

SC-514

IKK inhibitor

Selective, reversible and ATP-competitive inhibitor of IKK2 (IC_{50} =11.2 μ M). Does not inhibit other IKK isoforms or other serine/threonine or tyrosine kinases. Inhibits transcription of NF- κ B dependent genes in IL-1 β -induced rheumatoid arthritis-derived synovial fibroblasts in a dose-dependent fashion.

Citations: 6

[View Online »](#)

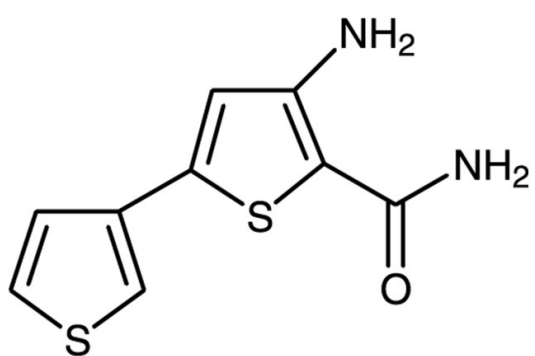
Ordering Information

[Order Online »](#)

BML-EI343-0010	10mg
BML-EI343-0050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	If resuspended in solution, store at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Yellow solid.
CAS	354812-17-2
Couple Target	IKK, NF-kappaB
Couple Type	Inhibitor
Formula	$C_9H_8N_2OS_2$
MW	224.3
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (>25mg/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-
fr@enzolifesciences.com](mailto:info-fr@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)