

CHIR 99021 (CT99021)

CHIR-99021 HCl is a synthetic small molecule inhibitor of GSK3 kinase, with IC₅₀ of 10 and 6.7nM for GSK3 α and GSK3 β , respectively. Has >500-fold selectivity for GSK3 over closely related CDC2 and ERK2 kinases. Functions as a WNT activator.

Materials Provided

Catalog Number:	92-1116
Quantity Supplied:	1 vial(s), 2 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:	501.8 g/mol
Source:	Synthetic
Purity:	≥ 99%
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:	5 mM
Reconstitution Volume:	0.797 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