

SB203580 4(4-Fluorophenyl)-2-(4-methylsulfinylphenyl)-5-(4-pyridyl) 1 H-imidazole

Publication Number MAN0003446

Rev. 2.00








Catalog Number:	PHZ1253
Quantity:	1.0 mg
Lot Number:	See product label.
Appearance:	White solid
Molecular Formula:	C ₂₁ H ₁₆ N ₃ FOS
Molecular Weight:	377.4
Purity:	98%, as assessed by TLC.
Summary:	SB203580 is a selective inhibitor of p38 mitogen-activated protein (MAP) kinase, which acts by competitively inhibiting ATP binding. This compound inhibits p38 MAPK stimulation of MAPKAPK2 (IC ₅₀ = 70 nM) and is a useful tool for assessing the physiological roles of p38.
Biological Activity:	K _i = 21 nM, determined by measuring competitive inhibition of ATP binding.
Solubility:	Soluble in DMSO at a concentration of 50 mg/mL.
Sterility:	This product is not sterile.
Storage:	Store, as supplied, at -20°C. Upon solubilization, apportion into working aliquots and store at -20°C. Avoid repeated freeze/thaw cycles. Solutions are stable at -20°C for up to three months.
Expiration Date:	Expires one year from date of receipt when stored as instructed.
References:	<p>Cuenda, A., et al. (1995) SB203580 is a specific inhibitor of a MAP kinase homologue which is stimulated by cellular stresses and interleukin-1. <i>FEBS Lett.</i> 364(2) 229-233.</p> <p>Young, P.R., et al. (1997) Pyridinyl imidazole inhibitors of p38 mitogen-activated protein kinase bind in the ATP site. <i>J. Biol. Chem.</i> 272(18)12116-12121.</p> <p>Ko, B.C., et al. (2002) Fyn and p38 signaling are both required for maximal hypertonic activation of the osmotic response element-binding protein/tonicity-responsive enhancer-binding protein (OREBP/TonEBP). <i>J. Biol. Chem.</i> 277(48):46085-46092.</p> <p>Kogut, M., et. al. (2002) Selective pharmacological inhibitors reveal the role of Syk tyrosine kinase, phospholipase C, phosphatidylinositol-3'-kinase, and p38 mitogen-activated protein kinase in Fc receptor-mediated signaling of chicken herophil degranulation. <i>Int. Immunopharmacol.</i> 2(7):963-973.</p> <p>Knebel, A., et al. (2002) Stress-induced regulation of eukaryotic elongation factor 2 kinase by SB203580-sensitive and -insensitive pathways. <i>Biochem. J.</i> 367(Pt 2):525-532.</p>







Related Products:

Product	Catalog no.
p38 [pTpY ^{180/182}] antibody	44-684
MEK3/6 [pSpT ^{189/193}]/[pSpT ^{207/211}] antibody	44-470

Explanation of Symbols

The symbols present on the product label are explained below:

Symbol	Description
	Catalog Number
	Research Use Only
	Use by
	Manufacturer
	Without, does not contain
	Protect from light
	Directs the user to consult instructions for use (IFU), accompanying the product.

Symbol	Description
	Batch code
	In vitro diagnostic medical device
	Temperature limitation
	European Community authorized representative
	With, contains
	Consult accompanying documents

Limited Use Label License: Research Use Only

The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial applications of any kind, including, without limitation, quality control and commercial services such as reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact outlicensing@lifetech.com or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.

For Research Use Only. Caution: Not for human or animal therapeutic or diagnostic use.

Manufacturing site: 7335 Executive Way | Frederick, MD 21704 | Toll Free in USA 800.955.6288

© 2011 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.

For support visit www.lifetechnologies.com/support or email techsupport@lifetech.com

www.lifetechnologies.com

Revision Date 13 May 2011

 life
technologies™