

NF-κB (p65) (human), (recombinant) (His- tag)

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-201-284-C002	2μg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store in liquid nitrogen.
Short Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Concentration	0.5mg/ml
Formulation	Liquid. In 50mM TRIS-HCl, pH 7.5, containing 100mM sodium chloride, 0.2% NP-40, 50mM imidazole and 10% glycerol.
Purity	≥95% (SDS-PAGE)
Source	Produced in SF21 insect cells. Fused to a His-tag. Produced in a baculovirus expression system.
Specific Activity	~1ng is required for high mobility shift assay. ~5ng are required for reconstituted transcription assays.
UniProt ID	Q04206

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