

NF-κB p50 subunit (human) (35-381), (recombinant) (GST- tag)

NF-κB p50 subunit acts as the DNA binding domain of the major NF-κB heterodimeric complex p50/p65. NF-κB plays an important role in inflammation, autoimmune response, cell proliferation, and apoptosis by regulating the expression of genes involved in these processes. NF-κB pathway studies, in conjunction with the p65 subunit.

Ordering Information

[Order Online »](#)

BML-UW9980-0050	50μg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Notes

Uses:

For NF-κB pathway studies.

Investigate DNA binding of p50.

Investigate inhibition by IκBa.

GST pull down of interacting partners.

Formulation

Liquid. In 50mM HEPES, pH 7.5, 50mM NaCl, 1mM DTT.

MW

65.0 kDa

Purity

≥95% (SDS-PAGE)

Purity Detail

Purified by glutathione affinity chromatography.

Source

Produced in *E. coli*.

UniProt ID

P19838

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)