

# PRODUCT INFORMATION



**CP 424,174**

Item No. 29675

**CAS Registry No.:** 210825-31-3

**Formal Name:** N-[[[4-chloro-2,6-bis(1-methylethyl)phenyl]amino]carbonyl]-3-(1-hydroxy-1-methylethyl)benzenesulfonamide

**MF:** C<sub>22</sub>H<sub>29</sub>ClN<sub>2</sub>O<sub>4</sub>S

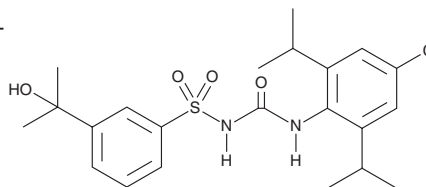
**FW:** 453.0

**Purity:** ≥98%

**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

CP 424,174 is supplied as a crystalline solid. A stock solution may be made by dissolving the CP 424,174 in the solvent of choice, which should be purged with an inert gas. CP 424,174 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of CP 424,174 in these solvents is approximately 30 mg/ml.

CP 424,174 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CP 424,174 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. CP 424,174 has a solubility of approximately 0.16 mg/ml in a 1:5 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

CP 424,174 is an inhibitor of stimulus-coupled IL-1 $\beta$  post-translational processing.<sup>1</sup> It inhibits ATP-induced proteolytic cleavage of pro-IL-1 $\beta$ , measured by secretion of the 17 kDa mature form of IL-1 $\beta$ , in LPS-primed human blood-derived monocytes when used at concentrations of 0.5 and 2.5  $\mu$ M. CP 424,174 inhibits ATP-induced IL-1 $\alpha$  and IL-1 $\beta$  production in LPS-primed mouse peritoneal macrophages (ED<sub>50</sub>s = 15 and 14 mg/kg, respectively).

## Reference

1. Perregaux, D.G., McNiff, P., Lalibertie, R., *et al.* Identification and characterization of a novel class of interleukin-1 post-translational processing inhibitors. *J. Pharmacol. Exp. Ther.* **299**(1), 187-197 (2001).

### 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