

QNZ

NF-κB inhibitor

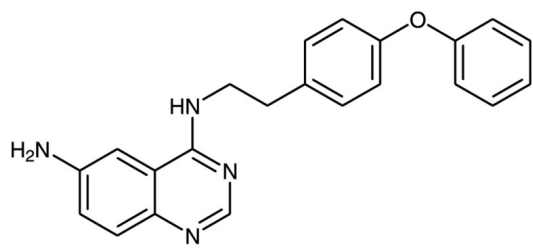
Inhibitor of NF-κB which displays potent inhibitory effects on both NF-κB transcriptional activation (IC_{50} = 11nM) and TNF-α production (IC_{50} = 7nM). A useful tool to explore the role of NF-κB in cellular signaling pathways.

Citations: 3

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance White solid.

CAS 545380-34-5

Couple Target NF-kappaB

Couple Type Inhibitor

Formula $C_{22}H_{20}N_4O$

MW 356.4

Purity ≥98% (TLC)

Solubility Soluble in DMSO (20 mg/ml); DMF (30mg/ml); 100% ethanol (10 mg/ml); not soluble in water.

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)