

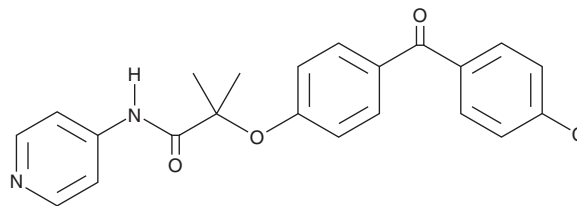
PRODUCT INFORMATION



HR68

Item No. 39855

CAS Registry No.: 2454582-64-8
Formal Name: 2-[4-(4-chlorobenzoyl)phenoxy]-2-methyl-N-4-pyridinyl-propanamide
Synonym: PP21
MF: C₂₂H₁₉ClN₂O₃
FW: 394.9
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥3 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

HR68 is supplied as a solid. A stock solution may be made by dissolving the HR68 in the solvent of choice, which should be purged with an inert gas. HR68 is soluble in methanol, chloroform, and acetonitrile.

Description

HR68 is an anticancer agent and a derivative of the peroxisome proliferator-activated receptor (PPAR) agonist fenofibric acid (Item No. 19262).¹ It decreases the viability of LN-229 glioblastoma cells (IC₅₀ = 1.17 μM). HR68 crosses the blood-brain barrier in a temozolomide-resistant orthotopic patient-derived xenograft (PDX) mouse model of glioblastoma.

Reference

- Ingraham, C.H., 4th., Stalinska, J., Carson, S.C., *et al.* Computational modeling and synthesis of pyridine variants of benzoyl-phenoxy-acetamide with high glioblastoma cytotoxicity and brain tumor penetration. *Sci. Rep.* **13**(1), 12235 (2023).

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, 11/17/2023

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