

Product Contents

Anti- β III Tubulin mAb:

Part No.	Size
G712A	100 μ g

Description: Anti- β III Tubulin mAb is a mouse monoclonal antibody raised against a peptide (EAQGPK) corresponding to the c-terminus of β III tubulin and is from clone 5G8.

Storage Buffer: Anti- β III Tubulin is supplied in PBS. There are no preservatives added.

Protein Concentration: 1mg/ml.

Storage Conditions: See the Product Information Label for storage recommendations. Avoid multiple freeze-thaw cycles and exposure to frequent temperature changes. See the expiration date on the Product Information Label.

Part# 9PIG712

Revised 2/08



Quality Control Assays

Immunohistochemistry: Anti- β III Tubulin mAb labels cell bodies, including axons and dendrites, but not nuclei of neurons in paraffin-embedded rat brain sections. The antibody is used at a concentration of 0.5 μ g/ml in conjunction with an anti-mouse Cy³ detection system.

Purity: Anti- β III Tubulin mAb is >90% pure, as determined by SDS-PAGE analysis followed by Coomassie[®] blue staining.



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.

© 1998–2008 Promega Corporation. All Rights Reserved.

Coomassie is a registered trademark of Imperial Chemical Industries, Ltd. Cy is a registered trademark of GE Healthcare Bio-sciences.

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# 9PIG712
Printed in USA. Revised 2/08