

Product Contents

Anti-TGF β ₁ pAb:

Part No.	Size
G122A	100 μ g

Description: Anti-TGF β ₁ is a polyclonal antibody generated in rabbits. The immunogen is biologically active, recombinant TGF β ₁. The antibody is supplied as purified IgG.

Formulation: Anti-TGF β ₁ pAb is supplied in phosphate buffered saline (PBS) containing 0.02mg/ml gentamicin as a preservative.

Protein Concentration: 1mg/ml.

Storage Conditions: Store at 4°C for short-term use. For long-term storage or infrequent use (1-2 times per month), store working aliquots at -20°C. Avoid multiple freeze-thaw cycles and exposure to frequent temperature changes. Performance of this product is guaranteed for at least six months from the date of purchase if stored and handled properly.

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC OR THERAPEUTIC PROCEDURES.

Part# 9PIG122

Revised 11/04



Quality Control Assays

Sensitivity: Using a Western blot assay, a 1:1,000 dilution of this antibody must detect a minimum of 500pg of TGF β ₁ standard.

Specificity: Using a Western blot assay, a 1:1,000 dilution of this antibody shows less than 5% cross-reactivity to TGF β ₂ and TGF β ₃.

Purity: The purity is >90% as determined by SDS-PAGE analysis on a 4-20% reducing gel.



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, 2004 Promega Corporation. All Rights Reserved.

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# 9PIG122
Printed in USA. Revised 11/04