

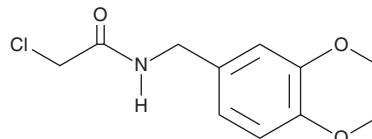
PRODUCT INFORMATION



DKM 2-93

Item No. 37293

CAS Registry No.: 65836-72-8
Formal Name: 2-chloro-N-[(3,4-dimethoxyphenyl)methyl]-acetamide
MF: C₁₁H₁₄ClNO₃
FW: 243.7
Purity: ≥98%
UV/Vis.: λ_{max}: 230 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

DKM 2-93 is supplied as a crystalline solid. A stock solution may be made by dissolving the DKM 2-93 in the solvent of choice, which should be purged with an inert gas. DKM 2-93 is soluble in chloroform and methanol.

Description

DKM 2-93 is a cysteine-reactive covalent inhibitor of ubiquitin-like modifier activating enzyme 5 (UBA5; IC₅₀ = 450 μM).¹ It decreases survival of PaCa-2 pancreatic cancer cells without affecting human pancreatic ductal epithelial cells (HPDEs) when used at a concentration of 50 μM. *In vivo*, DKM 2-93 (50 mg/kg per day) reduces tumor volume without reducing body weight in a PaCa-2 pancreatic cancer mouse xenograft model. It has also been used as an intermediate in the synthesis of topoisomerase I inhibitors and compounds active against *M. tuberculosis*.^{2,3}

References

1. Roberts, A.M., Miyamoto, D.K., Huffman, T.R., *et al.* Chemoproteomic screening of covalent ligands reveals UBA5 as a novel pancreatic cancer target. *ACS Chem. Biol.* **12(4)**, 899-904 (2017).
2. Kadagathur, M., Patra, S., Devabattula, G., *et al.* Design, synthesis of DNA-interactive 4-thiazolidinone-based indolo-/pyrroloazepinone conjugates as potential cytotoxic and topoisomerase I inhibitors. *Eur. J. Med. Chem.* **238**, 114465 (2022).
3. Lee, R.E., Protopopova, M., Crooks, E., *et al.* Combinatorial lead optimization of [1,2]-diamines based on ethambutol as potential antituberculosis preclinical candidates. *J. Comb. Chem.* **5(2)**, 172-187 (2003).

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, 10/05/2023

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