



## Digestion Module

- Rapid, non-selective release and recovery of intact N-Glycans from up to 50 µg of glycoprotein
- Flexible, high-throughput format
- Compatible with microplate liquid handling on a broad range of automation platforms

Product Code: GS24-RX

**NOTICE:** ProZyme was purchased by Agilent in July 2018. Documents for products and product lots manufactured before August 2019 will contain references to ProZyme. For more information about these products and support, go to: [www.agilent.com/en/contact-us](http://www.agilent.com/en/contact-us).



## TABLE OF CONTENTS

|                                         | page |
|-----------------------------------------|------|
| Kit Contents. . . . .                   | 2    |
| Storage Requirements                    |      |
| Additional Required Reagents/Equipment  |      |
| Optional Reagents and Supplies          |      |
| Safety and Handling. . . . .            | 3    |
| Introduction. . . . .                   | 4    |
| Using the Kit. . . . .                  | 5    |
| Preparation of Samples                  |      |
| Digest Time and Temperature             |      |
| Finishing Options                       |      |
| Protocol. . . . .                       | 9    |
| Getting Started                         |      |
| Preparation of Reagents                 |      |
| Digest                                  |      |
| Analysis of Released N-Glycans. . . . . | 19   |
| Tips & Hints. . . . .                   | 19   |
| References. . . . .                     | 20   |
| Technical Assistance. . . . .           | 21   |
| License to Use. . . . .                 | 22   |
| Trademarks and Tradenames. . . . .      | 23   |
| Product Use and Warranty. . . . .       | 24   |
| Other ProZyme Products & Kits. . . . .  | 24   |
| Ordering Information. . . . .           | 25   |

This product is intended for in vitro research use only.

NOTE: The following suggestions and data are based on information we believe to be reliable. They are offered in good faith, but without guarantee, as conditions and methods of use of our products are beyond our control. We recommend that the prospective user determine the suitability of our materials and suggestions before adopting them on a commercial scale. Suggestions for use of our products or the inclusion of descriptive material from patents and the citation of specific patents in this publication should not be understood as recommending the use of our products in violation of any patent or as permission to license to use any patents of ProZyme, Inc

## KIT CONTENTS

NOTE: We want successful results for our customers, so please read this entire booklet before starting the procedure.

| Item                                  | Qty             |
|---------------------------------------|-----------------|
| WS0283 Digestion (RX) Cartridges      | (24 Cartridges) |
| WS0256-24 Immobilization Reagent Set  | 1 ea            |
| WS0226-24 Denaturation Reagent (8 ml) |                 |
| WS0255-24 Blocking Reagent (1.5 ml)   |                 |
| WS0259-24 Digestion Reagent Set       | 1 ea            |
| WS0278-24 N-Glycanase® (75 µl)        |                 |
| WS0276 25x Digestion Buffer (700 µl)  |                 |
| WS0229 Finishing Reagent (1 ml)       |                 |
| Aluminum Sealing Film (1)             |                 |

## Storage Requirements

This kit is shipped with cold packs and should be stored at 2–8°C on receipt. For best results, equilibrate to ambient temperature prior to use.

## Additional Required Reagents/Equipment

Heater and Incubation Blocks capable of 50–100°C, available from ProZyme as product code GS150  
AssayMAP® Labware: Racks, Receiver Plates and Lids, available from ProZyme as a complete Starter Set (Product Code AM200) or separately in sets of 10.  
Labware: Gilson D-200 tips.  
Centrifuge (capable of 50–1000 x g) and microplate rotor with a height capacity of 44 mm  
Ultrapure, deionized water (Milli-Q® or equivalent)

Acetonitrile (100%, HPLC-grade)  
Pipettors & disposable tips (P5/P10, P200 and P1000)

## Optional Reagents and Supplies

Multichannel pipettors & disposable tips (P5/P10 and P200) (Gilson or equivalent, compatible with Gilson Diamond® D200 pipette tips)  
Pipette basins  
Microplate-compatible, centrifugal evaporator (e.g., SpeedVac® or similar) for optional Finishing of released N-Glycans.  
A rotor that holds two shallow-well microplates and fits into a small-capacity rotary evaporator may be purchased from Thermo Fisher Scientific (Savant Model RH2MP, Thermo Scientific Catalog number 20-548-127).

## SAFETY AND HANDLING

Please refer to the Safety Data Sheets (SDS) posted on ProZyme's website under the component name or Product Code.

<http://www.prozyme.com>

### General Laboratory Procedures

Use powder-free gloves for all sample handling procedures. Ensure that all glass, plasticware and solvents are free of glycosidases and environmental carbohydrates.

## INTRODUCTION

The GlykoPrep® Sample Preparation Platform (GlykoPrep) dramatically streamlines glycoanalysis by facilitating optional protein purification, quantitative deglycosylation and separation of N-Glycans, complete fluorescent labeling and efficient cleanup to reduce excess reagent peaks.

GlykoPrep is modular and can be integrated into any workflow, regardless of throughput or sample type. In order to match any standard sample preparation, Kit components are also available individually as the AssayMAP PA50 (for purification of Fc-containing antibodies only), Digestion Module and dye-specific Labeling & Cleanup Modules.

GlykoPrep is built on AssayMAP technology, microchromatography in a 96-well format, capable of automated high throughput. GlykoPrep may be performed using centrifugation to move liquid through the Cartridges (spin format), or with the Syringe Head on the Agilent AssayMAP Bravo Liquid Handling Workstation (GlykoPrep-plus). Using the spin format with a microplate centrifuge, up to 192 samples can be processed simultaneously with 2 Kits (96-Ct). Important general information for achieving success with the spin format, as well as special tips particular to individual Modules, may be found in the GlykoPrep Guidebook:

<http://www.prozyme.com/documents/TNGP100.pdf>

We also provide modified Microfuge Method instructions useful for those interested in using the spin format to run only a handful of samples with a benchtop microfuge and a PCR heater. Microfuge Method Instruction Manuals may be found on our website:

[http://www.prozyme.com/tech\\_notes.html](http://www.prozyme.com/tech_notes.html)

## USING THE KIT

### Preparation of Samples

#### Sample Quantities

The quantitative binding for the RX Cartridge is 50 µg of most standard glycoproteins.

NOTE: The binding capacity for your specific glycoprotein may need to be determined.

Cartridges are capable of binding more target, but will do so with increasing breakthrough, making the process non-quantitative.

For quantitative loading, prepare an excess of 10% or more sample; denature replicates together and load them individually.

Less than the maximum quantity may be processed, for example, when the sample is available only in limited amounts. The smallest amount of sample that will give good results depends on the sensitivity requirements of the analytical methods and the specific application (e.g., screening vs. QC release).

#### Sample Denaturation

Prior to deglycosylation, the samples are denatured by pre-mixing with Denaturation Reagent. The suggested sample concentration prior to deglycosylation is 1–5 mg/ml, and sufficient reagents have been provided for the standard sample concentration range.

NOTE: If quantitation is desired, pipetting less than 10 µl is not recommended; pipetting smaller volumes introduces variability, especially when samples are highly concentrated. If necessary, dilute the sample to within the 1-5 mg/ml range with Digestion Buffer before starting.

The Kit is useful for very dilute samples without requiring further concentration, by expanding this load step to multiple spins. See the GlykoPrep Guidebook section “Loading.”

When performed in a single spin, the amount loaded to each RX Cartridge should be 10–100 µl. The recommended starting ratio of Denaturation Reagent to sample is 1:1 (v/v).

NOTE: The mixture must be 50% Denaturation Reagent or more.

Example 1:

Sample concentration 1 mg/ml

Sample amount needed: 50 µg

50 µl (50 µg) Sample + 50 µl Denaturation Reagent  
= 100 µl denatured sample

Example 2:

Sample concentration 5 mg/ml

Sample amount needed: 50 µg

10 µl (50 µg) Sample + 90 µl Denaturation Reagent  
= 100 µl denatured sample

The current protocol employs a 5-minute, relatively gentle denaturation, but any custom denaturation may be performed and the subsequent protocol followed as described, as long as no SDS or other detergents are used. Please see the GlykoPrep Guidebook under Digestion Modules or contact us to discuss custom denaturation conditions for your

glycoprotein.

## Digest Time & Temperature

### Time

The Digest procedure has been optimized to deliver deglycosylation of most N-Glycans in 15-60 minutes. The optimal incubation time will vary depending on the specific glycoprotein; those which have proven to be resistant to deglycosylation via conventional enzymatic methods may require longer incubation times (up to 60 minutes). For glycoproteins that are comparatively easy to deglycosylate, such as monoclonal antibodies, a 15-minute incubation is generally sufficient. The selected Incubation Time will be used on page 16.

NOTE: It is critical not to exceed a 60-minute incubation, as the Cartridge resin bed may dry out, yielding uncertain results.

### Temperature

The GS150 Heater and Incubation Blocks are specially designed to provide rapid heat transfer through the Receiver Plate and into the packed bed of each Cartridge. The Incubation Blocks are sold separately (ProZyme Product Code WS0272) and can be used in any standard dry-block heater of the proper size, or pre-heated and used in an oven. Custom Incubation Blocks compatible with robotic systems are also available from ProZyme.

NOTE: If using the Microfuge Method format, PCR heat block substitutes for the GS150 Heater (with Incubation Blocks).

The GS150 Heater (with Incubation Blocks) is set to 50°C for the Digest procedure (deglycosylation and optional Finishing). Please allow a minimum of 1 hour to equilibrate the Blocks

before use. The Incubation Blocks have been designed with a thermometer well in the corner. We have verified that when the thermometer reads 50°C, the temperature in the Cartridge is ~37°C, the optimal temperature for deglycosylation. If using a different heater, confirm the block temperature.

## Finishing Options

The Elution step will differ depending on the dye used to label the N-Glycans. If using InstantDyes™, which require the glycosylamine produced by N-Glycanase digestion, no Finishing is necessary. Elute in Digestion Buffer following Finishing Option A (No Finishing, steps 9a–9d).

If proceeding to a Labeling Module using reductive amination (e.g., the GlykoPrep Rapid-Reductive-Amination™ 2-AB Labeling Module) or another label requiring a free reducing end, or for label-free analysis, elute in Finishing Reagent following Finishing Option B (steps 9–11). Finishing Reagent converts the glycosylamine produced by N-Glycanase digestion to a free reducing end.

## PROTOCOL

### Getting Started

#### Heater Setting

Turn on the GS150 Heater (with 2 Incubation Blocks). Set to 50°C and allow to equilibrate for a minimum of 1 hour.

#### Centrifuge Settings

If the centrifuge does not have x g settings, determine the setting and the specific microplate rotor combination by consulting the operation manual or the manufacturer's website:

\_\_\_\_\_ rpm = 1000 x g

\_\_\_\_\_ rpm = 300 x g

\_\_\_\_\_ rpm = 50 x g

### Preparation of Reagents

#### Digestion Buffer

NOTE: May be prepared up to one week before use. Store at 2–8°C.

25x Digestion Buffer (supplied with the Kit)  
Ultrapure water

Dilute one volume of 25x Digestion Buffer with twenty-four volumes of ultrapure water to obtain Digestion Buffer. Specifically, add 0.2 ml of 25x Digestion Buffer to 4.8 ml of ultrapure water to make 5 ml of Digestion Buffer.

For fewer samples, prepare 100 µl of Digestion Buffer for each sample to be processed.

Cap tightly and vortex on high for 10 seconds to mix.

## Enzyme Solution

NOTE: Should be prepared on the day of use. Store at RT.

N-Glycanase (supplied with the Kit)

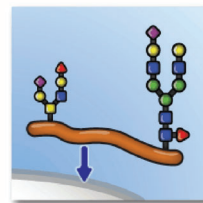
Digestion Buffer (prepared previously)

Spin the N-Glycanase briefly to collect the contents in the base of the vial, and pipette up and down several times to mix prior to use.

In a separate vial, prepare a mixture of 3  $\mu$ l of N-Glycanase and 9  $\mu$ l of Digestion Buffer for each sample to be processed. Pipet up and down several times to mix.

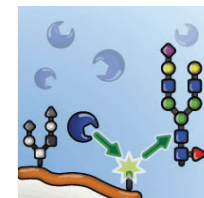
To prepare 24 samples, add 72  $\mu$ l of N-Glycanase to 216  $\mu$ l of 1x Digestion Buffer in a pipette basin. Mix the solution several times by pipette action.

NOTE: The pipette basin requires a minimum of ~100  $\mu$ l volume, so for fewer than 8 samples, do not use a basin.



## Digest

Samples (antibodies or other glycoproteins) are denatured and immobilized.



N-Glycans are released by N-Glycanase and eluted.

## Overview

---

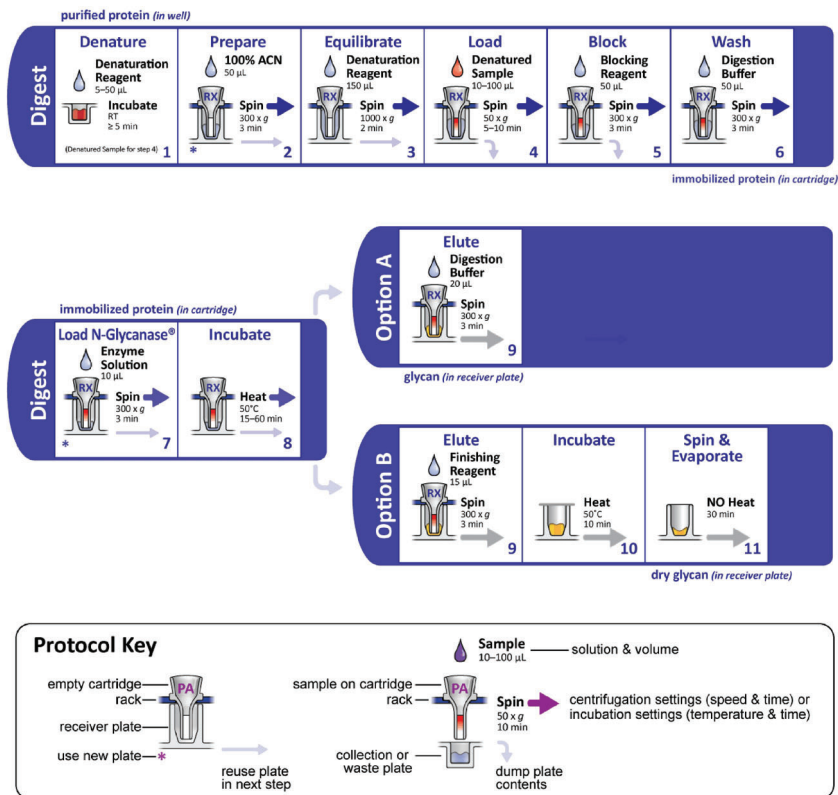
- 1 Denature
- 2 Prepare
- 3 Equilibrate
- 4 Load
- 5 Block
- 6 Wash
- 7 Load N-Glycanase
- 8 Incubate

### Finishing Option A (No Finishing)

- 9 Elute

### Finishing Option B

- 9 Elute (and Finish)
- 10 Incubate
- 11 Spin & Evaporate



## Reagents and other Supplies

### Glycoprotein Samples

NOTE: The quantity of purified sample loaded to the RX Cartridge may contain NO MORE than 50  $\mu$ g total protein, the quantitative binding capacity of the RX Cartridge.

RX Cartridges (supplied with the Kit, 1 per sample)

Prepare two balanced RX Cartridge assemblies  
(Cartridges on Racks on Receiver Plates with Lids).

Denaturation Reagent (supplied with the Kit)

Acetonitrile (100%, HPLC-grade), 50  $\mu$ L/sample

Blocking Reagent (supplied with the Kit)

Digestion Buffer (prepared previously)

Enzyme Solution (prepared previously)

Finishing Reagent (supplied with the Kit)

Aluminum Sealing Film (supplied with the Kit; film may be  
with scissors if using less than the full Kit)

## Procedure

NOTE: An incubation at elevated temperature is required for full deglycosylation. Before beginning, be sure each Incubation Block has equilibrated to 50°C; a thermometer may be placed in the corner well of the Incubation Block to monitor the temperature.

### Denature

- 1.a Add Denaturation Reagent to each sample as described in Sample Denaturation (page 5).
- 1.b Pipet up and down to mix.
- 1.c Incubate at room temperature for at least 5 minutes.

NOTE: Proceed through the Prepare, Equilibrate and Load steps without interruption, as evaporation can lead to airlock.

### Prepare

- 2.a Pipet 50 µl of 100% acetonitrile into the Sample Cup of each RX Cartridge.
- 2.b Spin at 300 x g for 3 minutes; do not empty the Receiver Plates.

### Equilibrate

- 3.a Pipet 150 µl of Denaturation Reagent into the Sample Cup of each RX Cartridge.
- 3.b Spin at 1000 x g for 2 minutes into the same Receiver Plate used for Step 2.b.

NOTE: Do not empty the Receiver Plate prior to loading the

denatured sample.

### Load

- 4.a Load each Denatured Sample into the Sample Cup of an RX Cartridge (see Sample Loading Technique in the GlykoPrep Guidebook).
- 4.b Empty the Receiver Plate and blot with a paper towel.

NOTE: Discard waste Acetonitrile Solution according to waste disposal procedures.

- 4.c Spin at 50 x g until all Sample Cups are empty. The estimated spin time is 5 minutes for volumes between 10 and 50 µl or 10 minutes for volumes up to 100 µl.

### Block

- 5.a Pipet 50 µl of Blocking Reagent into the Sample Cup of each RX Cartridge.
- 5.b Empty the Receiver Plate and blot with a paper towel.
- 5.c Spin at 300 x g for 3 minutes; do not remove the Receiver Plate.

### Wash

- 6.a Pipet 50 µl of Digestion Buffer into the Sample Cup of each RX Cartridge.
- 6.b Spin at 300 x g for 3 minutes.

### Load N-Glycanase

- 7.a Transfer RX Cartridges to fresh Receiver Plates.
- 7.b Pipet 10 µl of Enzyme Solution into the Sample Cup of

each RX Cartridge.

- 7.c Spin at 300 x g for 3 minutes;  
DO NOT DISCARD FLOW-THROUGH.

### Incubate

8. Incubate RX Cartridge assemblies on the equilibrated Incubation Blocks (Heater setting 50°C) for the chosen Incubation Time (not to exceed 60 minutes; see Time, page 7).

NOTE: If N-Glycans are to be labeled using InstantDyes, now would be a good time to prepare the Labeling Reagent.

### Finishing Options

If proceeding to labeling using InstantDyes (e.g., InstantAB™), elute in Digestion Buffer following Option A.

If proceeding to labeling using reductive amination (e.g., Rapid N-Glycan Preparation with 2-AB) or another label that requires a free reducing end, or for label-free analysis, elute in Finishing Reagent following Option B.

#### Option A (no Finishing)

### Elute

- 9.a Remove the RX Cartridge assemblies from the Incubation Blocks.

NOTE: If condensation is apparent, spin at 300 x g for 3 minutes and tap dish gently on the bench top to release Cartridges that may be stuck to the Lid.

- 9.b Pipet 20 µl of Digestion Buffer into the Sample Cup of each RX Cartridge; do not remove Rack from Receiver Plate.

- 9.c Spin at 300 x g for 3 minutes.

- 9.d Remove Cartridges and Rack from the Receiver Plate. The eluted N-Glycans are in the Receiver Plate; DO NOT DISCARD.

Proceed immediately to Labeling.

NOTE: Labeling with InstantDye™ requires the availability of reactive glycosylamine ends, such as those resulting from rapid digestion with N-Glycanase®. Glycosylamine ends spontaneously hydrolyze over time to reducing ends which are incompatible with InstantDye chemistry. To maximize labeling efficiency, Labeling should be performed immediately following collection of the N-Glycans from the GlykoPrep Digestion Module.

NOTE: Retain the RX Cartridges to recover the deglycosylated protein for further analysis (see Tips & Hints).

#### Option B

### Elute (and Finish)

- 9.a Remove the RX Cartridge assemblies from the Incubation Blocks. Keep the heat on for Incubate step.

NOTE: If condensation is apparent, spin at 300 x g for 3 minutes and tap dish gently on bench top to release Cartridges that may be stuck to the Lid.

- 9.b Pipet 15 µl of Finishing Reagent into the Sample Cup of each RX Cartridge; do not remove Rack from Receiver Plate.

- 9.c Spin at 300 x g for 3 minutes.

- 9.d Remove Cartridges and Rack from the Receiver Plate.

The eluted N-Glycans are in the Receiver Plate; DO NOT DISCARD.

#### Incubate

10. Cover the Receiver Plate with Aluminum Sealing Film and incubate on the equilibrated Incubation Block (with Heater still set at 50°C) for 10 minutes.

#### Spin & Evaporate

11. Remove the Aluminum Sealing Film and dry the N-Glycans in a centrifugal evaporator (SpeedVac, heat setting turned to the off position) for 30 minutes or until fully dry. This condenses N-Glycans into a pellet small enough to be dissolved by 5 µl of Labeling Reagent if proceeding to labeling using Rapid Reductive Amination (e.g., the GlykoPrep Rapid-Reductive-Amination 2-AB Labeling Module).

NOTE: The Finishing Reagent must be removed promptly by drying. Store N-Glycans at -20 or -80°C in the dark if further processing will not take place immediately.

NOTE: Retain the RX Cartridges to recover the deglycosylated protein for further analysis (see Tips & Hints).

## ANALYSIS OF RELEASED N-GLYCANS

Use standard techniques, such as LC, CE and MS, to analyze the released N-Glycans.

## TIPS & HINTS

Recovery of the Deglycosylated Protein from the Digestion (RX) Cartridge

Often, the deglycosylated protein is analyzed to evaluate the completeness of deglycosylation using such electrophoretic methods as SDS-PAGE or microfluidic lab-on-a-chip technology. Please contact us for guidelines for eluting your glycoprotein from the RX Cartridge.

## REFERENCES

Visit ProZyme's website for additional information and instructional videos:

<http://www.prozyme.com/glykoprep>

TechNote TNGP100 GlykoPrep Guidebook - General tips, tricks and troubleshooting suggestions when using kits or modules:

<http://www.prozyme.com/documents/TNGP100.pdf>

## TECHNICAL ASSISTANCE

ProZyme is committed to developing rapid, automatable methods for glycan analysis. Call us to discuss products in development.

If you have any questions or experience difficulties regarding any aspect of our products, please contact us:

TOLL FREE **(800) 457-9444** (US & CANADA)  
PHONE **(510) 638-6900**  
FAX **(510) 638-6919**  
E-MAIL **info@prozyme.com**  
WEB **www.prozyme.com**

ProZyme values customer opinions and considers customers an important source for information regarding advanced or specialized uses of our products. We encourage you to contact us. We welcome your suggestions about product performance or new applications and techniques.

## LICENSE TO USE

Purchase of this ProZyme product (the "Product") from ProZyme or one of its authorized distributors comes with a limited, non-transferable Use License to the purchaser ("Purchaser") under ProZyme patents (the "ProZyme Patents"), which Use License is subject to the following terms and conditions.

1. The Use License is only for Purchaser's in-house, in-vitro use of the purchased Product. Purchaser is not licensed to use, and shall not use the Product in any in vivo application and, for avoidance of doubt, is specifically prohibited from any use of the Product in humans.
2. The Product shall not be made available by Purchaser to any third party in any form, separately or in combinations, whether for any monetary or other consideration or at no charge, with the following exception: the Product may be made available to a third party who (a) agrees to be bound by all the terms and restrictions of this Agreement and (b) agrees to use it only for purposes of evaluation.
3. This Use License does not include the right to use the Product to perform for, or offer commercial services of any kind to, a third party, such as, without limitation, reporting the results of analyses or other activities to the third party for a fee or other commercial consideration. Persons wishing to perform such services must obtain a separate Commercial Services License from ProZyme, Inc.
4. Purchaser is responsible for qualification of the Product for the Purchaser's specific use.

Certain ProZyme products include components or reagents obtained under license from other companies. Purchase of these ProZyme products from ProZyme or one of its authorized distributors includes a license from the originating company to use the component(s) or reagent(s), but only (1) with the other materials included in the ProZyme product, and (2) for the purposes specified on ProZyme's product, on its packaging, or in the accompanying literature.

For a complete list of ProZyme Patents, please visit our website at <http://www.prozyme.com/patent.html>

The AssayMAP technology is licensed from Agilent Technologies.

## TRADEMARKS AND TRADENAMES

ProZyme, Glyko, GlykoPrep, InstantDye, Rapid-Reductive-Amination, InstantAB and N-Glycanase are trademarks or registered trademarks of ProZyme, Inc. in the United States and other countries.

Diamond is a registered trademark of Gilson, Inc, Middleton, WI, USA.

Milli-Q is a registered trademark of Millipore Corporation in the United States and/or other countries.

SpeedVac is a registered trademark of Thermo Fisher Scientific or its affiliates.

## PRODUCT USE AND WARRANTY

Terms and conditions of sale may be found at:

<http://www.prozyme.com/terms.html>

## OTHER PROZYME PRODUCTS & KITS

A wide variety of glycoanalysis products are available from ProZyme. A complete listing is accessible on our website by clicking on Glyko® Tools for Glycobiology:

<http://www.prozyme.com>

## ORDERING INFORMATION

For North American destinations:

Telephone orders may be placed between 8:00 am and 5:00 pm Pacific Time. Telefax or e-mail orders may be sent or messages recorded anytime.

TOLL FREE (800) 457-9444 (US & CANADA)

PHONE (510) 638-6900

FAX (510) 638-6919

E-MAIL [info@prozyme.com](mailto:info@prozyme.com)

Outside North America:

A list of ProZyme's distributors, with contact information, may be found at:

<http://www.prozyme.com/distributors.html>

If there is no distributor in your area, instructions for placing an international order may be found at:

<http://www.prozyme.com/ordering.html>



3832 Bay Center Place  
Hayward, CA 94545-3619  
USA

TOLL FREE (800) 457-9444  
PHONE (510) 638-6900  
FAX (510) 638-6919

E-MAIL [info@prozyme.com](mailto:info@prozyme.com)  
WEB [www.prozyme.com](http://www.prozyme.com)  
Rev. AD