

Thapsigargin

Thapsigargin is a naturally occurring, cell-permeable plant sesquiterpene lactone that non-competitively inhibits the sarco/endoplasmic reticulum calcium ATPase (SERCA, IC₅₀ ~2-80nM). Effectively increases intracellular calcium levels upon SERCA inhibition, leading to cell activation, increased histamine release from mast cells, and increased cancer cell proliferation. Acts as an anti-inflammatory and induces autophagy in-vivo.

Materials Provided

Catalog Number:	92-1148
Quantity Supplied:	1 vial(s), 1 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:	650.76 g/mol
Source:	Synthetic
Purity:	≥ 97%
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:	0.307 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