

GW-441756

TrkA inhibitor

Potent and selective inhibitor of NGFR kinase (TrkA, $IC_{50}=2$ nM). Blocks TrkA-mediated γ H2AX production and apoptosis in TrkA-overexpressing cells.

Citations: 1

[View Online »](#)

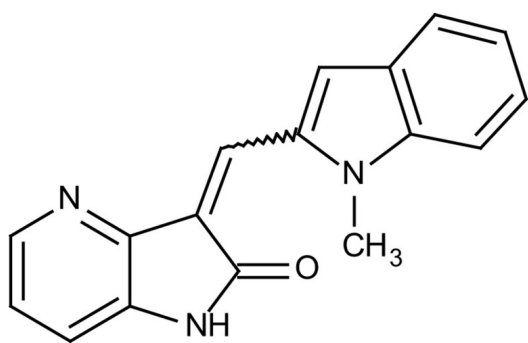
Ordering Information

[Order Online »](#)

BML-EI364-0010	10mg
BML-EI364-0050	50mg

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	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Orange solid.
CAS	504433-23-3
Couple Target	Trk
Couple Type	Inhibitor
Formula	$C_{17}H_{13}N_3O$
MW	275.3
Purity	≥98% (TLC); mixture of E and Z isomers
Solubility	Soluble in DMSO (>25 mg/ml).

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-
be@enzolifesciences.com](mailto:info-be@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)