

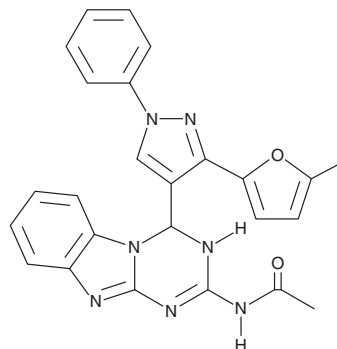
PRODUCT INFORMATION



PRGL493

Item No. 32748

CAS Registry No.: 2479378-45-3
Formal Name: N-(4-(3-(5-methylfuran-2-yl)-1-phenyl-1H-pyrazol-4-yl)-3,4-dihydrobenzo[4,5]imidazo[1,2-a][1,3,5]triazin-2-yl)acetamide
MF: C₂₅H₂₁N₇O₂
FW: 451.5
Purity: ≥98%
UV/Vis.: λ_{max}: 305 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

PRGL493 is supplied as a crystalline solid. A stock solution may be made by dissolving the PRGL493 in the solvent of choice, which should be purged with an inert gas. PRGL493 is soluble in the organic solvent chloroform at a concentration of approximately 1 mg/ml.

Description

PRGL493 is an inhibitor of long-chain acyl-CoA synthetase 4 (ACSL4).¹ It inhibits formation of arachidonoyl-CoA (AA-CoA) from arachidonic acid (Item Nos. 90010 | 90010.1 | 10006607) in, as well as the proliferation and migration of, PC3 and MDA-MB-231 cancer cells when used at a concentration of 50 μM. PRGL493 (5 μM) reduces production of progesterone (Item No. 15876) induced by 8-Br-cyclic AMP (8-Br-cAMP) in MA-10 Leydig tumor cells. It inhibits tumor growth in a PC3 mouse xenograft model when administered at a dose of 0.25 mg/kg.

Reference

1. Castillo, A.F., Orlando, U.D., Maloberti, P.M., et al. New inhibitor targeting Acyl-CoA synthetase 4 reduces breast and prostate tumor growth, therapeutic resistance and steroidogenesis. *Cell. Mol. Life Sci.* (2020).

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, 03/26/2024

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