

Recombinant Human TGF- β 1 (HEK 293 derived)

Transforming growth factor beta 1 (TGF- β 1) is a cytokine peptide responsible for controlling proliferation, differentiation and other functions in cells. Also involved in hematopoiesis, wound healing and immune system control, endothelial differentiation, stimulating osteoblast bone formation, causing chemotaxis, and in the proliferation and differentiation of osteoblasts. Synthesized in many cell types, and binds many different receptors.

Materials Provided

Catalog Number:	92-1222
Quantity Supplied:	1 vial(s), 2 μ g 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:	25 kg/mol
Source:	Recombinant
Purity:	\geq 98% by HPLC
Endotoxin Level:	<0.01 EU per 1 μ g
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:	5 μ g/mL
Reconstitution Volume:	0.400 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