

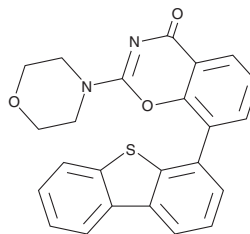
# PRODUCT INFORMATION



## LTURM 34

Item No. 35571

**CAS Registry No.:** 1879887-96-3  
**Formal Name:** 8-(4-dibenzothieryl)-2-(4-morpholinyl)-4H-1,3-benzoxazin-4-one  
**MF:** C<sub>24</sub>H<sub>18</sub>N<sub>2</sub>O<sub>3</sub>S  
**FW:** 414.5  
**Purity:** ≥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

LTURM 34 is supplied as a solid. A stock solution may be made by dissolving the LTURM 34 in the solvent of choice, which should be purged with an inert gas. LTURM 34 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of LTURM 34 in these solvents is approximately 2 and 5 mg/ml, respectively.

### Description

LTURM 34 is an inhibitor of DNA protein kinase (DNA-PK; IC<sub>50</sub> = 0.034 μM).<sup>1</sup> It is selective for DNA-PK over PI3Kα, PI3Kβ, PI3Kγ, and PI3Kδ at 10 μM. LTURM 34 (10 μM) inhibits the growth of HOP-92, SNB-75, and UO-31 cancer cells.

### Reference

1. Morrison, R., Al-Rawi, J.M.A., Jennings, I.G., *et al.* Synthesis, structure elucidation, DNA-PK and PI3K and anti-cancer activity of 8- and 6-aryl-substituted-1-3-benzoxazines. *Eur. J. Med. Chem.* **110**, 326-339 (2016).

#### 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, 11/23/2022

#### 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](http://WWW.CAYMANCHEM.COM)