

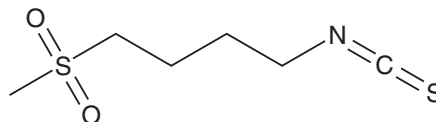
PRODUCT INFORMATION



Erysolin

Item No. 39661

CAS Registry No.: 504-84-7
Formal Name: 1-isothiocyanato-4-(methylsulfonyl)-butane
MF: $C_6H_{11}NO_2S_2$
FW: 193.3
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 204, 241 nm
Supplied as: A solid
Storage: $-20^\circ C$
Stability: ≥ 4 years
Item Origin: Synthetic



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

Laboratory Procedures

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

Description

Erysolin is an isothiocyanate that has been found in *E. perofskianum* and has anticancer activity.¹ It inhibits the proliferation of MCF-7, HeLa, A549, HepG2, and SW480 cells (IC_{50} s = 1.62, 1.12, 5.22, 14, and 8.14 μM , respectively). Erysolin (3 μM) enhances arsenic trioxide-induced apoptosis and cytotoxicity in HL-60, U937, and K562 cells.²

References

1. Li, B.-L., Zhao, S.-Z., Zhou, H., *et al.* Synthesis and antitumor activity of erysolin and its metabolites. *J. Asian Nat. Prod. Res.* **25**(4), 369-378 (2023).
2. Doudican, N.A., Bowling, B., and Orlow, S.J. Enhancement of arsenic trioxide cytotoxicity by dietary isothiocyanates in human leukemic cells via a reactive oxygen species-dependent mechanism. *Leuk. Res.* **34**(2), 229-234 (2010).

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, 10/16/2023

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