

# 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](#).

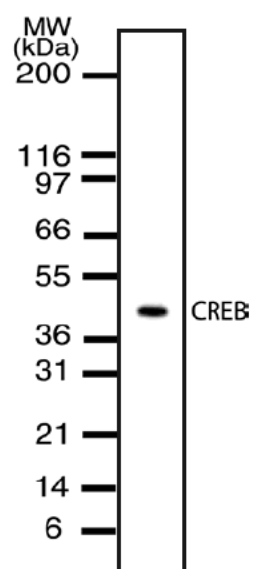
## Ordering Information

[Order Online »](#)

ADI-KAM-TF030-E	100µg
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of  
HeLa cell lysate, probed  
with KAM-TF030.

## 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](#).



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)