

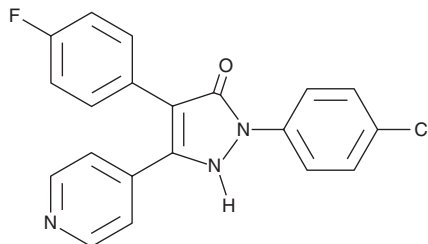
# PRODUCT INFORMATION



## p38 MAPK Inhibitor

Item No. 17650

**CAS Registry No.:** 219138-24-6  
**Formal Name:** 2-(4-chlorophenyl)-4-(4-fluorophenyl)-1,2-dihydro-5-(4-pyridinyl)-3H-pyrazol-3-one  
**Synonyms:** p38 MAP Kinase Inhibitor, p38 Mitogen-activated Protein Kinase Inhibitor  
**MF:** C<sub>20</sub>H<sub>13</sub>ClFN<sub>3</sub>O  
**FW:** 365.8  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 260, 282 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

p38 MAPK inhibitor is supplied as a crystalline solid. A stock solution may be made by dissolving the p38 MAPK inhibitor in the solvent of choice. p38 MAPK inhibitor is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of p38 MAPK inhibitor in ethanol is approximately 0.3 mg/ml and approximately 5 mg/ml in DMSO and DMF.

p38 MAPK inhibitor is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, p38 MAPK inhibitor should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. p38 MAPK inhibitor has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

p38 MAPK inhibitor is a potent inhibitor of p38 MAP kinase (IC<sub>50</sub> = 35 nM).<sup>1</sup> It inhibits senescence induced by the oncogene RAS.<sup>2</sup>

### References

1. de Laszlo, S.E., Visco, D., Agarwal, L., *et al.* Pyrroles and other heterocycles as inhibitors of p38 kinase. *Bioorg. Med. Chem. Lett.* **8**(19), 2689-2694 (1998).
2. Zhang, R., Bitler, B.J., and Peterson, J. Methods and compositions for screening modulators of oncogene-induced senescence. United States Patent Application Publication US20140272947A1 (2014), 1-10, US 14/202,909.

#### 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, 11/02/2022

#### 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