CREB monoclonal antibody (82B514)

The transcription factor CREB (cAMP response element binding) binds the palindromic octanucleotide TGACGTCA. Members of the ATF/CREB family include CREB-1, CREB-2 (also designated ATF-4), ATF-1, ATF-2, and ATF-3. Phosphorylation of CREB allows recruitment of CREB-binding protein (CBP), a large co-activator that contacts the general transcriptional machinery. CREB-mediated transcription regulates diverse cellular responses, including intermediary metabolism, neuronal signaling, cell proliferation, and apoptosis.

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

ADI-KAM-TF030-E

100µg

Manuals, SDS & CofA

View Online »



Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name cAMP responsive element binding protein

Application WB

Clone 82B514

Formulation Liquid. In PBS containing 0.2% gelatin and 0.05% sodium

azide.

Host Mouse

Immunogen Synthetic peptide corresponding to a portion of human

CREB.

Isotype IgG1

Purity Detail Protein G affinity purified.

Recommendation Dilutions/Conditions Western Blot (2-4µg/ml)Suggested dilutions/conditions

may not be available for all applications. Optimal conditions

must be determined individually for each application.

Species Reactivity Human

UniProt ID P16220

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

