

PRODUCT INFORMATION

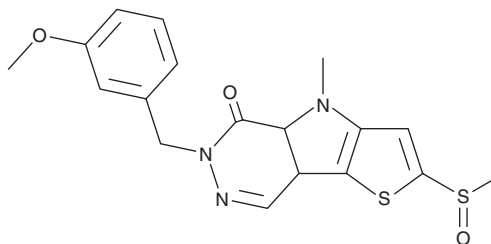


ML-202

Item No. 37326

CAS Registry No.: 1221186-52-2
Formal Name: 4,6-dihydro-6-[(3-methoxyphenyl)methyl]-4-methyl-2-(methylsulfinyl)-5H-thieno[2',3':4,5]pyrrolo[2,3-d]pyridazin-5-one

MF: C₁₈H₁₇N₃O₃S₂
FW: 387.5
Purity: ≥98%
UV/Vis.: λ_{max}: 270, 305 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

ML-202 is supplied as a solid. A stock solution may be made by dissolving the ML-202 in the solvent of choice, which should be purged with an inert gas. ML-202 is soluble in the organic solvent DMSO.

Description

ML-202 is an activator of pyruvate kinase M2 (PKM2).¹ It induces PKM2 activation in a cell-free assay with a 50% activation concentration (AC₅₀) value of 73 nM.

Reference

1. Jiang, J., Boxer, M.B., Heiden, M.G.V., *et al.* Evaluation of thieno[3,2-b]pyrrole[3,2-d]pyridazinones as activators of the tumor cell specific M2 isoform of pyruvate kinase. *Bioorg. Med. Chem. Lett.* **20(11)**, 3387-3393 (2010).

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, 10/02/2023

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