

Tyramine

Tyramine is a naturally occurring monoamine derivative of tyrosine that acts as a catecholamine releasing agent. Has nonpsychoactive peripheral sympathomimetic effects, and can lead to a hypertensive crisis if foods rich in tyramine are ingested in conjunction with monoamine oxidase inhibitors. Occurs widely in foods as a result of fermentation or decay via tyrosine decarboxylation.

Materials Provided

Catalog Number:	92-1223
Quantity Supplied:	1 vial(s), 10 mg (re-packaged) 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:	137.18 g/mol
Source:	Synthetic
Purity:	≥ 98.5% by GC
Endotoxin Level:	N/A
Formulation:	N/A
Storage:	RT Please avoid multiple freeze/thaw cycles.

Reagent Preparation

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

Stock Concentration:	100 mM
Reconstitution Volume:	0.729 mL
Reconstitution Solvent:	DMSO

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