

Zaprinast

Zaprinast is a synthetic cGMP-specific phosphodiesterase (PDE) inhibitor from which sildenafil (Viagra) was developed. It is selective for PDE5, PDE6, PDE9 and PDE11 (IC50s of 0.76, 0.15, 29, and 12 μM , respectively), and is also a GPR35 agonist. It was an unsuccessful clinical drug candidate.

Materials Provided

Catalog Number:	92-1048
Quantity Supplied:	1 vial(s), 10 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:	271.28 g/mol
Source:	Synthetic
Purity:	$\geq 99\%$
Endotoxin Level:	N/A
Formulation:	N/A
Storage:	23-25°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

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

Stock Concentration:	20 mM
Reconstitution Volume:	1.84 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