Provided

Catalog Number:	92-1294
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:	25700 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