

AG-879

NGF receptor inhibitor

AG-879, a member of the tyrphostin family of tyrosine kinase inhibitors, inhibits NGF receptor (pp140^{C-trk}) autophosphorylation selectively with no inhibition of EGF or PDGF receptor phosphorylation (IC₅₀=10 μM). AG-879 inhibits NGF-induced neurite outgrowth in PC12 cells.

Citations: 1

[View Online »](#)

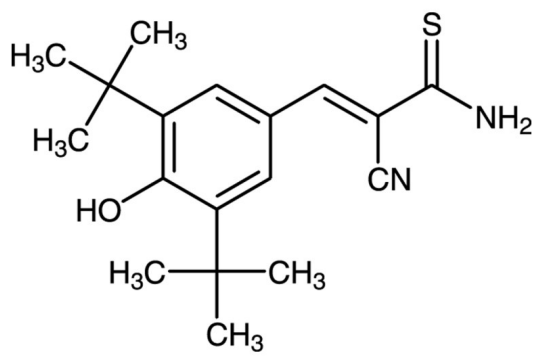
Ordering Information

[Order Online »](#)

BML-EI258-0005	5mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Yellow solid.
CAS	148741-30-4
Couple Target	Trk
Couple Type	Inhibitor
Formula	$C_{18}H_{24}N_2OS$
MW	316.5
MeltingPoint	223-224°C
Purity	≥99% (HPLC)
Solubility	Soluble in 100% ethanol (25mg/ml) or DMSO (25mg/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