

IKK γ (human) monoclonal antibody (46B844)

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 1

[View Online »](#)

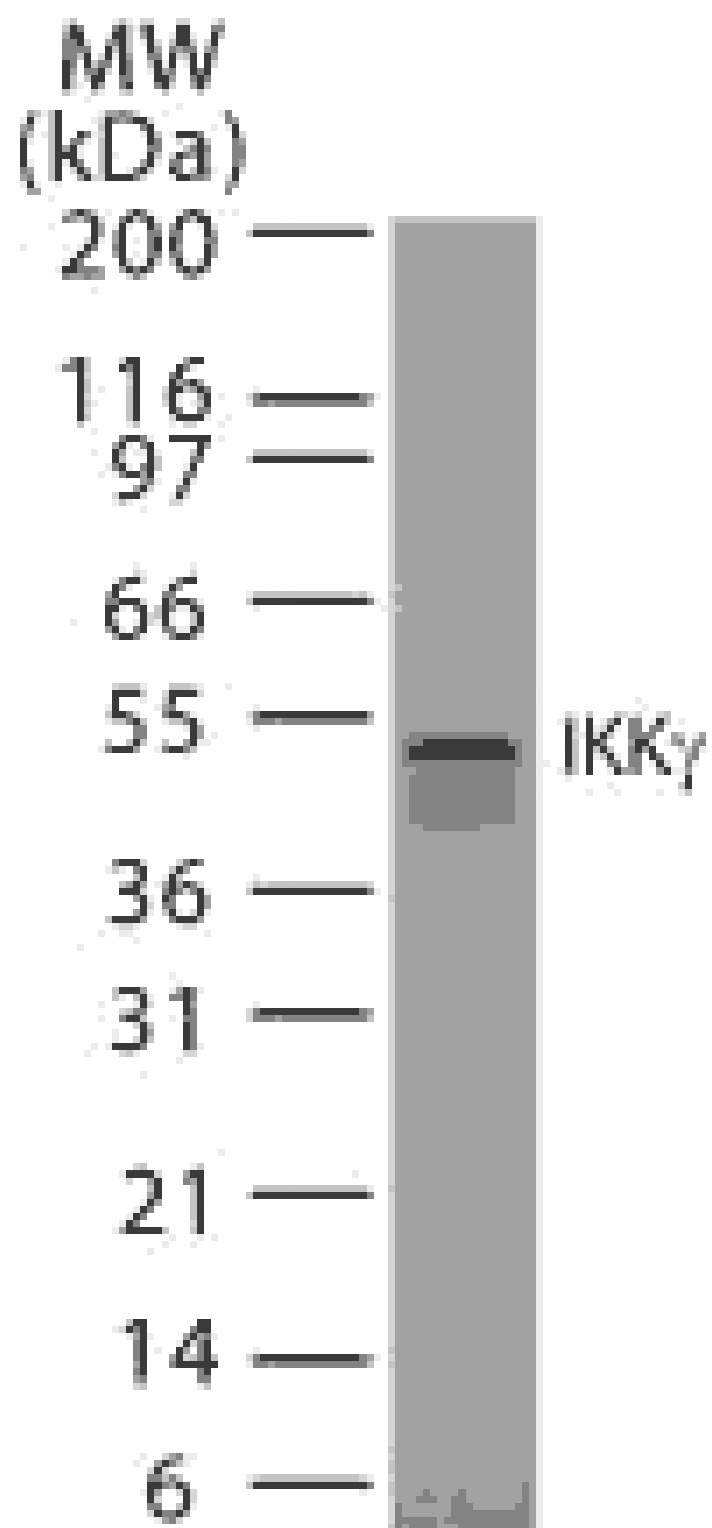
Ordering Information

[Order Online »](#)

ALX-804-424-C100	100 μ g
------------------	-------------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	Stable for 6 months when stored at +4°C.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Inhibitor of nuclear factor κ -B kinase γ , IKK3, NEMO
Application	WB
Application Notes	Detects a band of ~48kDa by Western blot.
Clone	46B844
Crossreactivity	Does not cross-react with mouse IKK γ .
Formulation	Liquid. In PBS containing 0.2% gelatin and 0.05% sodium azide.
Host	Mouse
Immunogen	Recombinant human IKK γ (inhibitor of NF- κ B kinase γ ; IKK3; NEMO).
Isotype	IgG1
Purity Detail	Protein G-affinity purified.
Recommendation Dilutions/Conditions	Western Blot (2 μ g/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human

UniProt ID

Q9Y6K9



Worry-free Guarantee

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-fr@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

This antibody is covered by our **Worry-Free Guarantee**.