

(-)-p-Bromotetramisole . oxalate

Alkaline phosphatase inhibitor

Potent alkaline phosphatase inhibitor. Mimics the action of orthovanadate in the potentiation of fluorouracil antiproliferative activity. Cell permeable.

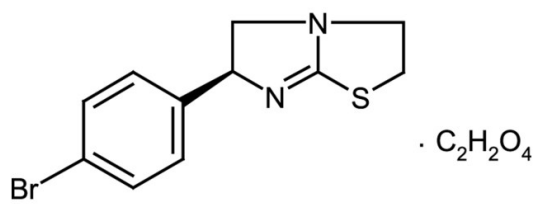
Ordering Information

[Order Online »](#)

ALX-430-025-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name (-)-(S)-6-(4-Bromophenyl)-2,3,5,6-tetrahydroimidazo[2,1-b]thiazole . oxalate, L-p-Bromotetramisole . oxalate

Appearance White powder.

CAS 62284-79-1

Couple Target Alkaline phosphatase

Couple Type Inhibitor

Formula $C_{11}H_{11}BrN_2S \cdot C_2H_2O_4$

MW 283.2 . 90.0

Purity ≥98% (HPLC)

Solubility Soluble in water, methanol, DMF or DMSO.

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)