

Melanin Concentrating Hormone (6-17)

Melanin Concentrating Hormone (6-17) is a hypothalamic neuropeptide involved in feeding, mood, sleep-wake cycle, and energy balance. Overexpression is associated with obesity. Agonist of the melanin-concentrating hormone receptor (MCHR1).

Materials Provided

Catalog Number:	92-1173
Quantity Supplied:	1 vial(s), 500 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:	1537.88 g/mol
Source:	Synthetic
Purity:	≥ 97%
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:	0.325 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