

Wiskostatin

Blocks actin filament assembly

A selective, reversible inhibitor of N-WASP (neural Wiskott-Aldrich syndrome protein) that selectively blocks actin filament assembly. Appears to bind to N-WASP, stabilize the autoinhibited conformation and prevent the activation of Arp2/3 (actin-related protein 2/3) complex.

Citations: 2

[View Online »](#)

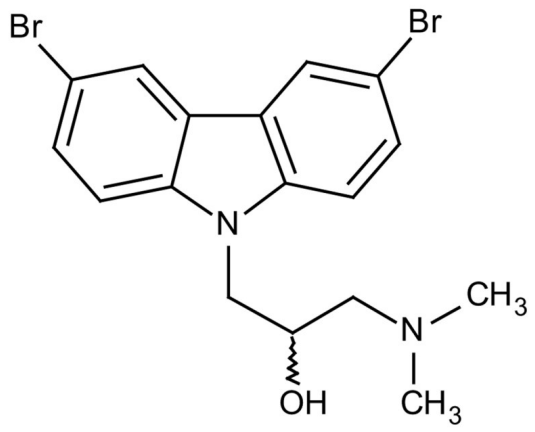
Ordering Information

[Order Online »](#)

BML-T126-0001	1mg
---------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at room temperature for up to 1 year. Store solutions at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance	White to off-white solid.
CAS	253449-04-6
MW	426.15
MeltingPoint	148-150°C
Purity	≥98% (HPLC)
Solubility	Soluble in ethanol (10 mg/ml with warming) or DMSO (25 mg/ml).

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