

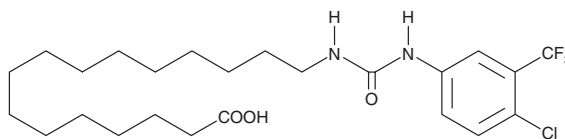
PRODUCT INFORMATION



CAY10726

Item No. 23859

CAS Registry No.: 1611446-66-2
Formal Name: 16-[[[4-chloro-3-(trifluoromethyl)phenyl]amino]carbonyl]amino]hexadecanoic acid
MF: C₂₄H₃₆ClF₃N₂O₃
FW: 493.0
Purity: ≥98%
UV/Vis.: λ_{max}: 253 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

CAY10726 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10726 in the solvent of choice, which should be purged with an inert gas. CAY10726 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CAY10726 in these solvents is approximately 12, 5, and 20 mg/ml, respectively.

CAY10726 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CAY10726 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. CAY10726 has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CAY10726 is an arylurea fatty acid.¹ It decreases ATP production by 28% in MDA-MB-231 breast cancer cells when used at a concentration of 10 μM. CAY10726 decreases proliferation and initiates apoptosis in MDA-MB-231 cells *via* depletion of the phospholipid cardiolipin and its precursor phosphatidylglycerol from the mitochondrial membrane. *In vivo*, CAY10726 (2.5-40 mg/kg) reduces tumor volume and increases tumor cell apoptosis in a mouse MDA-MB-231 breast cancer xenograft model in a dose-dependent manner.

Reference

1. Rawling, T., Choucair, H., Koolaji, N., *et al.* A novel arylurea fatty acid that targets the mitochondrion and depletes cardiolipin to promote killing of breast cancer cells. *J. Med. Chem.* **60**(20), 8661-8666 (2017).

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, 02/07/2025

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