

PRODUCT INFORMATION



ML-265

Item No. 13942

CAS Registry No.: 1221186-53-3
Formal Name: 6-[(3-aminophenyl)methyl]-4,6-dihydro-4-methyl-2-(methylsulfinyl)-5H-thieno[2',3':4,5]pyrrolo[2,3-d]pyridazin-5-one

Synonyms: CID-44246499, NCGC00186528, TEPP-46

MF: C₁₇H₁₆N₄O₂S₂

FW: 372.5

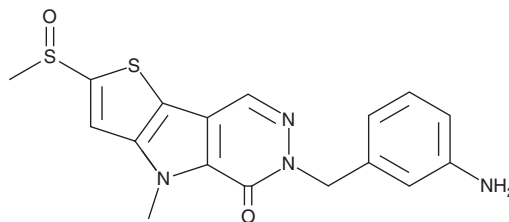
Purity: ≥95%

UV/Vis.: λ_{max}: 269, 304 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

ML-265 is supplied as a crystalline solid. A stock solution may be made by dissolving the ML-265 in the solvent of choice, which should be purged with an inert gas. ML-265 is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml.

Description

ML-265 activates tumor-specific PKM2 (EC₅₀ = 92 nM) by binding to the dimer-dimer interface between two subunits of PKM2 and inducing tetramerization, which is the most active form of the enzyme.¹ It demonstrates >100-fold selectivity for PKM2 over the related PKM1, PKR, and PKL isoforms.¹ At 50 mg/kg, ML-265 has been shown to reduce tumor size, weight, and occurrence in mice bearing H1299 cell xenografts in a model of human non-small cell lung carcinoma.¹

Reference

1. Walsh, M.J., Brimacombe, K.R., Anastasiou, D., *et al.* ML265: A potent PKM2 activator induces tetramerization and reduces tumor formation and size in a mouse xenograft model. *Probe Reports from the NIH Molecular Libraries Program* 1-21 (2012).

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, 11/18/2022

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