

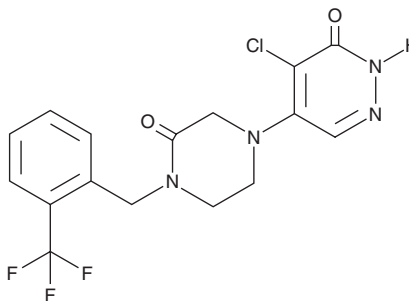
PRODUCT INFORMATION



GFB-8438

Item No. 43035

CAS Registry No.: 2304549-73-1
Formal Name: 4-chloro-5-[3-oxo-4-[[2-(trifluoromethyl)phenyl]methyl]-1-piperazinyl]-3(2H)-pyridazinone
MF: C₁₆H₁₄ClF₃N₄O₂
FW: 386.8
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

GFB-8438 is supplied as a solid. A stock solution may be made by dissolving the GFB-8438 in the solvent of choice, which should be purged with an inert gas. GFB-8438 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

GFB-8438 is an inhibitor of transient receptor potential canonical 5 (TRPC5; IC₅₀ = 0.18 μM).¹ It is selective for TRPC5 over TRPC6 and voltage-gated sodium channel 1.5 (Na_v1.5; IC₅₀s = >30 μM for both) and a panel of 100 receptors and kinases at 10 μM but does inhibit TRPC4 (IC₅₀ = 0.29 μM). GFB-8438 (30 mg/kg per day) decreases urinary excretion of total protein and albumin in a rat model of high salt- and unilateral nephrectomy-induced renal hypertension.

Reference

1. Yu, M., Ledebor, M.W., Daniels, M., *et al.* Discovery of a potent and selective TRPC5 inhibitor, efficacious in a focal segmental glomerulosclerosis model. *ACS Med. Chem. Lett.* **10(11)**, 1579-1585 (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, 04/09/2025

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