

# (-)-p-Bromotetramisole . oxalate

## Alkaline phosphatase inhibitor

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

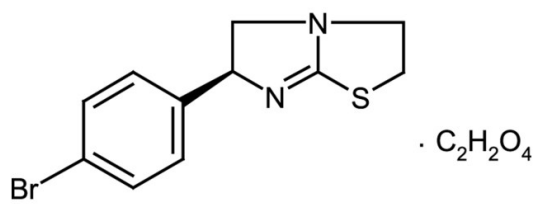
### Ordering Information

[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)