



Anti- p27Kip1, monoclonal

CATALOG NO.: 53744

CLONE SPM348

BACKGROUND:

p27Kip1, a candidate tumor suppressor gene, functions as a negative regulator of G1 progression and has been proposed to function as a possible mediator of TGF- β induced G1 arrest.

SOURCE & REACTIVITY:

Mouse anti-p27Kip1 monoclonal antibody was raised against a mouse recombinant p27 protein with the Ig isotype being IgG1 / κ . Anti-p27Kip1 reacts with p27Kip1 at the molecular weight of 27 kDa on western blot. Species reactivity includes human, mouse, and rat, but does not react with cow, while others are not known.

APPLICATION:

The following concentration ranges are recommended starting points for this product.

FC

IFC

WB (Ab 1 μ g/ml for 2hrs at room temperature (RT))

IP (Ab 2 μ g/mg protein lysate)

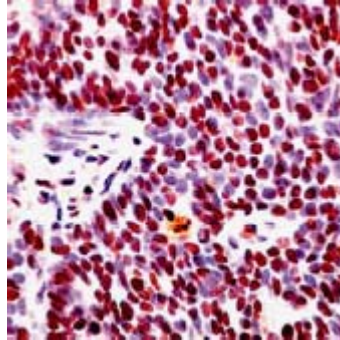
IHC (Formalin/paraffin)

Working dilution: 1:50 for 30 min at RT

Staining of formalin-fixed tissues requires boiling tissue sections in 10mM citrate buffer, pH 6.0 for 10 min followed by cooling at RT for 20 min.

Positive Control: Colon carcinoma, MAD109, Ls174T cells.

Staining Pattern: Nuclear



Human colon ca.
stained with
anti-p27Kip1

This product is for in vitro research purposes only.

STORAGE:

0.25 ml antibody (at 200 μ g/ml), purified from ascites fluid by Protein G, in 1X PBS (pH 7.4) with 0.2% BSA and 0.05% sodium azide. Store at 2-8 $^{\circ}$ C for up to 1 year. Avoid repeated freeze thaw cycles.

