

# PRODUCT INFORMATION



## JX-401

Item No. 16898

CAS Registry No.: 349087-34-9

Formal Name: [2-methoxy-4-(methylthio)phenyl]  
[4-(phenylmethyl)-1-piperidinyl]-methanone

Synonyms: p38 MAPK Inhibitor VI,  
p38 Mitogen-activated Protein Kinase Inhibitor VI

MF: C<sub>21</sub>H<sub>25</sub>NO<sub>2</sub>S

FW: 355.5

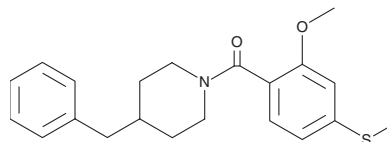
Purity: ≥95%

UV/Vis.: λ<sub>max</sub>: 210, 265 nm

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥4 years



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

## Laboratory Procedures

JX-401 is supplied as a crystalline solid. A stock solution may be made by dissolving the JX-401 in the solvent of choice. JX-401 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of JX-401 in these solvents is approximately 1, 10, and 20 mg/ml, respectively.

JX-401 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, JX-401 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. JX-401 has a solubility of approximately 0.1 mg/ml in a 1:5 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

JX-401 is a potent, reversible inhibitor of the p38α isoform of MAP kinase (IC<sub>50</sub> = 32 nM).<sup>1</sup> It does not inhibit the p38γ isoform.<sup>1</sup> JX-401 is effective in mammalian cells, as it blocks the differentiation of myoblasts into myotubes.<sup>1</sup>

## Reference

1. Friedmann, Y., Shriki, A., Bennett, E.R., *et al.* JX401, a p38α inhibitor containing a 4-benzylpiperidine motif, identified *via* a novel screening system in yeast. *Mol. Pharmacol.* **70**(4), 1395-1405 (2006).

### 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, 11/30/2022

## 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