

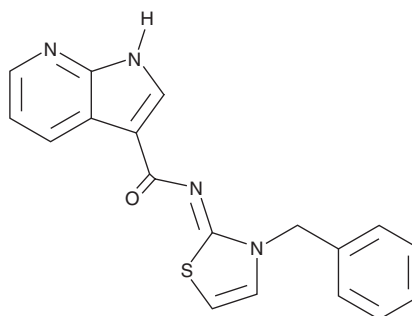
PRODUCT INFORMATION



LATS-IN-1

Item No. 36623

CAS Registry No.: 1424635-83-5
Formal Name: N-[3-(phenylmethyl)-2(3H)-thiazolylidene]-1H-pyrrolo[2,3-b]pyridine-3-carboxamide
Synonyms: LATS Inhibitor 1, TRULI
MF: C₁₈H₁₄N₄OS
FW: 334.4
Purity: ≥98%
UV/Vis.: λ_{max}: 330 nm
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

LATS-IN-1 is supplied as a solid. A stock solution may be made by dissolving the LATS-IN-1 in the solvent of choice, which should be purged with an inert gas. LATS-IN-1 is soluble in the organic solvent DMSO at a concentration of approximately 50 mg/ml.

Description

LATS-IN-1 is an inhibitor of large tumor suppressor kinase 1 (LATS1) and LATS2 (IC₅₀ = 2 nM for both), key regulators of Hippo regenerative signaling.¹ It inhibits phosphorylation of YES1-associated transcriptional regulator (YAP1) in serum-starved HEK293A cells (EC₅₀ = 510 nM). LATS-IN-1 (10 μM) increases the proliferation of Müller glial cells and isolated mouse cardiomyocytes.

Reference

1. Kastan, N., Gnedeva, K., Alisch, T., *et al.* Small-molecule inhibition of Lats kinases may promote Yap-dependent proliferation in postmitotic mammalian tissues. *Nat. Commun.* **12**(1), 3100 (2021).

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, 11/30/2022

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