

AG-1288

Guanylyl cyclase inhibitor

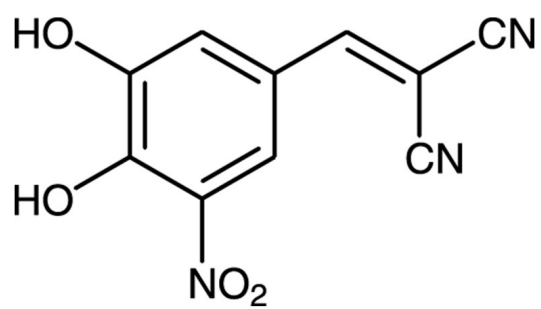
Tyrosine kinase inhibitor. It has been shown to block TNF α induced cytotoxicity and to block glucocorticoid-induced COX-2 activity in human amnion cells (IC_{50} =21 μ M). It is also reported to inhibit guanylyl cyclase.

Citations: 3

[View Online »](#)

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 1 year. 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	α -Cyano-(3,4-dihydroxy-5-nitro)cinnamionitrile
Appearance	Pale orange solid.
CAS	116313-73-6
Couple Target	Guanylyl cyclase
Couple Type	Inhibitor
Formula	$C_{10}H_5N_3O_4$
MW	231.2
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (35mg/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-fr@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