



Anti-Cyclin D2

CATALOG NO: 53422

BACKGROUND:

Cyclin D2 is a G₁ cyclin required for G₁-phase progression and is a strong candidate for a proto-oncogene. Cyclin D2 can phosphorylate pRB when associated with cdk4 and/or cdk6.

SOURCE AND REACTIVITY:

Mouse anti-cyclin D2 monoclonal antibody was raised against purified human recombinant full length cyclin D2 protein with the Ig Isotype being IgG / κ . Anti-cyclin D2 reacts with cyclin-D2 at the molecular weight of 34 kDa on western blot. Species reactivity includes human, mouse, and rat, while others are not known.

APPLICATION:

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

WB

IFC

FC

IP (2 μ g/mg protein lysate)

Positive Control: U-2-OS, C82, RD, Bristol-8, LoVo36 cells.

Cellular localization: Nuclear

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 °C for up to 1 year. Avoid repeated freeze thaw cycles.