

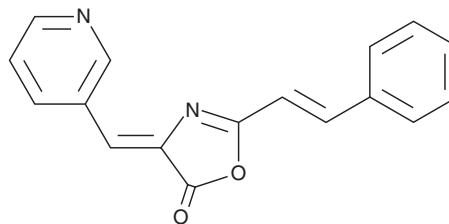
PRODUCT INFORMATION



TC-DAPK6

Item No. 42378

CAS Registry No.: 315694-89-4
Formal Name: 2-(2-phenylethenyl)-4-(3-pyridinylmethylene)-5(4H)-oxazolone
MF: $C_{17}H_{12}N_2O_2$
FW: 276.3
Purity: $\geq 98\%$
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

TC-DAPK6 is supplied as a solid. A stock solution may be made by dissolving the TC-DAPK6 in the solvent of choice, which should be purged with an inert gas. TC-DAPK6 is slightly soluble (0.1-1 mg/ml) in acetonitrile and DMSO.

Description

TC-DAPK6 is an inhibitor of death-associated protein kinase 1 (DAPK1) and DAPK3 ($IC_{50}S = 69$ and 225 nM, respectively).¹ It is selective for DAPK1 and DAPK3 over 47 other kinases at 10 μM . Hollow mesoporous silica nanocarriers (HMSNs) encapsulating TC-DAPK6 increase the latency to fall and distance traveled on the rod in the rotarod test and decrease the time spent in the center in the open field test in a mouse model of kainic acid-induced epilepsy.²

References

1. Okamoto, M., Takayama, K., Shimizu, T., *et al.* Identification of death-associated protein kinases inhibitors using structure-based virtual screening. *J. Med. Chem.* **52**(22), 7323-7327 (2009).
2. Geng, C., Ren, X., Cao, P., *et al.* Macrophage membrane-biomimetic nanoparticles target inflammatory microenvironment for epilepsy treatment. *Theranostics* **14**(17), 6652-6670 (2024).

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, 01/29/2025

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