

PRODUCT INFORMATION



T-26c

Item No. 19962

CAS Registry No.: 869296-13-9

Formal Name: 4-[[[1,4-dihydro-2-[[[(3-methoxyphenyl)methyl]amino]carbonyl]-4-oxothieno[2,3-d]pyrimidin-5-yl]methoxy]methyl]-benzoic acid

MF: C₂₄H₂₁N₃O₆S

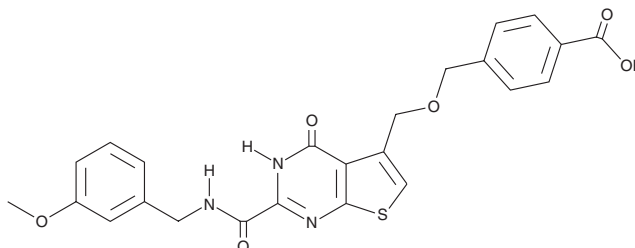
FW: 479.5

Purity: ≥98%

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

T-26c is supplied as a crystalline solid. A stock solution may be made by dissolving the T-26c in the solvent of choice, which should be purged with an inert gas. T-26c is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml.

Description

T-26c is a selective inhibitor of matrix metalloproteinase-13 (MMP-13; IC₅₀ = 0.0069 nM).¹ It is selective for MMP-13 over MMP-1, -2, -3, -7, -8, -9, -10, and -14 (IC₅₀s = ≥18 nM). T-26c inhibits IL-1β- and oncostatin M-induced collagen degradation in bovine nasal septum cartilage explants by 87.4% when used at a concentration of 0.1 μM.

Reference

1. Nara, H., Sato, K., Naito, T., *et al.* Thieno[2,3-d]pyrimidine-2-carboxamides bearing a carboxybenzene group at 5-position: Highly potent, selective, and orally available MMP-13 inhibitors interacting with the S1'' binding site. *Bioorg. Med. Chem.* **22**(19), 5487-5505 (2014).

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, 11/01/2022

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