

BAY 11-7085

NF- κ B inhibitor

An irreversible inhibitor of cytokine-inducible I κ B α phosphorylation (IC_{50} = 10 μ M). In intact HUVECs it inhibits TNF α -induced phosphorylation and degradation of I κ B α without affecting constitutive phosphorylation. Inhibition of I κ B α phosphorylation and degradation results in the inactivation of NF- κ B and NF- κ B-dependent expression of endothelial-leukocyte adhesion molecules, E-selectin, VCAM-1 and ICAM-1. BAY 11-7085 displayed potent anti-inflammatory activity in the rat carrageenan paw edema assay and adjuvant arthritis model.

Citations: 10

[View Online »](#)

Ordering Information

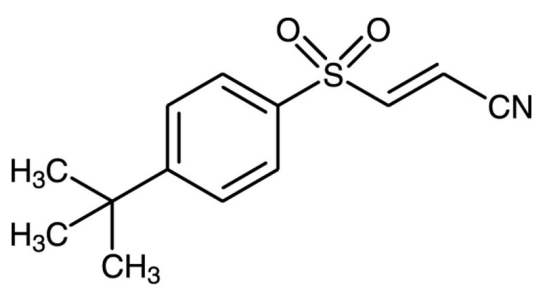
[Order Online »](#)

BML-EI279-0010

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

Appearance	White solid.
CAS	196309-76-9
Couple Target	NF-kappaB
Couple Type	Inhibitor
Formula	$C_{13}H_{15}NO_2S$
MW	249.3
Purity	≥99% (TLC)
Solubility	Soluble in DMSO (>25mg/ml) or 100% ethanol (>25mg/ml). 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