

Cyk-4 phospho Ser170 antibody (pAb)

Catalog Nos: 39265, 39266

RRID: AB_2793211

Isotype: IgG

Application(s): ICC, IF, WB

Reactivity: Human

Volumes: 200 µl, 10 µl

Purification: Protein A Chromatography

Host: Rabbit

Molecular Weight: 32 kDa

Background: Cyk-4 is a member of a GTPase activating protein family that regulates the catalytic function of Rho and Ras small G-proteins. Cyk-4 regulates the initiation of cytokinesis through its interaction with a variety of microtubule organizing proteins. The activity of Cyk-4 is regulated by phosphorylation of serine 170 by Polo-like Kinase 1.

Reference:

Burkard, ME *et al* (2009) PLoS Biol. 2009 7:e1000111.

Immunogen: This Cyk-4 phospho Ser170 antibody was raised against a peptide containing phospho Ser170 of human Cyk-4.

Buffer: PBS pH 7.5 containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

Applications Validated by Active Motif:

ICC/IF: 1:200 - 1:500 dilution

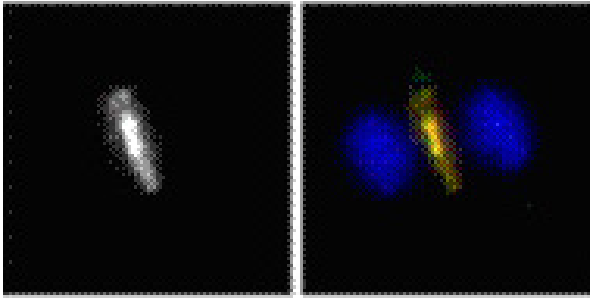
WB: 1:1,000 - 1:2,500 dilution

DB: 1:2000

This Cyk-4 antibody must first be immunoprecipitated and then run on a gel before Western blotting.

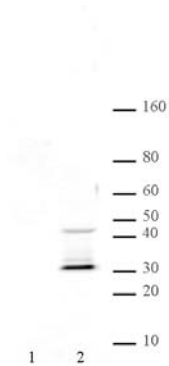
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

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



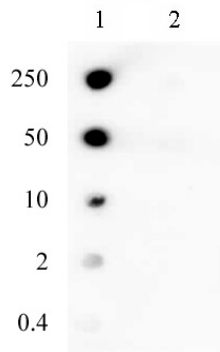
Cyk-4 phospho Ser170 pAb tested by immunofluorescence.

Normal human retinal pigment epithelial cells stained with 39265 (1:300). Left: 39265. Right: merge with DAPI.



Cyk-4 phospho Ser170 Rabbit pAb tested by Western Blot.

Recombinant Cyk-4 fragment (100 ng) untreated (lane 1) and incubated with the kinase Plk-1 (lane 2) was probed with 39265 at 1:2,000.



Cyk-4 phospho Ser170 pAb tested by dot blot to show specificity for serine 170.

Dot blot analysis was used to confirm the specificity of 39265 for phospho Ser170 Cyk-4. The modified and unmodified peptides to the immunogen were spotted onto PVDF and probed with 39265 at a dilution of 1:2,000. The amount of peptide (picomoles) spotted is 250,50,10 and 2 from top to bottom.

Lane 1: phospho Ser170 peptide.
Lane 2: unmodified Ser170 peptide.