

BAY 11-7082

NF- κ B inhibitor

Inhibitor of cytokine induced I κ B α phosphorylation resulting thereby in a decreased expression of NF- κ B and adhesion molecules. Potential anti-inflammatory agent.

Citations: 27

[View Online »](#)

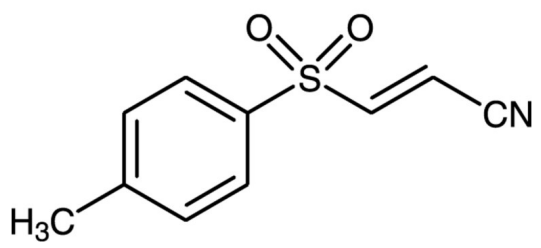
Ordering Information

[Order Online »](#)

BML-EI278-0010	10mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light. Keep under inert gas.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	(E)-3-[(4-methylphenylsulfonyl)-2-propenenitrile
Appearance	White to off-white solid.
CAS	19542-67-7
Couple Target	NF-kappaB
Couple Type	Inhibitor
Formula	$C_{10}H_9NO_2S$
MW	207.2
Purity	≥99% (HPLC)
Solubility	Soluble in DMSO, 100% ethanol, methanol or methylene dichloride; insoluble in water.

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-be@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