

Dolastatin 15

Microtubule inhibitor

Marine natural product from the sea hare *Dolabella auricularia*. Noncompetitively inhibits binding of vincristine to tubulin but has been shown to bind to the RZX/MAY region. Inhibits microtubule assembly and induces apoptosis and Bcl-2 phosphorylation in several malignant cell types.

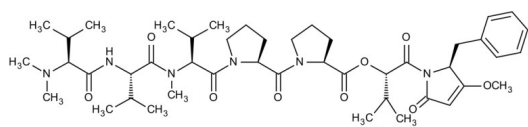
Ordering Information

[Order Online »](#)

BML-T124-0001	1mg
---------------	-----

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	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White solid.
CAS	123884-00-4
Couple Target	Microtubule
Couple Type	Inhibitor
MW	837.1
Purity	≥98% (HPLC)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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