

Instruction Manual

Catalog Number	D-64410
Description	Recombinant Murine Endothelial Cell Growth Factor (rMu VEGF) is a homodimeric protein secreted by a variety of vascularized tissues. The reported activities of VEGF include stimulation of endothelial cell growth, angiogenesis, and capillary permeability. Murine VEGF is a 39.036 kDa homodimeric, non-glycosylated protein consisting of two 165 amino acid polypeptide chains.
Quantity	10 µg
Molecular Mass	39 kDa
Specific Activity	see below
Source	E. coli
Purity	Greater than 95% (determined by SDS-PAGE, FPLC and HPLC analysis).
Endotoxin Level	< 0.1 ng per µg of rMu VEGF (1EU/µg).
Biological Activity	Recombinant murine VEGF165 is fully biologically active when compared to standards. The biological activity of PromoKine's murine VEGF was determined by the dose-dependent stimulation of the proliferation of human umbilical vein endothelial cells (HUVEC) using a concentration range of 1.0 to 7.0 ng/ml.
Formulation	Lyophilized from a sterile-filtered, concentrated (1 mg/ml) solution in PBS (pH 7.4) with no additives.
Reconstitution	Please Note: Always centrifuge product briefly before opening the vial. The lyophilized protein should be reconstituted in sterile, ultra-pure water to a concentration of 0.1 - 1.0 mg/ml. This solution can then be diluted into other aqueous buffers and stored at -20°C for future use.
Storage & Stability	The lyophilized rMu VEGF, though stable at room temperature for up to 3 weeks, is best stored desiccated at -20°C. Reconstituted rMu VEGF should be used immediately or stored long-term in undiluted working aliquots at -20°C. For long-term storage, it is recommended to add a carrier protein (0.1% endotoxin-free HSA or BSA; e.g. Cat.No. C-69500A). Avoid repeated freeze / thaw cycles.

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.