

# Stattic

## STAT inhibitor

Small molecule STAT3-specific inhibitor which blocks activation, dimerization and nuclear translocation via binding to the STAT3 SH2 domain. It is a very useful tool for exploring STAT3-mediated signaling pathways. It blocks LPS-mediated STAT3 tyrosine phosphorylation which leads to inhibition of LPS-mediated IL-1 $\beta$  and IL-6 production. In primary cultures of rat neural stem cells it blocks IL-15-induced differentiation.

Citations: 10

[View Online »](#)

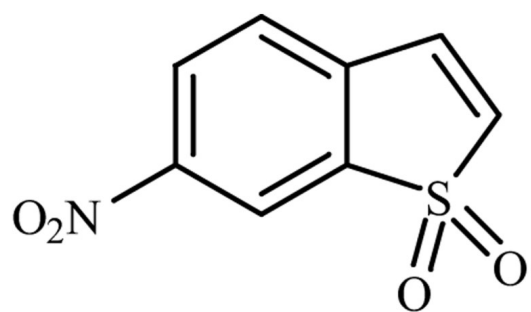
## Ordering Information

[Order Online »](#)

BML-EI368-0010	10mg
----------------	------

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Store solutions at -20°C for up to 3 months.
<b>Long Term Storage</b>	Ambient
<b>Shipping</b>	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	6-Nitro[b]thiophene-1,1-dioxide
<b>Appearance</b>	Pale yellow solid.
<b>CAS</b>	19983-44-9
<b>Couple Target</b>	STAT
<b>Couple Type</b>	Inhibitor
<b>Formula</b>	$C_8H_5NO_4S$
<b>MW</b>	211.2
<b>Purity</b>	≥98% (TLC)
<b>Solubility</b>	Soluble in DMSO (>25 mg/ml), 100% ethanol (10 mg/ml).

**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)