

Gambogic amide

TrkA ligand

A novel small molecule NGF mimic discovered using a cell-based chemical genetic screen for TrkA agonists. Gambogic amide selectively binds to TrkA, triggers its tyrosine phosphorylation, elicits PI3-kinase/Akt and MAPK activation, provokes neurite outgrowth in PC12 cells and prevents neuronal cell death. EC₅₀ for antiapoptotic activity is 5nM (see reference S.-W. Jang, et al. 2007).

Citations: 3

[View Online »](#)

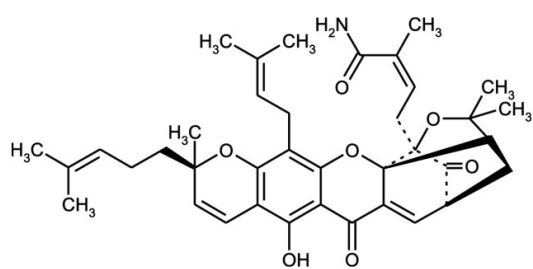
Ordering Information

[Order Online »](#)

BML-N159-0001	1mg
BML-N159-0005	5mg

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 Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance Yellow-orange solid.

CAS 286935-60-2

Couple Target Trk

Couple Type Ligand

Formula $C_{38}H_{45}NO_7$

MW 627.8

Purity ≥95% (HPLC)

Solubility Soluble in DMSO (20mg/ml) or ethanol (20mg/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-
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)