

# PRODUCT INFORMATION



## STD1T

Item No. 37791

CAS Registry No.: 893075-58-6

Formal Name: 4,5,6,7-tetrahydro-2-[[[2-[[5-(2-thienyl)-3-isoxazolyl]methoxy]acetyl]amino]-benzo[b]thiophene-3-carboxamide

MF:  $C_{19}H_{19}N_3O_4S_2$

FW: 417.5

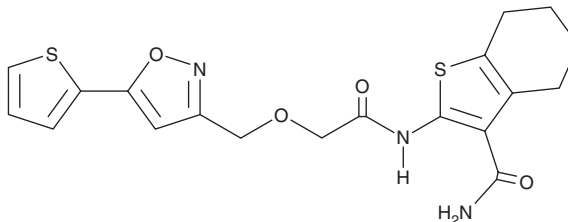
Purity:  $\geq 98\%$

UV/Vis.:  $\lambda_{max}$ : 225, 256, 262, 292 nm

Supplied as: A solid

Storage:  $-20^{\circ}\text{C}$

Stability:  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

STD1T is supplied as a solid. A stock solution may be made by dissolving the STD1T in the solvent of choice, which should be purged with an inert gas. STD1T is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of STD1T in these solvents is approximately 16 and 20 mg/ml, respectively.

### Description

STD1T is an inhibitor of ubiquitin-specific protease 2a (USP2a;  $IC_{50} = 3.3 \mu\text{M}$ ).<sup>1</sup> It is selective for USP2a over USP7 at 10  $\mu\text{M}$ . STD1T (20  $\mu\text{M}$ ) decreases cyclin D1 protein levels in HCT116 colon and MCF-7 breast cancer cells.

### Reference

1. Tomala, M.D., Magiera-Mularz, K., Kubica, K., *et al.* Identification of small-molecule inhibitors of USP2a. *Eur. J. Med. Chem.* **150**, 261-267 (2018).

#### 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, 10/19/2022

#### 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