

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

Cat. No.	Amount
ABD-019	100 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 145 μ l ddH₂O
Western Blot: 1 μ g/ml
Immunocytochemistry: 25 μ g/ml

Description

anti-CK2 α is a murine monoclonal antibody (MAb) that was produced against the recombinant human α -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 (319)MEH-PYF(324).

Immunogen

SDS-denatured, purified recombinant human CK2 α -subunit.

Cross-reactivity

Bovine, mouse, rat, hamster (so far tested).

Selected References:

Schmidt-Spaniol et al. (1992) Isolation and characterization of a monoclonal anti CK-2 alpha subunit antibody of the IgG1 subclass. *Hybridoma*. 11:53.
Nastainczyk et al. (2002) Epitope analysis of the MAb 1AD9 antibody detection site in human protein kinase CK2alpha-subunit. *Hybrid. Hybridomics*. 22:87.