

Phospho-RAF1 (S43) Recombinant Monoclonal Antibody [1F7]

(Catalog #A74178)

Background

Serine/threonine-protein kinase linking Ras GTPases and MAPK/ERK cascade. Controls cell fate decisions like proliferation, differentiation, survival, and apoptosis. Initiates MAPK cascade. Phosphorylates key targets including BAD, adenylyl cyclases, PPP1R12A, and TNNT2. Regulates NF- κ B, motility, apoptosis, proliferation, and angiogenesis. Involved in mitochondria-mediated cell protection, Rho signaling, wound healing, and oncogenic transformation.

Description

Phospho-RAF1 (S43) Recombinant Monoclonal Antibody [1F7]. Unconjugated. Raised in: HEK293F Cell.

Formulation

Buffer: Rabbit IgG in phosphate buffered saline, pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.

Specificity

Human

Isotype

Rabbit IgG

Uniprot ID

P04049

Purification

Affinity Chromatography

Immunogen

A synthesized peptide derived from human Phospho-RAF1 (S43)

Storage

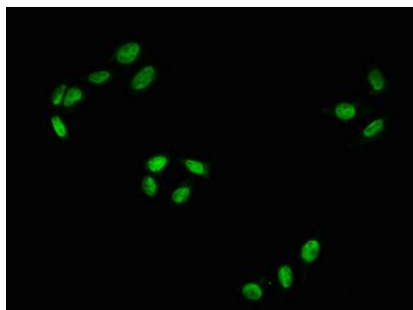
Upon receipt, store at -20°C or -80°C. Avoid repeated freeze.

Alternative Names

RAF proto-oncogene serine/threonine-protein kinase, Proto-oncogene c-RAF, cRaf, Raf-1, RAF1, RAF

Application

ELISA, IF; Recommended dilution: IF:1:20 - 1:200



Immunofluorescence staining of HeLa with p-RAF1 rMab at 1:100, counter-stained with DAPI. The cells were fixed in 4% formaldehyde, permeabilized using 0.2% Triton X-100 and blocked in 10% normal Goat Serum. The cells were then incubated with the antibody overnight at 4°C. The secondary antibody was Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L).