

Humanin, HN

Humanin is a 21 or 24 amino acid mitochondrial-derived peptide that has neuroprotective, cytoprotective, and anti-apoptotic effects. May be useful in protecting against alzheimer's disease, huntington's disease, prion disease, stroke, oxidative stress, atherosclerosis, type I diabetes, and heart attack.

Materials Provided

Catalog Number:	92-1221
Quantity Supplied:	1 vial(s), 1 mg per vial

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:	2687.3 g/mol
Source:	Synthetic
Purity:	≥ 95% by 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:	1 mM
Reconstitution Volume:	0.372 mL
Reconstitution Solvent:	H ₂ O

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