

S-Trityl-L-cysteine

Cell cycle blocker

Potent, cell permeable cysteine thioether displaying anti-mitotic and anti-tumor properties. Inhibits mitosis by blocking basal and microtubule-activated ATPase activities of the mitotic kinesin Eg5 ($IC_{50}=1\mu M$ and $140nm$). Induces mitotic arrest in HeLa cells ($IC_{50}=700nM$) and inhibits the growth of NCI tumor cell line panel ($GI_{50}=1.31\mu M$).

Citations: 12

[View Online »](#)

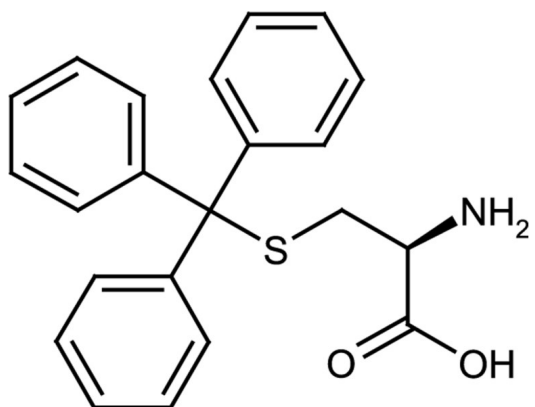
Ordering Information

[Order Online »](#)

ALX-105-011-M500	500mg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name NSC 83265, STLC

Appearance White to off-white solid.

CAS 2799-07-7

Couple Target Kinesin

Couple Type Inhibitor

Formula $C_{22}H_{21}NO_2S$

MW 363.5

Purity ≥98% (TLC)

RTECS AY7710000

Solubility Soluble in DMSO or acetic acid.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)