

Melanotan II

Melanotan II is a synthetic analog of the melanocortin peptide hormone α -MSH, developed at the University of Arizona. Effects of its use include melanogenesis and aphrodesia. Not FDA approved, but Melanotan II products are currently sold and purchased online.

Materials Provided

Catalog Number:	92-1122
Quantity Supplied:	1 vial(s), 1 mg 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:	1024.18 g/mol
Source:	Synthetic
Purity:	$\geq 95\%$
Endotoxin Level:	N/A
Formulation:	N/A
Storage:	$\leq -20^{\circ}\text{C}$ Please avoid multiple freeze/thaw cycles.

Reagent Preparation

To avoid condensation, equilibrate the vial to ambient temperature before opening.

Stock Concentration:	2 mM
Reconstitution Volume:	0.488 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