

Recombinant Human BMP-6

Bone morphogenetic protein 6 (BMP-6) is a signaling protein and member of the TGF β superfamily, responsible for inducing bone and cartilage growth, particularly prominent in early development. Induces the osteogenic markers in mesenchymal stem cells.

Materials Provided

Catalog Number:	92-1269
Quantity Supplied:	1 vial(s), 10 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:	26200 g/mol
Source:	Recombinant
Purity:	\geq 95%
Endotoxin Level:	< 1EU/ 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:	100 μ g/mL
Reconstitution Volume:	0.1 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