

PRODUCT INFORMATION



SMPT

Item No. 44829

CAS Registry No.: 112241-19-7

Formal Name: 4-[1-(2-pyridyldithio)ethyl]-benzoic acid, 2,5-dioxo-1-pyrrolidinyl ester

Synonym: 4-Succinimidylloxycarbonyl- α -methyl- α -(2-pyridyldithio)toluene

MF: $C_{18}H_{16}N_2O_4S_2$

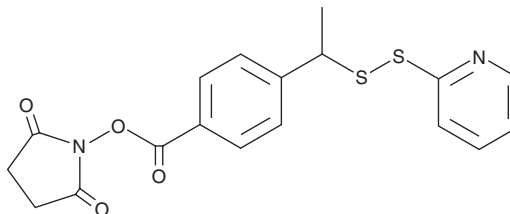
FW: 388.5

Purity: $\geq 90\%$

Supplied as: A 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

SMPT is supplied as a solid. A stock solution may be made by dissolving the SMPT in the solvent of choice, which should be purged with an inert gas. SMPT is soluble (≥ 10 mg/ml) in DMSO.

Description

SMPT is an antibody-toxin cross-linking agent that contains an amine-reactive N-hydroxysuccinimide (NHS) group and a sulfhydryl-reactive pyridyldithio group.¹ SMPT has been used in the formation of antibody-toxin conjugates to induce toxicity in lymphoma and leukemia cells. SMPT-crosslinked immunotoxins have increased stability *in vivo* compared to those cross-linked using succinimidyl 3-(2-pyridyldithio) propionate (SPDP) or 2-iminothiolane (2-IT).

Reference

1. Thorpe, P.E., Wallace, P.M., Knowles, P.P., *et al.* New coupling agents for the synthesis of immunotoxins containing a hindered disulfide bond with improved stability *in vivo*. *Cancer Res.* **47(22)**, 5924-5931 (1987).

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, 03/05/2026

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