

SB203580

p38 MAP Kinase Inhibitor

Catalog # tlr1-sb20

For research use only

Version # 15K04-MM

PRODUCT INFORMATION

Content:

- 5 mg SB203580

Storage and stability:

- SB203580 is provided as a translucent film and shipped at room temperature. Store at -20°C in the dark. Solid product is stable 1 year at -20°C.

- Upon resuspension, prepare aliquots of SB203580 and store at -20°C. Avoid freeze-thaw cycles. Protect from light. Resuspended product is stable 3 month at -20°C when properly stored.

Quality control:

- Purity: ≥98% (UHPLC)
- The absence of bacterial contamination (e.g. lipoproteins and endotoxins) has been confirmed using HEK-Blue™ TLR2 and HEK-Blue™ TLR4 cells.

DESCRIPTION

SB203580 is a pyridinyl imidazole inhibitor widely used to elucidate the roles of p38 mitogen-activated protein kinase (MAPK)¹. It inhibits p38 catalytic activity by binding to the ATP binding pocket. SB203580 also inhibits the phosphorylation and activation of protein kinase B (PKB, also known as Akt)². Both kinases are involved in a wide array of signaling pathways, including the TLR signaling pathway³. Moreover, several studies suggest that p38 MAPKs regulate distinct phases of autophagy. p38 can elicit autophagy via Beclin1. Contrarily, p38 α has also been reported to inhibit autophagy by interfering with the trafficking of Atg9. SB203580 is generally used at low concentrations (1-10 μ M), as it has been reported that at high concentrations SB203580 (> 20 μ M) can induce the activation of the serine/threonine kinase Raf-1⁴.

CHEMICAL PROPERTIES

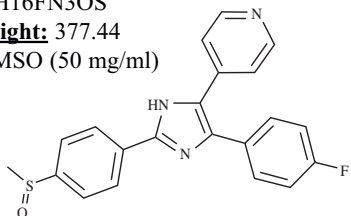
Synonym: 4-(4'-Fluorophenyl)-2-(4'-methylsulfinylphenyl)-5-(4'-pyridyl)-imidazole

Formula: C₆H₁₆FN₃O₃S

Molecular weight: 377.44

Solubility: DMSO (50 mg/ml)

Structure:



METHODS

Preparation of sterile stock solution (25 mM)

1. Add 530 μ l DMSO to 5 mg of SB203580.
2. Vortex until complete solubilization.
3. Aliquote and store at -20°C.
4. Once SB203580 is solubilized, dilutions can be prepared using aqueous buffers.

Working concentration: 1-10 μ M

PROTOCOLS

For reference only; as described in the indicated publications.

Cell Culture Assay³

Cells: Mouse bone-marrow derived immature dendritic cells

Working concentration: 1-10 μ M

Incubation time: 1-2 h

Methods: RT-PCR and Western blot

Cell Culture Assay⁵

Cells: Primary cardiac fibroblasts

Working concentration: 10 μ M

Incubation time: 45 min

Methods: Western blot and assays to measure collagen synthesis

1. **Cuenda A. et al., 1995.** SB 203580 is a specific inhibitor of a MAP kinase homologue which is stimulated by cellular stresses and interleukin-1. FEBS Lett. 364:229-233. 2. **Lali FV. et al., 2000.** The pyridinyl imidazole inhibitor SB203580 blocks phosphoinositide-dependent protein kinase activity, protein kinase B phosphorylation, and retinoblastoma hyperphosphorylation in interleukin-2-stimulated T cells independently of p38 mitogen-activated protein kinase. J Biol Chem. 275(10):7395-402. 3. **Jarnicki AG. et al., 2008.** Attenuating Regulatory T Cell Induction by TLR Agonists through Inhibition of p38 MAPK Signaling in Dendritic Cells Enhances Their Efficacy as Vaccine Adjuvants and Cancer Immunotherapeutics. J. Immunol., 180: 3797 - 3806. 4. **Kalmes A et al., 1999.** Raf-1 is activated by the p38 mitogen-activated protein kinase inhibitor, SB203580. FEBS Lett. 444(1):71-4. 5. **Mir SA. et al., 2012.** Inhibition of signal transducer and activator of transcription 3 (STAT3) attenuates interleukin-6 (IL-6)-induced collagen synthesis and resultant hypertrophy in rat heart. J Biol Chem. 287(4):2666-77.

RELATED PRODUCTS

Product	Description	Cat. Code
AZD6244	MEK1/2 Inhibitor	inh-ad62
PD0325901	MEK Inhibitor	inh-pd32
SB202190	MAP Kinase Inhibitor	tlr1-sb90

TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Hong Kong: +852 3-622-34-80

E-mail: info@invivogen.com

 **InvivoGen**
www.invivogen.com