## IKKγ (human) monoclonal antibody (46B844)

This antibody is covered by our Worry-Free Guarantee.

Citations: 1

View Online »

**Ordering Information** 

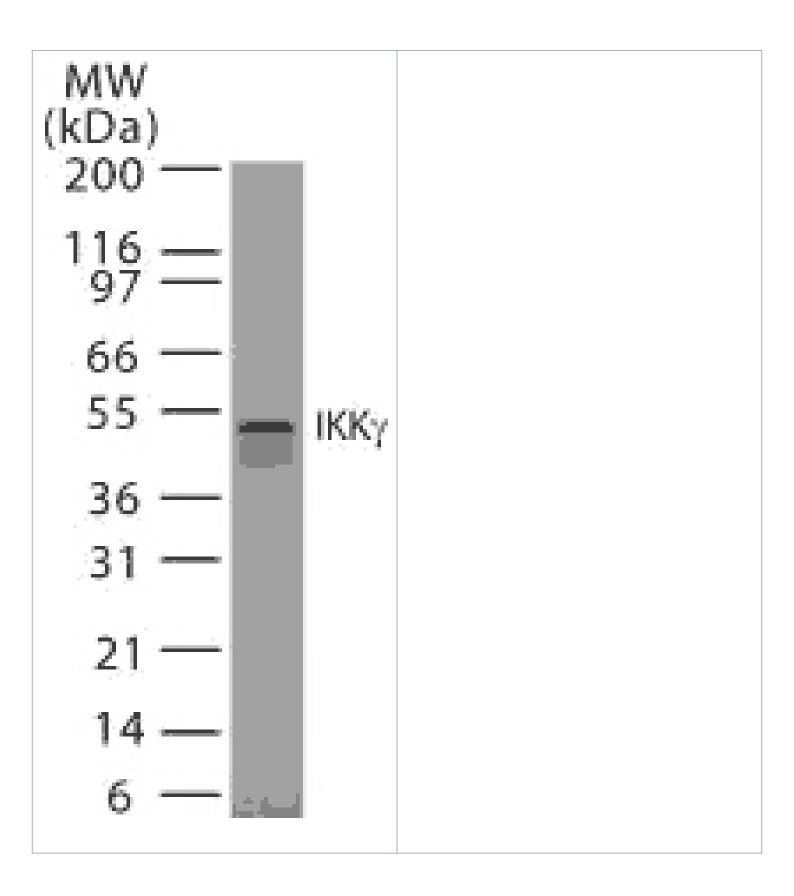
Order Online »

ALX-804-424-C100

100µg

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 IKKy.

Formulation Liquid. In PBS containing 0.2% gelatin and 0.05% sodium

azide.

**Host** Mouse

**Immunogen** Recombinant human IKKγ (inhibitor of NF-κB kinase γ;

IKK3; NEMO).

lsotype lgG1

Purity Detail Protein G-affinity purified.

Recommendation Dilutions/Conditions Western Blot (2µg/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

Last modified: May 29, 2024



info-

eu@enzolifesciences.com