

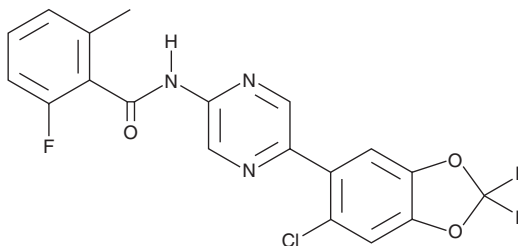
# PRODUCT INFORMATION



## CM-4620

Item No. 41977

**CAS Registry No.:** 1713240-67-5  
**Formal Name:** N-[5-(6-chloro-2,2-difluoro-1,3-benzodioxol-5-yl)-2-pyrazinyl]-2-fluoro-6-methyl-benzamide  
**Synonyms:** Zegrocraetin  
**MF:** C<sub>19</sub>H<sub>11</sub>ClF<sub>3</sub>N<sub>3</sub>O<sub>3</sub>  
**FW:** 421.8  
**Purity:** ≥98%  
**Supplied as:** A 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

CM-4620 is supplied as a solid. A stock solution may be made by dissolving the CM-4620 in the solvent of choice, which should be purged with an inert gas. CM-4620 is sparingly soluble (1-10 mg/ml) in acetonitrile and DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

## Description

CM-4620 is an inhibitor of Orai1/STIM1-containing store-operated calcium channels (SOCs; IC<sub>50</sub> = 120 nM).<sup>1</sup> It is selective for Orai1/STIM1-containing SOCs over Orai2/STIM2-containing SOCs (IC<sub>50</sub> = 900 nM). CM-4620 (3 μM) decreases cholecystokinin- or tauro lithocholic acid 3-sulfate-induced cell death in isolated mouse, rat, and human acini. *In vivo*, CM-4620 (20 mg/kg) reduces blood amylase and lipase levels and pancreatic trypsin activity in a mouse model of cerulein-induced acute pancreatitis.

## Reference

1. Waldron, R.T., Chen, Y., Pham, H., et al. The Orai Ca<sup>2+</sup> channel inhibitor CM4620 targets both parenchymal and immune cells to reduce inflammation in experimental acute pancreatitis. *J. Physiol.* **597(12)**, 3085-3105 (2019).

### 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, 10/31/2024

## 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