

Nocodazole

Microtubule inhibitor

Synthetic reversible microtubule inhibitor.

Citations: 11

[View Online »](#)

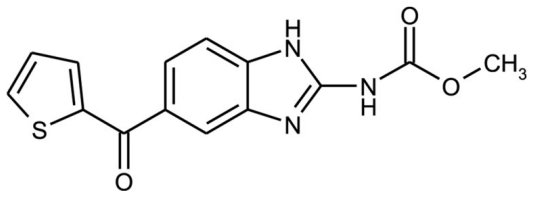
Ordering Information

[Order Online »](#)

BML-T101-0010	10mg
---------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored at +20°C. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Methyl[5-(2-thienylcarbonyl)-1H-benzimidazol-2-yl]carbamate, R 17934
Appearance	Solid. White, off-white, orange and possibly even a brown/pink color.
CAS	31430-18-9
Couple Target	Microtubule
Couple Type	Inhibitor
Formula	$C_{14}H_{11}N_3O_3S$
MW	301.3
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (10mg/ml) or dimethyl formamide; almost insoluble in water.

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