

Product Contents

Anti-Human GDNF pAb:

Part No.	Size
G279A	200µg

Description: Anti-Human GDNF pAb is generated in chickens. It 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.

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

Immunogen: Recombinant Human GDNF (rhGDNF) (Cat.# G2781).

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.

Note: The storage buffer for Anti-Human GDNF 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 rhGDNF at a concentration of 0.1µg/ml (10ng/well). Anti-Human GDNF pAb is added to the first row of wells at 20µg/ml, diluted serially to 0.154µg/ml and detected using Anti-Chicken IgY, HRP Conjugate (Cat.# G1351) and TMB substrate. Specifications require a half-maximal detection at a concentration of <2.5µg/ml under these conditions.

Cross-Reactivity: Anti-Human GDNF pAb cross-reacts with human and mouse GDNF and is presumed to cross-react with rat and Rhesus monkey GDNF based on factor sequence identity across species; it does not cross-react with BDNF, NGF, TGFα and TGFβ.

Purity: Anti-Human GDNF pAb is ≥90% pure as determined by SDS-PAGE.

Part# 9PIG279

Revised 6/08



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.

EGGstract is a registered trademark of Promega Corporation.

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# 9PIG279
Printed in USA. Revised 6/08