

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 »](#)

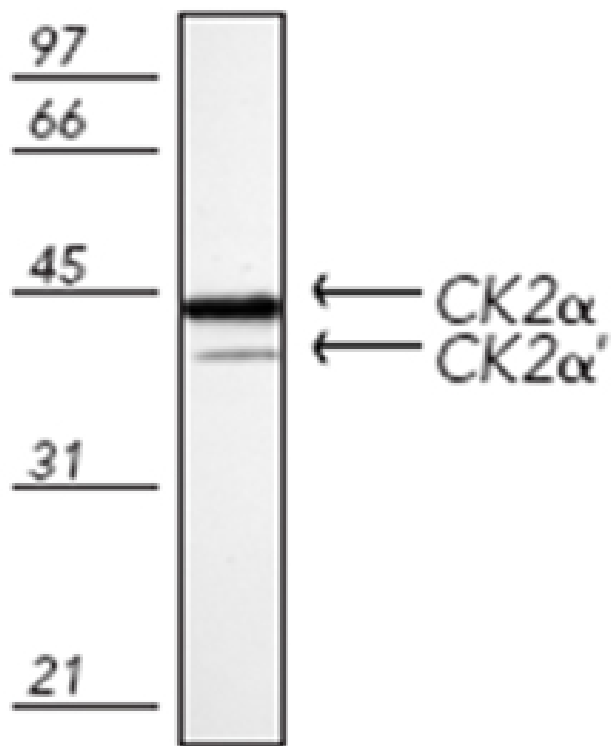
Ordering Information

[Order Online »](#)

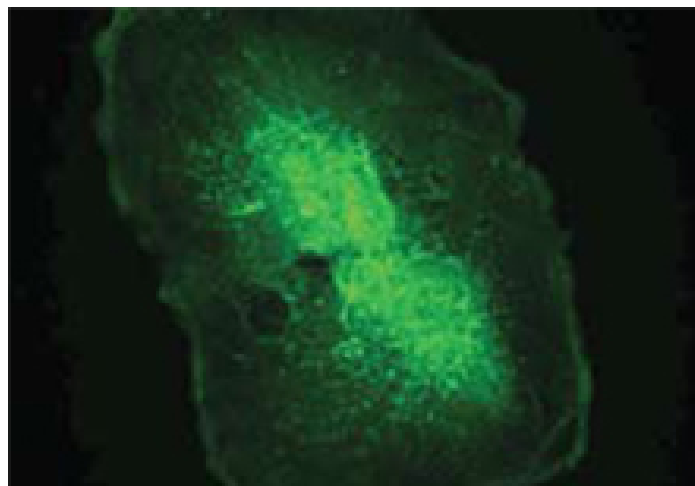
ADI-KAP-ST010-E	100µg
-----------------	-------

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 Immunohistochemistry (1.7-8.5 μ g/ml)Western Blot (4 μ g/ml, ECL)Suggested
Dilutions/Conditions 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')



ENZO LIFE SCIENCES,
INC.
Tel: +1 800.942.0430
info-
usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
Tel: +33 472 440 655
Phone: +41 61 926 8989
info-
eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Tel: +33 472 440 655
info-
be@enzolifesciences.com

France
Phone: +33 472 440 655
fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
de@enzolifesciences.com

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

Worry-Free Guarantee This antibody is covered by our **Worry-Free Guarantee**.