

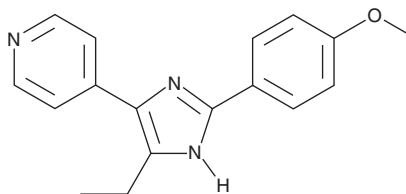
PRODUCT INFORMATION



SB 202474

Item No. 18749

CAS Registry No.: 172747-50-1
Formal Name: 4-[4-ethyl-2-(4-methoxyphenyl)-1H-imidazol-5-yl]-pyridine
MF: C₁₇H₁₇N₃O
FW: 279.3
Purity: ≥97%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Special Conditions: Protect from light



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

SB 202474 is supplied as a solid. A stock solution may be made by dissolving the SB 202474 in the solvent of choice, which should be purged with an inert gas. SB 202474 is soluble in organic solvents such as methanol and DMSO.

Description

SB 202190 (Item No. 10010399) and SB 203580 (Item No. 13067) are potent, selective inhibitors of the MAP kinases p38 α and p38 β .^{1,2} SB 202474 is a structural analog of SB 202190 and SB 203580 that is used as a negative control in studies of p38 inhibition.³⁻⁵

References

1. Jiang, Y., Chen, C., Li, Z., *et al.* Characterization of the structure and function of a new mitogen-activated protein kinase (p38 β). *J. Biol. Chem.* **271**(30), 17920-17926 (1996).
2. Davies, S.P., Reddy, H., Caivano, M., *et al.* Specificity and mechanism of action of some commonly used protein kinase inhibitors. *Biochem. J.* **351**(1), 95-105 (2000).
3. Lee, J.C., Laydon, J.T., McDonnell, P.C., *et al.* A protein kinase involved in the regulation of inflammatory cytokine biosynthesis. *Nature* **372**(6508), 739-746 (1994).
4. Sweeney, G., Somwar, R., Ramlal, T., *et al.* An inhibitor of p38 mitogen-activated protein kinase prevents insulin-stimulated glucose transport but not glucose transporter translocation in 3T3-L1 adipocytes and L6 myotubes. *J. Biol. Chem.* **274**(15), 10071-10078 (1999).
5. Chang, A.Y.W. Pro-life role for c-Jun N-terminal kinase and p38 mitogen-activated protein kinase at rostral ventrolateral medulla in experimental brain stem death. *J. Biomed. Sci.* **19**(1), (2012).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/02/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM