

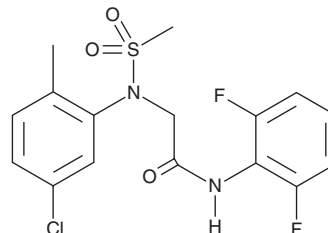
PRODUCT INFORMATION



BRD6125

Item No. 21804

CAS Registry No.: 708219-39-0
Formal Name: 2-[(5-chloro-2-methylphenyl)(methylsulfonyl)amino]-N-(2,6-difluorophenyl)-acetamide
Synonym: FPH1
MF: C₁₆H₁₅ClF₂N₂O₃S
FW: 388.8
Purity: ≥98%
UV/Vis.: λ_{max}: 261 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

BRD6125 is supplied as a crystalline solid. A stock solution may be made by dissolving the BRD6125 in the solvent of choice, which should be purged with an inert gas. BRD6125 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of BRD6125 in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

BRD6125 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BRD6125 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. BRD6125 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

BRD6125, also known as functional proliferation hits 1 (FPH1), is a small molecule that increases the number and activity of hepatocytes *in vitro*.¹ It dramatically increases expansion of primary human hepatocytes and induces maturation of hepatocyte-like cells (iHeps). It also induces maturation in two pluripotent cell lines. This product is also available as part of Cayman's Hepatocyte Differentiation Reagent Kit (Item No. 44737).

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, 10/27/2025

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