

Product Information

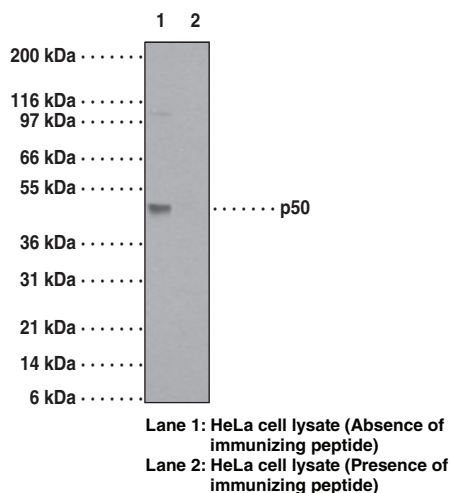


NF-κB (p50) Monoclonal Antibody (Clone 2J10D7)

Catalog No. 13755

- Contents:** This vial contains 100 µg of protein G-purified IgG in 200 µl PBS containing 0.2% gelatin and 0.05% sodium azide.
- Antigen:** A portion of amino acids 150-200 of human NF-κB (p50)
- Host:** Mouse, clone 2J10D7
- Isotype:** IgG_{1κ}
- Cross Reactivity:** (+) Human NF-κB (p50); predicted to react with murine and rat NF-κB (p50)
- Storage:** ≤6 months at 4°C; ≥6 months at -20°C
- Application:** Western blot (WB) and immunohistochemistry (IHC). The recommended starting concentration for WB is 2-5 µg/ml and 5 µg/ml for IHC.

NF-κB is a transcription factor activated by various extra and intracellular stimuli such as cytokines, UV radiation, stress, injury, and by bacterial and viral products. It is involved in regulation of various cellular events including cell growth, differentiation, proliferation, apoptosis, and inflammation. NF-κB1 (p50), a 50 kDa functional sub-unit of NF-κB, is a member of the Rel protein family. It is synthesized as a p105 precursor protein and consists of an N-terminal conserved RHD-region containing a nuclear localization signal, DNA-binding and dimerization domains. NF-κB (p50) forms homodimers or heterodimerizes with p65, forming the functional NF-κB factor. NF-κB (p50) directs the nuclear translocation of NF-κB and is instrumental in its DNA-binding. A pathological role of NF-κB has been suggested in AIDS, hematogenic cancer cell metastasis, and rheumatoid arthritis.



Related Products

NF-κB (p50) Polyclonal Antibody - Cat. No. 13754 • NF-κB (p50-NLS) Inhibitory Peptide Set - Cat. No. 13760 • NF-κB (human p50) Transcription Factor Assay Kit - Cat. No. 10006912 • NF-κB p50 (human recombinant) - Cat. No. 10009818 • NF-κB (human p50/p65) Combo Transcription Factor Assay Kit - Cat. No. 10011223

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 03/15/2010