

Galanin

Galanin is a neuropeptide that acts as a cellular messenger in the nervous system, regulating physiologic functions related to endocrine, metabolism, and behavior. Predominantly involved in inhibition of action potentials in neurons, and has been implicated in functions related to nociception, sleep, cognition, appetite, mood, and blood pressure.

Materials Provided

Catalog Number:	92-1107
Quantity Supplied:	1 vial(s), 0.1 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:	3157.41 g/mol
Source:	Synthetic
Purity:	≥ 97% 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:	500 µM
Reconstitution Volume:	0.063 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