

# AG-126

## Inhibits tyrosine kinase-dependent B cell receptor signaling

AG-126 prevents lipopolysaccharide-induced lethal toxicity (septic shock) in mice. Administered at 1.5 mg/mouse, it reduced LPS-induced lethal toxicity from 95% to 10%. AG-126 greatly reduces LPS-induced tyrosine phosphorylation of p42<sup>MAPK</sup> in murine peritoneal macrophages at 50  $\mu$ M. EGFR, Her-2/neu receptor or PDGF-R are not inhibited at 100  $\mu$ M. Inhibits tyrosine kinase-dependent B cell receptor signaling.

Citations: 9

[View Online »](#)

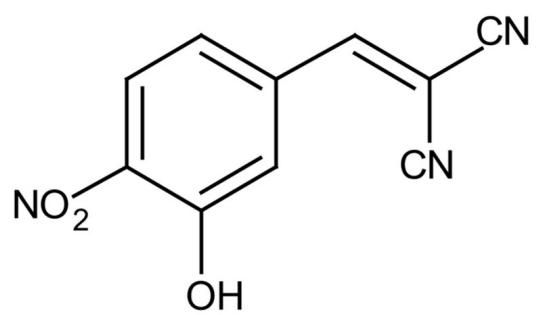
## Ordering Information

[Order Online »](#)

BML-EI263-0005	5mg
----------------	-----

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended.
<b>Handling</b>	Store tightly sealed in the dark. Protect from light.
<b>Long Term Storage</b>	Ambient
<b>Shipping</b>	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	3-Hydroxy-4-nitrobenzylidenemalononitrile
<b>Appearance</b>	Yellow solid.
<b>CAS</b>	118409-62-4
<b>Formula</b>	$C_{10}H_5N_3O_3$
<b>MW</b>	215.2
<b>MeltingPoint</b>	181-182°C
<b>Purity</b>	≥98% (HPLC)
<b>Solubility</b>	Soluble in DMSO (25mg/ml) or 100% ethanol (5mg/ml warm).

**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)