

Recombinant Human BMP - 7

Bone morphogenetic protein 7 (BMP-7) is a protein member of the TGFB superfamily that induces cartilage and bone formation and is expressed in the brain, kidneys and bladder. Has roles in calcium regulation, bone homeostasis, and epithelial osteogenesis.

Materials Provided

Catalog Number:	92-1225
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:	28800 g/mol
Source:	Recombinant
Purity:	≥ 98% by HPLC
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:	5 µM
Reconstitution Volume:	0.069 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