

NAK1/Nur77 antibody (pAb)

Catalog No: 40982

Isotype: IgG

Application(s): WB

Reactivity: Human, Mouse

Purification: Affinity Purified

Host: Rabbit

Concentration: 0.5 µg/µl

Quantity: 100 µg

Molecular Weight: 84 kDa

Background: NAK1 (Nur77, NR4A1, NP10 or GFRP1) is a member of the steroid-thyroid hormone-retinoid receptor superfamily. NAK1 (Nur77, NR4A1, NP10 or GFRP1) expression is induced by phytohemagglutinin in human lymphocytes and by serum stimulation of arrested fibroblasts. The encoded protein acts as a nuclear transcription factor. Translocation of the protein from the nucleus to mitochondria induces apoptosis. Multiple alternatively spliced variants, encoding the same protein, have been identified.

Immunogen: This NAK1/Nur77 antibody was raised against a peptide corresponding to amino acid residues 251-266 of human NAK1/Nur77.

Buffer: PBS containing 0.05% sodium azide and 0.05% BSA. Sodium azide is highly toxic.

Application Notes:

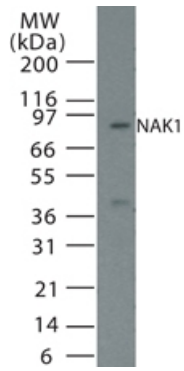
Validated Applications:

WB: 1:500 dilution

For optimal results, primary antibody incubations should be performed at room temperature. The addition of 0.1% Tween 20 to all Blotto solutions may also reduce background. Individual optimization may be required.

Storage and Guarantee: Antibodies in solution can be stored at -80°C for 2 years. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.

**NAK1/Nur77 pAb tested by Western blot.**

NAK1/Nur77 detection by Western blot. The analysis was performed using NIH/3T3 whole-cell extracts and NAK1/Nur77 pAb at a 1:500 dilution.