

RG-13022

Tyrosine kinase inhibitor

RG-13022 is a non-phenolic tyrphostin analog, which inhibits the EGF receptor ($IC_{50}=5\text{ }\mu\text{M}$) and PDGF receptor tyrosine kinases, and is long-acting. It inhibits EGF-stimulated HER14 cell proliferation ($IC_{50}=1\text{ }\mu\text{M}$) as well as tumor growth *in vivo*. It inhibits both EGF- and PDGF-stimulated DNA synthesis in human glioma cell lines. RG-13022 is non-phenolic and therefore possesses less antioxidant activity than other tyrphostins.

Citations: 2

[View Online »](#)

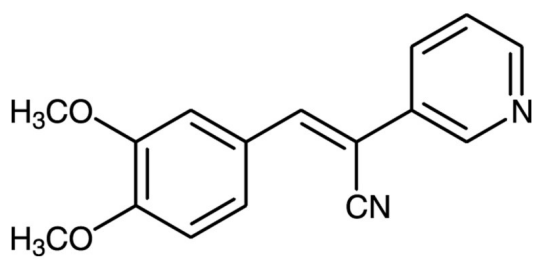
Ordering Information

[Order Online »](#)

BML-EI252-0005	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 room temperature for up to 2 years. Store solutions at -20°C for up to 3 months.
Handling	Store tightly sealed in the dark. Protect from light.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	3,4-Dimethoxy- α -(3-pyridyl)-(Z)-cinnamonnitrile
Appearance	White solid.
CAS	136831-48-6
Couple Target	EGFR, PDGF receptor
Couple Type	Inhibitor
Formula	$C_{16}H_{14}N_2O_2$
MW	266.3
Purity	$\geq 98\%$ (TLC)
Solubility	Soluble at 25mg/ml in DMSO or 100% ethanol.

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)