
Phospho-MAPK3-T202 Polyclonal Antibody

(Catalog #A71504)

Background

The protein encoded by this gene is a member of the MAP kinase family. MAP kinases, also known as extracellular signal-regulated kinases (ERKs), act in a signaling cascade that regulates various cellular processes such as proliferation, differentiation, and cell cycle progression in response to a variety of extracellular signals. This kinase is activated by upstream kinases, resulting in its translocation to the nucleus where it phosphorylates nuclear targets. Alternatively spliced transcript variants encoding different protein isoforms have been described.

Description

Phospho-MAPK3-T202 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

P27361

Purification

Affinity Purification

Immunogen

A synthetic phosphorylated peptide around T202 of human ERK1/ERK2 (NP_002737.2).

Storage

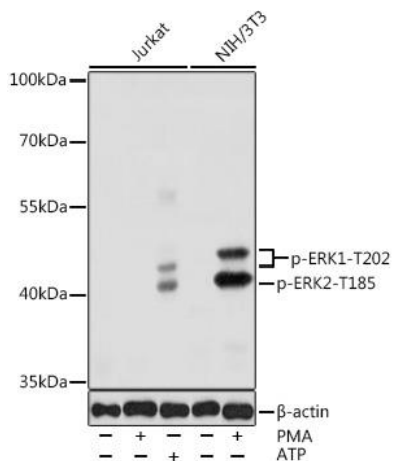
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

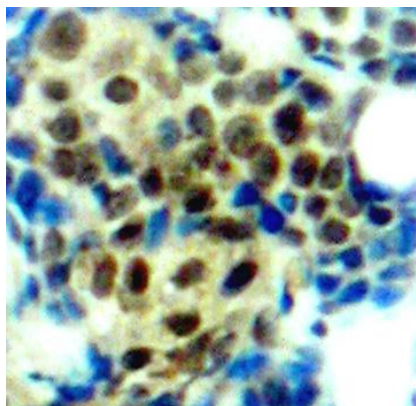
MAPK3; ERK-1; ERK1; ERT2; HS44KDAP; HUMKER1A; P44ERK1; P44MAPK; PRKM3; p44-ERK1; p44-MAPK; mitogen-activated protein kinase 3

Application

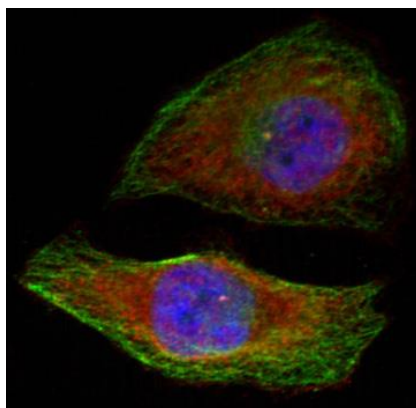
ELISA, WB, IHC, IF; Recommended dilution: WB 1:500 - 1:1000, IHC 1:50 - 1:100, IF 1:20 - 1:100



Western blot analysis of various lysates using Phospho-MAPK3-T202 Polyclonal Antibody at 1:1000 dilution. Jurkat and NIH/3T3 cells were treated by PMA/TPA (200 nM) at 37°C for 30 minutes after serum-starvation overnight. Jurkat cells were treated by ATP(5 mM) at 30°C for 1 hour. Secondary antibody: HRP-conjugated Goat anti-Rabbit IgG (H+L) at 1:10000 dilution. Lysates/proteins: 25µg per lane. Blocking buffer: 3% nonfat dry milk in TBST.



Immunohistochemistry of paraffin-embedded human breast carcinoma tissue, using Phospho-MAPK3-T202 Polyclonal Antibody.



Immunofluorescence analysis of methanol-fixed HeLa cells showing cytoplasmic, nuclear staining using Phospho-MAPK3-T202 Polyclonal Antibody.