

Product Contents

Anti- β -Galactosidase, Purified Monoclonal Antibody:

Part No.	Size
Z378A	100 μ g
Z378B	2mg

Description: This antibody (subclass IgG2a(k)) was purified from ascites of a mouse hybridoma and recognizes *E. coli* β -galactosidase.

Concentration: 2.0–2.5mg/ml in a solution containing 100mM Tris-HCl (pH 8.0), 150mM NaCl, 0.02% NaN₃.

Physical Purity: \geq 95% homogeneous as tested by SDS-polyacrylamide gel electrophoresis followed by Coomassie® blue staining.

Storage Conditions: Store Anti- β -Galactosidase at -20°C . Avoid multiple freeze-thaw cycles and exposure to frequent temperature changes. The performance of this product is guaranteed for at least twelve months from the date of purchase if stored and handled properly.

Part# 9PIZ378

Revised 6/04



Quality Control Assays

Activity Assays

Dot Blot: When this antibody is diluted 1:5,000, it will detect 100ng of both native and denatured β -galactosidase in a dot blot assay using anti-mouse IgG alkaline phosphatase.

Immunoprecipitation: This antibody has been tested for immunoprecipitation of β -galactosidase using Pansorbin® cells. Ten microliters of the monoclonal antibody at 25 μ g/ml will immunoprecipitate \geq 90% of β -galactosidase activity when added to 100 μ l of a 1 μ g/ml solution of β -galactosidase.



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.

© 1999, 2004 Promega Corporation. All Rights Reserved.

Coomassie is a registered trademark of Imperial Chemical Industries, Ltd. Pansorbin is a registered trademark of Calbiochem.

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# 9PIZ378
Printed in USA. Revised 6/04