

(+)-CGP-12177A

(+)-CGP-12177A is a synthetic hydrophilic inhibitor of the adrenergic receptors β_1 , β_2 , and β_3 ($K_i=0.9$, 4 , and 88 nM, respectively). Inhibits isoproterenol stimulation of adenylyl cyclase activity at low concentrations. Can also act as an agonist of these receptors at higher concentrations.

Materials Provided

Catalog Number:	92-1216
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:	315.8 g/mol
Source:	Synthetic
Purity:	$\geq 98\%$ by HPLC
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:	0.316 mL
Reconstitution Solvent:	H ₂ O

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