

Wnt3a protein

WNT3A is a signaling glycoprotein that is important for multiple developmental processes, including cell fate regulation and patterning during embryogenesis, and cell-cell signaling during morphogenesis of developing neural tube. Also implicated in oncogenesis and adipogenesis. Binds receptors of the Frizzled family of transmembrane receptors.

Materials Provided

Catalog Number:	92-1100
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:	37 kg/mol
Source:	Recombinant, Mouse
Purity:	≥ 75% by SDS-PAGE and HPLC
Endotoxin Level:	N/A
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:	60 µg/mL
Reconstitution Volume:	0.166 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