

# IKK $\gamma$ (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 $\mu$ g
------------------	-------------

Manuals, SDS & CofA

[View Online »](#)

MW  
(kDa)

200

116

97

66

55

36

31

21

14

6



IKKy

## Handling & Storage

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

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Inhibitor of nuclear factor $\kappa$ -B kinase $\gamma$ , IKK3, NEMO
<b>Application</b>	WB
<b>Application Notes</b>	Detects a band of ~48kDa by Western blot.
<b>Clone</b>	46B844
<b>Crossreactivity</b>	Does not cross-react with mouse IKK $\gamma$ .
<b>Formulation</b>	Liquid. In PBS containing 0.2% gelatin and 0.05% sodium azide.
<b>Host</b>	Mouse
<b>Immunogen</b>	Recombinant human IKK $\gamma$ (inhibitor of NF- $\kappa$ B kinase $\gamma$ ; IKK3; NEMO).
<b>Isotype</b>	IgG1
<b>Purity Detail</b>	Protein G-affinity purified.
<b>Recommendation Dilutions/Conditions</b>	Western Blot (2 $\mu$ g/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
<b>Species Reactivity</b>	Human
<b>UniProt ID</b>	Q9Y6K9

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)