

anti-hCK2, β -subunit
anti-human casein kinase 2, β -subunit
murine, monoclonal

Cat. No.	Amount
ABD-020	50 μ g

For *in vitro* use only
Quality guaranteed for 12 months
Store at -20°C

Avoid freeze / thaw cycles

Form

Liquid. Supplied lyophilized from PBS.

Handling and use

Reconstitute in 125 μ l ddH₂O
Western Blot: 1 μ g/ml
Immunocytochemistry: 12 μ g/ml

Description

anti-CK2 β is a murine monoclonal antibody (MAb) that was produced against the recombinant human β -subunit of protein kinase CK2 (casein kinase), a pleiotropic and constitutively active serine/threonine protein kinase. The antibody is an affinity-purified IgG (by protein A agarose), the epitope was mapped to (123)GLSDI(127).

Immunogen

Immunogen: Purified recombinant human CK2 β -subunit.

Cross-reactivity

Cross-reactivity: Bovine, mouse, rat, hamster (so far tested).

Selected References:

Nastainczyk et al. (1995) Isolation and characterization of a monoclonal anti-protein kinase CK2 beta-subunit antibody of the IgG class for the direct detection of CK2 beta-subunit in tissue cultures of various mammalian species and human tumors. *Hybridoma*. 14:335.
Sayed et al. (2000) Stress-induced activation of protein kinase CK2 by direct interaction with p38 mitogen-activated protein kinase. *J. Biol. Chem.* 275:16569.