

# 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 »](#)

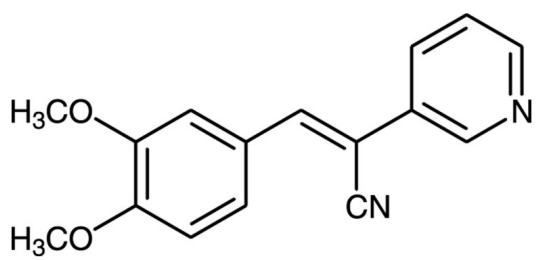
## Ordering Information

[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- $\alpha$ -(3-pyridyl)-(Z)-cinnamionitrile
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.



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)