

PRODUCT INFORMATION



IκBζ Polyclonal Antibody

Item No. 13926

Overview and Properties

Contents: This vial contains 100 μg of protein G-purified IgG in 200 μl PBS, with 0.2% gelatin and 0.05% sodium azide.

Synonyms: INAP, MAIL

Immunogen: Synthetic peptide corresponding to mouse IκBζ amino acids 684-699 and 285-298

Species Reactivity: (+) Mouse IκBζ

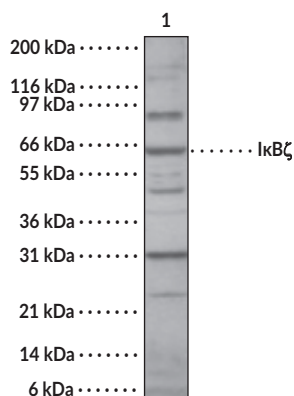
Storage: -20°C (as supplied)

Stability: ≥1 year

Host: Rabbit

Applications: Western blot (WB); the recommended starting concentration is 1-3 μg/ml. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Lane 1: RAW cell lysate

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/15/2024

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

I κ B ζ is an ankyrin repeat-containing nuclear protein that is highly homologous to the I κ B family member Bcl-3. Transcription of I κ B ζ is upregulated by stimulation with TLR ligands, IL-1, and IL-6 in cultured B-lymphocytes and monocytes/macrophages, but only faintly so in T-lymphocytes, fibroblasts, and endothelial cells. I κ B ζ preferentially associates with the NF- κ B subunit p50 rather than p65 and recombinant I κ B ζ proteins inhibit the DNA binding of the p65/p50 heterodimer and the p50/p50 homodimer.

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM