

[pThr²⁸⁶]Ca²⁺

/calmodulin-

dependent protein

kinase II (rat) (α

subunit) monoclonal

This antibody is covered by our [No-Fluorescence Guarantee](#).

antibody (B1)

[Citations](#)

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-802-0090

100 μ l

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name CaM kinase II α subunit, CAMK2 α subunit

Application ELISA, Flow Cytometry, ICC, IF, IP, WB

Clone 22B1

Formulation Liquid. In PBS containing 0.05% sodium azide.

Host Mouse

Immunogen Synthetic peptide corresponding to aa 281-294 (M281HRQET(PSO3)VDCLKKFN294) of the α -subunit of mouse and rat CaM kinase II (Ca²⁺/Calmodulin-dependent Protein Kinase II) thiophosphorylated at (Thr286).

Isotype IgG1

Recommendation Dilutions/Conditions Flow Cytometry (FACS, 1 μ g 106 cells) Immunocytochemistry (1:100) Immunofluorescence (1:100) Western Blot (1:2,000) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Species Reactivity Human, Mouse, Rat

Specificity Recognizes the phosphorylated α -subunit of CaM kinase II. This antibody does not detect the non-phosphorylated α -subunit of rat CaM kinase II nor the ~60kDa β -subunit in either phosphorylation state.

UniProt ID

P11798 (mouse), P11275 (rat)

Worry-free Guarantee

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

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)