

Temozolomide

AGT inhibitor

Antitumor and antiangiogenic compound. DNA alkylating agent. Induces G2/M arrest and apoptosis through adduction of a methyl group to O6 position of guanine in genomic DNA. Inactivates DNA repair protein O(6)-alkylguanine DNA alkyltransferase (AGT) in base excision repair (BER) pathway. Chemotherapeutic agent. Modifies drug resistance.

Citations: 17

[View Online »](#)

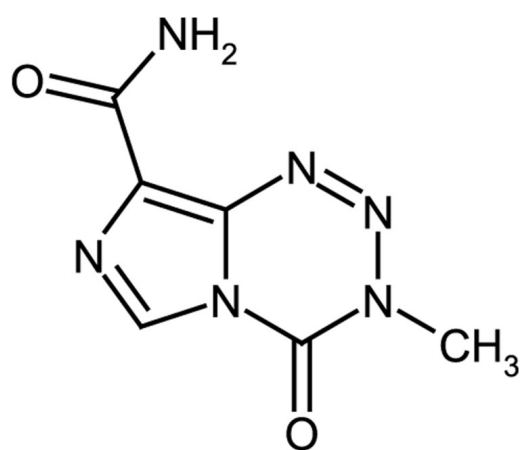
Ordering Information

[Order Online »](#)

ALX-420-044-M025	25mg
ALX-420-044-M100	100mg

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 8-Carbamoyl-3-methylimidazo[5,1-d]-1,2,3,5-tetrazin-4(3H)-one

Appearance White to off-white crystalline powder.

CAS 85622-93-1

Couple Target DNA alkyltransferase

Couple Type Inhibitor

Formula $C_6H_6N_6O_2$

MI **14: 9139**

MW 194.2

Purity ≥98% (HPLC)

RTECS NJ5927050

Solubility Soluble in DMSO; slightly soluble in water.

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-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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

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