

[pThr²⁸⁶]Ca²⁺ /calmodulin- dependent protein kinase II (rat) (α subunit) monoclonal antibody (B1)

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

Citations: 3 [View Online »](#)

Ordering Information [Order Online »](#)

| | |
|------------------|-------|
| ALX-802-009-R100 | 100μl |
|------------------|-------|

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 Flow Cytometry (FACS, 1 μ g 106 cells) Immunocytochemistry
Dilutions/Conditions (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)



ENZO LIFE SCIENCES,
INC.
Phone: +1 942 042 0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
& Luxembourg
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 266 6000
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

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