

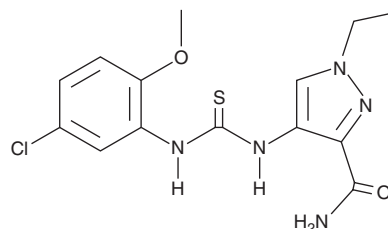
PRODUCT INFORMATION



BRD-9424

Item No. 29013

CAS Registry No.: 957485-64-2
Formal Name: 4-[[[(5-chloro-2-methoxyphenyl)amino]thioxomethyl]amino]-1-ethyl-1H-pyrazole-3-carboxamide
Synonyms: FPH2, Functional Proliferation Hit 2
MF: C₁₄H₁₆ClN₅O₂S
FW: 353.8
Purity: ≥98%
UV/Vis.: λ_{max}: 275 nm
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

BRD-9424 is supplied as a solid. A stock solution may be made by dissolving the BRD-9424 in the solvent of choice, which should be purged with an inert gas. BRD-9424 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of BRD-9424 in these solvents is approximately 0.25, 10, and 15 mg/ml, respectively.

BRD-9424 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BRD-9424 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. BRD-9424 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

BRD-9424, also known as functional proliferation hit 2 (FPH2), is a small molecule that induces proliferation of primary human hepatocytes *in vitro*.¹

Reference

1. Shan, J., Schwartz, R.E., Ross, N.T., *et al.* Identification of small molecules for human hepatocyte expansion and iPS differentiation. *Nat. Chem. Biol.* **9**(8), 514-520 (2013).

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, 12/06/2022

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