

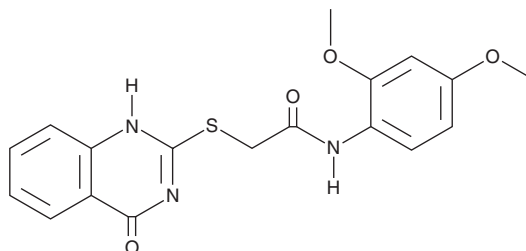
PRODUCT INFORMATION



Z16078526

Item No. 43072

CAS Registry No.: 852222-94-7
Formal Name: 2-[(3,4-dihydro-4-oxo-2-quinazolinyl)thio]-N-(2,4-dimethoxyphenyl)-acetamide
Synonym: Z160
MF: C₁₈H₁₇N₃O₄S
FW: 371.4
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

Z16078526 is supplied as a solid. A stock solution may be made by dissolving the Z16078526 in the solvent of choice, which should be purged with an inert gas. Z16078526 is soluble (≥10 mg/ml) in DMSO.

Description

Z16078526 is an inducer of mitochondrial uncoupling protein 1 (UCP1).¹ It increases levels of Ucp1 in primary mouse brown adipocytes when used at a concentration of 10 μM. Z16078526 (10 μM) increases oxygen consumption rate (OCR), Pka activity, and lipolysis in primary mouse brown adipocytes. It increases body temperature in mice when administered at a dose of 1.5 mg/kg.

Reference

1. Vergnes, L., Lin, J.Y., Davies, G.R., *et al.* Induction of UCP1 and thermogenesis by a small molecule via AKAP1/PKA modulation. *J. Biol. Chem.* **295**(44), 15054-15069 (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, 04/30/2025

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