

Product Contents

Anti-Rat CNTF pAb:

Part No.	Size
G163A	200µg

Description: Anti-Rat CNTF pAb is generated in chickens and purified using the EGGstract® IgY Purification System^(a) (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.

Formulation: Provided at 0.5mg/ml as frozen liquid in 0.1M NaCl, 0.01M K₂HPO₄ and 50µg/ml gentamicin.

Immunogen: Rat Recombinant CNTF (rrCNTF).

Storage Conditions: Store long-term (infrequent use or 1–2 times per month) at –20°C. For daily/weekly use, store at 4°C. 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.

Quality Control Assays

Activity Assays

Biological Activity: A multiwell plate is coated with rrCNTF at a concentration of 0.5µg/ml. Anti-Rat CNTF 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. Specifications require a detection level of twofold greater than background at a concentration of ≤0.4µg/ml under these conditions.

Cross-Reactivity: Anti-Rat CNTF pAb cross-reacts with human, mouse and rat CNTF.

Dot Blot: Anti-Rat CNTF pAb, at a concentration of 0.5µg/ml, must detect 100ng of rrCNTF. No detection is seen for equivalent amounts of interleukin-6 or nerve growth factor.

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

Part# 9PIG163

Revised 4/05



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.

© 1997–2005 Promega Corporation. All Rights Reserved.

^(a)U.S. Pat. Nos. 4, 357,272 and 4,550,019 and U.K. Pat. No. 2,057,451 for recovering purified antibodies from egg yolk.

EGGstract is a trademark of Promega Corporation and is registered with the U.S. Patent and Trademark Office.

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# 9PIG163

Printed in USA. Revised 4/05