

Anti-Human proBDNF Antibody

Mouse Monoclonal antibody

Catalog number

H10001G-MA (100 µg)

Description

Anti-Human proBDNF Antibody is a mouse monoclonal antibody that is specific for the prodomain of human brain-derived neurotrophic factor (proBDNF). It does not recognize mature BDNF.

Immunogen

Purified GST-prodomain protein was used to immunize mice as well as to screen hybridoma clones. Clone mAb287 was selected from the initial screening and further characterized using recombinant proteins, neurons, brain lysates and sections from *bdnf* KO and the wild type littermates as described¹.

Specificity

The anti-proBDNF antibody recognizes the prodomain of human BDNF and does not recognize mature BDNF.

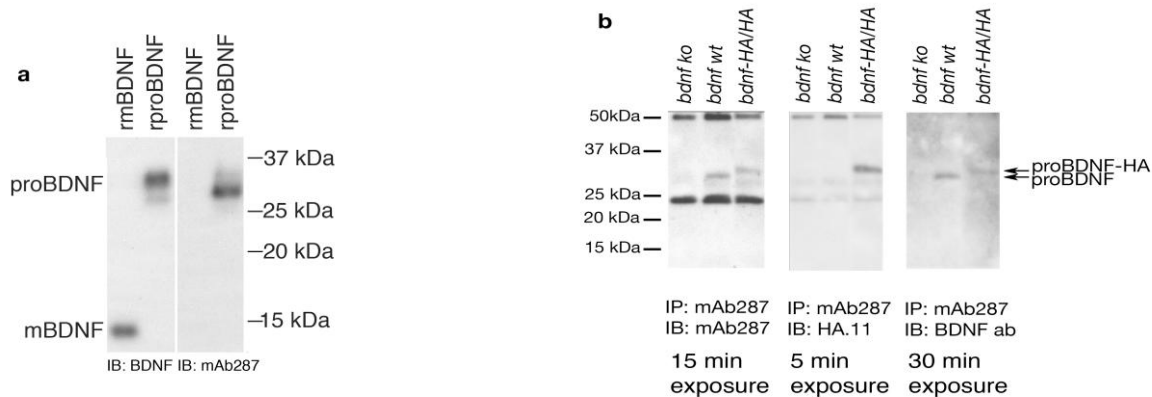


Figure 1. Characterization of anti-prodomain antibody. **(a)** Western blot analysis documenting immunospecificity of monoclonal mAb287 antibody. Purified recombinant mature BDNF (50 ng) and purified recombinant proBDNF (50 ng) were separated using SDS-PAGE and probed with anti-mature BDNF antibody (N-20, Santa Cruz) (left panel) or the monoclonal mAb287 antibody (right panel). **(b)** Detection of proBDNF from brain lysates using mAb287 antibody. The hippocampal lysates of *bdnf*^{-/-}, *bdnf*^{+/+} and *bdnf*-HA/HA mice were immunoprecipitated with mAb287 antibody. After the separation by SDS-PAGE, immunoblotting was performed using either anti-mAb287, anti-HA.11 or anti-mature BDNF antibodies as indicated. Note the lack of cleaved prodomain, which has predicted mass of 14-16 kDa.

Applications

Western blot, immunoprecipitation, immunostaining and affinity purification.

Storage/Stability

Stable for 1 year at -20°C from shipment date. Briefly centrifuge the original vial prior to opening the cap to collect the product at the bottom of the tube.

Starting dilution recommendations

Optimal working conditions should be determined by the user. General recommendations are western blot at 1:5000 to 1:10,000.

Reference

1. Yang, J. et al. *Nature Neuroscience* doi: 10.1038/nn.2244 (2009)

GeneCopoeia, Inc.
9620 Medical Center Drive, #101
Rockville, Maryland 20850
Tel: 301-762-0888 Fax: 301-762-8333
Email: inquiry@genecopoeia.com
Web: www.genecopoeia.com

GeneCopoeia Products are for Research Use Only
Trademarks: GeneCopoeia™ (GeneCopoeia Inc.)

Copyright © 2010 GeneCopoeia Inc.
mAB-proBDNF-122010