



Anti-CsK

CATALOG NO: 53412

BACKGROUND:

C-terminal Src kinase (Csk, p50Csk) is ubiquitously expressed in 50kDa cytoplasmic tyrosine kinase. Csk is involved in the regulation integrin control of cell attachment and shape. Over expression of Csk is localized in focal adhesions and developing embryonic brain, with especially high levels of expression in lymphoid tissues.

SOURCE AND REACTIVITY:

Rabbit anti-Csk polyclonal antibody was raised against a synthetic peptide derived from the C-terminus of human Csk protein. Anti-Csk reacts with Csk at the molecular weight of 50 kDa on western blot. Species reactivity includes human and mouse, while others are not known.

APPLICATION:

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

WB: 2.5 to 5 ug/ml for 2hr, room temperature

Positive Control: Jurkat cells

Cellular Localization: Cytoplasmic



Western blot analysis of Csk in Jurkat cell lysate with anti-Csk

This product is for in vitro research purposes only.

STORAGE:

0.25 ml of affinity purified antibody at 0.2 mg/ml 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.