

Recombinant Human EGF

Epidermal growth factor (EGF) is a 53 amino acid polypeptide growth factor that binds EGF receptor (EGFR) and stimulates cell growth, proliferation, differentiation, and survival. It is a potent mitogenic factor, and defects result in hypomagnesemia type 4.

Materials Provided

Catalog Number:	92-1113
Quantity Supplied:	1 vial(s), 100 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:	6200 g/mol
Source:	Recombinant
Purity:	≥ 98% by SDS PAGE
Endotoxin Level:	< 0.1 ng/µg (<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:	500 µg/mL
Reconstitution Volume:	0.200 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