

SB220025

SB220025 is a synthetic small molecule potent and selective inhibitor of the p38 mitogen-activated protein kinase (MAPK, IC₅₀=60nM), 50-1000-fold selective over other MAP kinases. Potent inhibitor of angiogenesis, inflammatory cytokine production, and TNF- α production.

Materials Provided

Catalog Number:	92-1241
Quantity Supplied:	1 vial(s), 0.5 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:	447.76 g/mol
Source:	Synthetic
Purity:	$\geq 98\%$
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:	5 mM
Reconstitution Volume:	0.223 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