

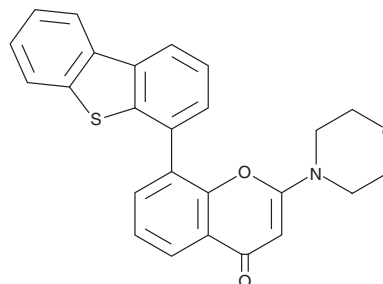
# PRODUCT INFORMATION



NU 7441

Item No. 14881

**CAS Registry No.:** 503468-95-9  
**Formal Name:** 8-(4-dibenzothieryl)-2-(4-morpholinyl)-4H-1-benzopyran-4-one  
**Synonym:** Ku-57788  
**MF:** C<sub>25</sub>H<sub>19</sub>NO<sub>3</sub>S  
**FW:** 413.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 234, 288 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

NU 7441 is supplied as a crystalline solid. A stock solution may be made by dissolving the NU 7441 in the solvent of choice, which should be purged with an inert gas. NU 7441 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of NU 7441 in these solvents is approximately 12 and 1 mg/ml, respectively.

NU 7441 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, NU 7441 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. NU 7441 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

DNA-dependent protein kinase (DNA-PK) catalyzes nonhomologous end-joining, which is required to repair lethal DNA double-strand breaks. Because cells that are defective in DNA double-strand break repair are highly sensitive to ionizing radiation and topoisomerase II poisons, modulating DNA-PK is one strategy to defer cancer cell resistance to radiation or chemotherapeutic treatments.<sup>1</sup> NU 7441 is a selective DNA-PK inhibitor with an IC<sub>50</sub> value of 14 nM.<sup>2</sup> It inhibits other members of the PI3K-related kinase family, including mTOR, PI3K, ataxia telangiectasia mutated (ATM), and ataxia telangiectasia and Rad3 related (ATR) with IC<sub>50</sub> values of 1.7, 5, >100, and >100 μM, respectively.<sup>2</sup> NU 7441 has been shown to increase the cytotoxicity of ionizing radiation and etoposide (Item No. 12092) in human colon cancer cell lines *in vitro* and to potentiate the effects of etoposide in mice bearing human colon cancer xenograft tumors *in vivo*.<sup>3</sup>

## References

1. Veuger, S.J., Curtin, N.J., Richardson, C.J., *et al.* Radiosensitization and DNA repair inhibition by the combined use of novel inhibitors of DNA-dependent protein kinase and poly(ADP-ribose) polymerase-1. *Cancer Res.* **63(18)**, 6008-6015 (2003).
2. Leahy, J.J.J., Golding, B.T., Griffin, R.J., *et al.* Identification of a highly potent and selective DNA-dependent protein kinase (DNA-PK) inhibitor (NU7441) by screening of chromenone libraries. *Bioorg. Med. Chem. Lett.* **14(24)**, 6083-6087 (2004).
3. Zhao, Y., Thomas, H.D., Batey, M.A., *et al.* Preclinical evaluation of a potent novel DNA-dependent protein kinase inhibitor NU7441. *Cancer Res.* **66(10)**, 5354-5362 (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, 12/12/2022

## 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