

[D-Ala2, D-Leu5]-Enkephalin (DADLE)

[D-Ala2, D-Leu5]-Enkephalin (DADLE) is a synthetic peptide that is a selective agonist of the delta opioid receptor, but also binds mu opioid receptors of subtype $\mu 1$. It has antinociceptive and analgesic properties via neuron protection, and in-vivo effects include reduction in arterial blood pressure and heart rate.

Materials Provided

Catalog Number:	92-1055
Quantity Supplied:	1 vial(s), 5 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:	568.7 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:	10 mM
Reconstitution Volume:	0.879 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