

α MSH

α -Melanocyte stimulating hormone (α -MSH) is a pituitary hormone that plays a key role in appetite control and the skins UV protection and development of pigmentation. Also has effects on energy homeostasis, sexual activity, and as an anti-inflammatory. Potent agonist of the melanocortin receptors MC1, MC3, MC4, and MC5 ($K_i=0.12, 31, 660,$ and 5700nM , respectively), but not MC2.

Materials Provided

Catalog Number:	92-1172
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:	1664.88 g/mol
Source:	Synthetic
Purity:	$\geq 97\%$
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:	1 mg/mL
Reconstitution Volume:	1.0 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