

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 β and IL-6 production. In primary cultures of rat neural stem cells it blocks IL-15-induced differentiation.

Citations: 10

[View Online »](#)

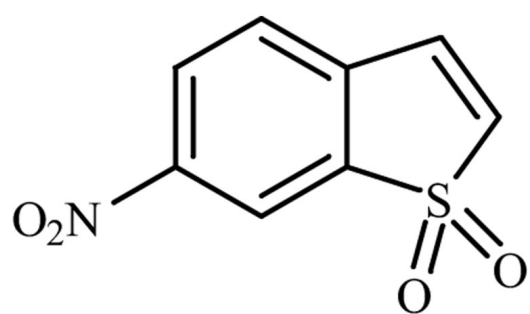
Ordering Information

[Order Online »](#)

BML-EI368-0010	10mg
BML-EI368-0050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	6-Nitro[b]thiophene-1,1-dioxide
Appearance	Pale yellow solid.
CAS	19983-44-9
Couple Target	STAT
Couple Type	Inhibitor
Formula	$C_8H_5NO_4S$
MW	211.2
Purity	≥98% (TLC)
Solubility	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

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