

Product Contents

Anti-Human BDNF pAb:

Part No.	Size
G164A	200µg

Description: Anti-Human BDNF pAb is generated in chickens. Anti-Human BDNF pAb is purified using the EGGstract® IgY Purification System (Cat.# G2610).

Applications: For Western blot assay, use antibody at 1µg/ml; for detection and quantitation by ELISA, use antibody at 1µg/ml; for immunocytochemistry, use antibody at 1–10µg/ml; for biological neutralization, use antibody at 10µg/ml.

Immunogen: Human Recombinant BDNF (rhBDNF) (Cat.# G1491).

Formulation: Provided at 0.5mg/ml as frozen liquid in 0.15M NaCl, 0.1M Tris Buffer and 50µg/ml gentamicin.

Storage Conditions: See product information label for storage recommendations. Avoid multiple freeze-thaw cycles and exposure to frequent temperature changes. The performance of this product is guaranteed for 6 months from date of purchase if stored and handled properly.

Note: The storage buffer for Anti-Human BDNF pAb has changed from 0.1M NaCl, 0.01M K₂HPO₄, 50µg/ml gentamicin to 0.15M NaCl, 0.1M Tris Buffer, 50µg/ml gentamicin.

Quality Control Assays

Activity Assays

Biological Activity: A multiwell plate is coated with rhBDNF at a concentration of 0.5µg/ml. Anti-Human BDNF pAb is added to the first row of wells at 10µg/ml, diluted serially to 0.01µg/ml and detected using Anti-Chicken IgY, HRP Conjugate (Cat.# G1351) and TMB substrate. Anti-Human BDNF pAb gives a signal of twofold greater than background at a concentration of ≤0.4µg/ml under these conditions.

Cross-Reactivity: In a sandwich ELISA where Anti-Human BDNF pAb is used as a detect antibody at a 1:500 dilution, the cross-reactivity with 100ng/ml of nerve growth factor, neurotrophin (NT)-3 or NT-4 is less than 3%. Anti-Human BDNF pAb cross-reacts with mouse, rat, monkey, goat, sheep, pig and horse BDNF.

Purity: Anti-Human BDNF pAb is ≥90% pure as determined by 4–20% SDS-PAGE.

Part# 9PIG164

Revised 8/07



AF9PI G164 0807G164



Promega

Promega Corporation

2800 Woods Hollow Road	
Madison, WI 53711-5399	USA
Telephone	608-274-4330
Toll Free	800-356-9526
Fax	608-277-2516
Internet	www.promega.com

PRODUCT USE LIMITATIONS, WARRANTY, DISCLAIMER

Promega manufactures products for a number of intended uses. Please refer to the product label for the intended use statements for specific products. Promega products contain chemicals which may be harmful if misused. Due care should be exercised with all Promega products to prevent direct human contact.

Each Promega product is shipped with documentation stating specifications and other technical information. Promega products are warranted to meet or exceed the stated specifications. Promega's sole obligation and the customer's sole remedy is limited to replacement of products free of charge in the event products fail to perform as warranted. Promega makes no other warranty of any kind whatsoever, and SPECIFICALLY DISCLAIMS AND EXCLUDES ALL OTHER WARRANTIES OF ANY KIND OR NATURE WHATSOEVER, DIRECTLY OR INDIRECTLY, EXPRESS OR IMPLIED, INCLUDING, WITHOUT LIMITATION, AS TO THE SUITABILITY, PRODUCTIVITY, DURABILITY, FITNESS FOR A PARTICULAR PURPOSE OR USE, MERCHANTABILITY, CONDITION, OR ANY OTHER MATTER WITH RESPECT TO PROMEGA PRODUCTS. In no event shall Promega be liable for claims for any other damages, whether direct, incidental, foreseeable, consequential, or special (including but not limited to loss of use, revenue or profit), whether based upon warranty, contract, tort (including negligence) or strict liability arising in connection with the sale or the failure of Promega products to perform in accordance with the stated specifications.

© 1997–2007 Promega Corporation. All Rights Reserved.

EGGstract is a registered trademark of Promega Corporation.

Products may be covered by pending or issued patents or may have certain limitations. Please visit our Web site for more information.

All specifications are subject to change without prior notice.

Product claims are subject to change. Please contact Promega Technical Services or access the Promega online catalog for the most up-to-date information on Promega products.

Part# 9PIG164

Printed in USA. Revised 8/07