

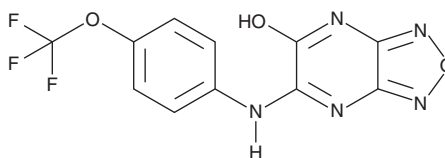
PRODUCT INFORMATION



SHS4121705

Item No. 31053

CAS Registry No.: 2379550-82-8
Formal Name: 6-[[4-(trifluoromethoxy)phenyl]amino]-[1,2,5]oxadiazolo[3,4-b]pyrazin-5(3H)-one
MF: C₁₁H₆F₃N₅O₃
FW: 313.2
Purity: ≥98%
UV/Vis.: λ_{max}: 231, 331 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

SHS4121705 is supplied as a crystalline solid. A stock solution may be made by dissolving the SHS4121705 in the solvent of choice, which should be purged with an inert gas. SHS4121705 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of SHS4121705 in these solvents is approximately 30, 25, and 10 mg/ml, respectively.

Description

SHS4121705 is an orally bioavailable mitochondrial uncoupler.¹ It increases oxygen consumption rate in L6 rat myoblast cells with an EC₅₀ value of 4.3 μM. SHS4121705 (25 mg/kg per day in the diet) reduces hepatic steatosis, liver triglyceride levels, and plasma alanine aminotransferase (ALT) levels in Stelic animal model (STAM) mice, a model of non-alcoholic steatohepatitis (NASH).

Reference

1. Salamoun, J.M., Garcia, C.J., Hargett, S.R., *et al.* 6-Amino[1,2,5]oxadiazolo[3,4-b]pyrazin-5-ol derivatives as efficacious mitochondrial uncouplers in STAM mouse model of nonalcoholic steatohepatitis. *J. Med. Chem.* **63**(11), 6203-6224 (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, 11/07/2025

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