

Casein Kinase 2 α polyclonal antibody

The casein kinases are serine/threonine protein kinases involved in the regulation of multiple cell signaling pathways. There are at least seven identified mammalian isoforms in the casein kinase I (CKI) family, which are monomeric and function in cell growth and the replication and repair of DNA. Casein kinase II (CK2) is a ubiquitous tetrameric holoenzyme comprised of two catalytic alpha subunits and two regulatory beta subunits. CK2 enzymes are known to regulate metabolic signal transduction, transcription, and translation.

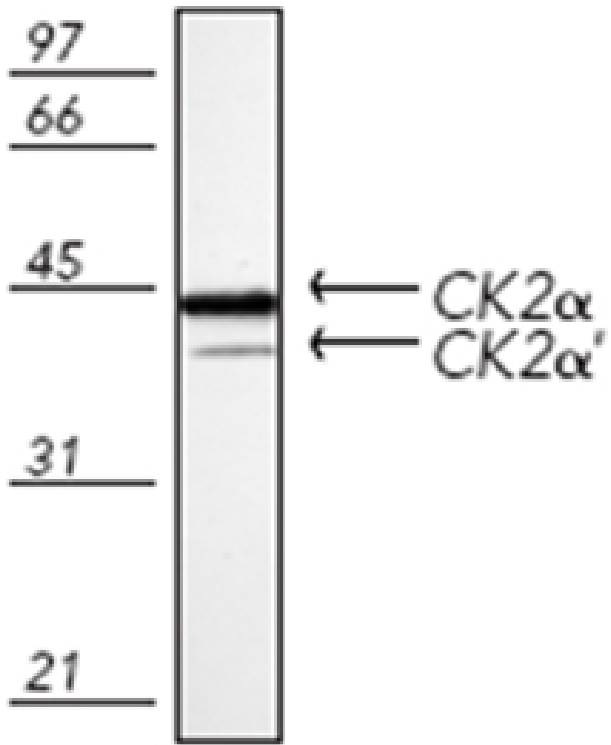
This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 5

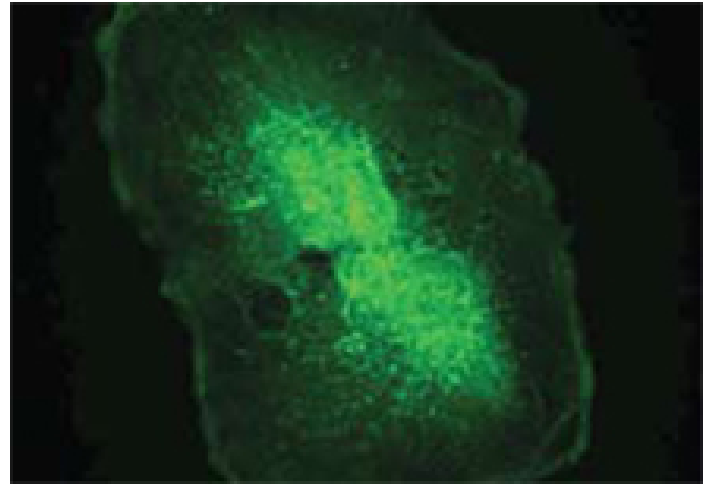
[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of mouse spleen tissue extract, probed with Casein Kinase 2 alpha polyclonal antibody (Prod. No. ADI-KAP-ST010).



Proliferating human Caco-2 cells stained with Casein Kinase 2 alpha polyclonal antibody (Prod. No. ADI-KAP-ST010).

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name

CSNK, CK

Application

IF, IHC, WB

Application Notes

Detects band of ~41kDa and ~45kDa by Western blot. An additional band with an apparent molecular mass of ~38 kDa is detected in brain and heart tissue extracts.

Formulation

Liquid. In borate buffered saline, pH 8.2, containing 25mM sodium borate, 100mM boric acid, 75mM sodium chloride, and 5mM EDTA (disodium salt).

GenBank ID

X70251 (CK2 alpha), M55268 (CK2 alpha')

Host

Rabbit

Immunogen

Synthetic peptide corresponding to a portion of human CK2 α (casein kinase 2 α). The sequence is completely conserved in mouse, rat, rabbit, chicken, Xenopus, zebra fish, Drosophila and C. elegans.

Purity Detail

Immunoaffinity purified.

Recommendation Dilutions/Conditions

Immunohistochemistry (1.7-8.5 μ g/ml)Western Blot (4 μ g/ml, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Source

Purified from rabbit serum.

Species Reactivity

Bovine, Chinook salmon, Drosophila, Guinea pig, Human, Monkey, Mouse, Rabbit, Rat, Sheep

UniProt ID

P68400 (CK2 alpha), P19784 (CK2 alpha')

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)