

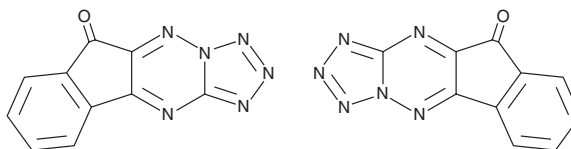
PRODUCT INFORMATION



KP372-1

Item No. 17808

CAS Registry No.: 1374996-60-7
Formal Name: 6H-indeno[1,2-ε]tetrazolo[1,5-b]
[1,2,4]triazin-6-one, compd. with
10H-indeno[2,1-ε]tetrazolo[1,5-b]
[1,2,4]triazin-10-one (1:1)
MF: C₁₀H₄N₆O • C₁₀H₄N₆O
FW: 224.2
Purity: ≥95% (sum of isomers)
UV/Vis.: λ_{max}: 218, 267 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

KP372-1 is supplied as a crystalline solid. A stock solution may be made by dissolving the KP372-1 in the solvent of choice, which should be purged with an inert gas. KP372-1 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of KP372-1 in these solvents is approximately 20 mg/ml.

KP372-1 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, KP372-1 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. KP372-1 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

KP372-1 is a specific Akt inhibitor that demonstrates at least 10-fold selectivity against a panel of additional kinase targets, including CDK1, ERK1, GSK3β, LCK, MEK1, PKA, PKC, and S6K.¹ In acute myelogenous leukemia cells, KP372-1 is reported to inhibit the kinase activity of Akt, PDK1, and FLT3, decreasing the colony-forming ability of these cells with an IC₅₀ value less than 200 nM.¹

Reference

1. Zeng, Z., Samudio, I.J., Zhang, W., *et al.* Simultaneous inhibition of PDK1/AKT and Fms-like tyrosine kinase 3 signaling by a small-molecule KP372-1 induces mitochondrial dysfunction and apoptosis in acute myelogenous leukemia. *Cancer Res.* **66**(7), 3737-3746 (2006).

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, 04/06/2023

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