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 (IC $_{50}$ =10 μ M). AG-879 inhibits NGF-induced neurite outgrowth in PC12 cells.

Citations: 1

View Online »

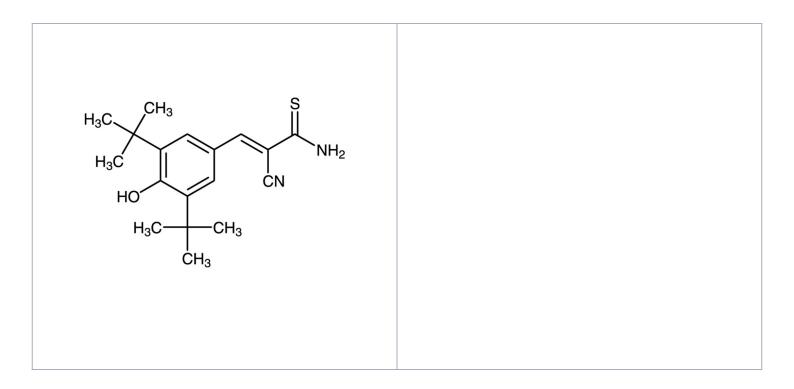
Ordering Information

Order Online »

BML-EI258-0005	5mg
BML-EI258-0025	25mg

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

