

Product Information

GlycoProfile[®] 2-AB Labeling Kit

Product Code **PP0520**
 Store at Room Temperature

TECHNICAL BULLETIN

Product Description

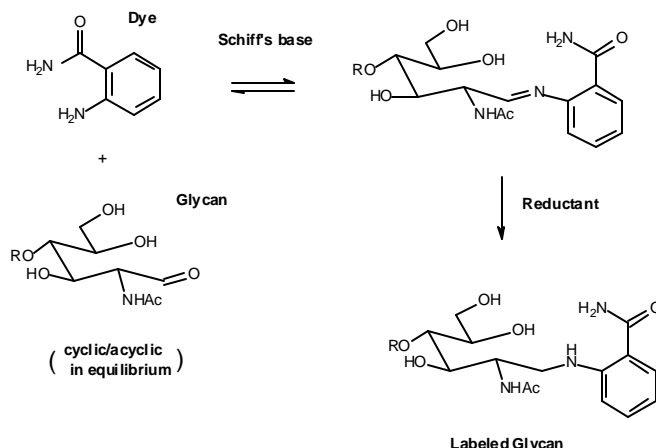
Glycan analysis has become an increasingly critical aspect of glycomics and proteomics, as the role of glycoproteins in cell signaling, cell adhesion, immune response, and disease states is made clear through ongoing research. Because glycans tend to have low spectral activity in both UV and visible light, it is often necessary to label them in order to enhance detection.

The fluorophores 2-AB (2-aminobenzamide) and 2-AA (anthranilic acid or 2-aminobenzoic acid) provide valuable tools for glycan analysis, due to their sensitivity and stability when bound to glycans.¹

Samples containing mixed pools of glycans can often be detected at picomolar concentrations. While 2-AB is somewhat less sensitive than 2-AA, it is suitable for downstream glycan analysis by HPAE (high-performance anion-exchange), HPLC, and ESI-MS (electrospray mass spectroscopy) methods.²⁻⁸ Although suitable for the same applications, 2-AA is better suited for SDS-PAGE.

Glycan binding to both dyes is very robust with no degradation during post-cleanup analysis. Labeling can be performed on either purified or pooled samples, including a variety of sources, such as N-linked, O-linked, and GPI anchored glycans. For samples containing sialated oligosaccharides, sialic acid loss is negligible. Other commonly used methods such as radiolabels, antibody labels, and various probes do not display the stability, flexibility, and ease of use observed with 2-AA and 2-AB.

Glycans with a free reducing sugar exist in equilibrium between the cyclic (closed ring) and acyclic (open ring) structures. A stable Schiff's base is formed when the carbonyl carbon of an acyclic reducing sugar is linked to the amine moiety of the dye in a nucleophilic manner. Following formation of the Schiff's base, the resulting imine group is reduced using sodium cyanoborohydride, resulting in a stable labeled glycan.



Components

The kit contains sufficient reagents for labeling up to 36 samples. Two sets of components have been provided. Each set is sufficient for labeling up to 18 samples based on a 5 µl reaction volume (to accommodate use with the GlycoProfile Glycan Cleanup Cartridges, Product Code G 8169). Prior to the clean up step, larger samples can be split accordingly.

Mixed glycan samples should contain between 100 picomoles to 50 nanomoles of purified glycans. With a single pure glycan, as little as 5 picomoles may be labeled and detected in subsequent HPLC analysis.

2-AB (2-Aminobenzamide) (Product Code A 9478)	2 X 5 mg
DMSO (Dimethyl sulfoxide) (Product Code D 4942) 350 µl per vial	2 X 1 vial
Acetic Acid, Glacial (Product Code A 9353)	2 X 200 µl
Reductant (Sodium Cyanoborohydride) (Product Code R 5153)	2 X 6 mg

Equipment Recommended but Not Provided

GlycoProfile Glycan Clean-Up Cartridges
(Product Code G 8169)

Precautions and Disclaimer

This product is for R&D use only, not for drug, household or other uses. For in vitro research use only. Please consult the Material Safety Data Sheet for information regarding hazards and safe handling practices.

Labeling of glycans with 2-AB is covered under US Patent No 5,747,347 and its foreign equivalents.

It is recommended that the entire Technical Bulletin be read prior to starting the procedure.

Storage/Stability

The kit, as supplied, is stable for at least one year when stored unopened at room temperature. The 2-AB reagent (Product Code A 9478) is light sensitive and must be stored in the dark. Labeled glycans should be stored at -20°C in the dark. For optimal results, the labeling reagent should be used within one hour of component reconstitution and preparation.

Procedure

Glycan samples may be prepared by enzymatic deglycosylation or chemical deglycosylation via hydrazinolysis. For mixed pools of glycans, the optimal sample size is 100 pmole to 50 nmole. For purified samples of a single glycan, much lower amounts may be labeled. As little as 5 pmole of a single purified glycan be detected in post-labeling analysis by HPLC.

1. Prior to labeling, glycan samples should be purified to remove protein, peptides, salts, detergents, and any additional contaminating substances.
2. Transfer the glycan sample solutions into the reaction vials and dry completely. Excess moisture will have negative effects on labeling and stability.
 - a. Optimal results are obtained when samples are dried by centrifugal evaporation. Use caution with lyophilization. Ensure that each sample dries as a concentrated mass at the bottom of the vials.
 - b. Do not expose samples to extreme pH or high temperatures.
 - i. Temperatures greater than 28°C and/or low pH will result in desialylation.
 - ii. Exposure to high pH may result in epimerization of reducing sugars.
3. Prepare the Labeling Solution immediately prior to labeling.
 - a. Tap or briefly spin down component vials to avoid loss of reagent in the cap or on the walls of ampules. To open ampules, hold both the body and the top of the ampule, then gently, but firmly, snap open at the colored break-ring, directing the break away from your body.
 - b. Add $150\ \mu\text{l}$ of Acetic Acid, Glacial (Product Code A 9353) to the entire ampule of DMSO (Product Code D 4942). Note that DMSO forms a semi-solid at cooler temperatures, hence warming may be required.
 - c. Add $100\ \mu\text{l}$ of the DMSO/Acetic Acid mixture (step 3b) to the entire vial of 2-AB (Product Code A 9478). Mix until dissolved; vortexing may be required.
 - d. Add the entire volume from step 3c to the vial of Reductant (Product Code R 5153). Mix until completely dissolved. If insoluble particulates remain, the solution may be heated to 65°C for up to 3 minutes. Any remaining particulates can be dissolved by adding $10\ \mu\text{l}$ of water, (HPLC grade or better). This is the Labeling Solution and should be protected from light. It is stable for 1 hour.
4. Add $5\ \mu\text{l}$ of the Labeling Solution (step 3d) to each dried glycan sample. Cap and mix thoroughly. Tap or spin down to collect the dissolved sample in the bottom of each vial.
5. Incubate glycan samples for 3 hours at 65°C in a heating block, dry oven, or sand tray. Avoid moist heat, if possible. If a water bath or moist heat is used, the vials must be tightly sealed.

Note: Incubation periods of between 2 and 4 hours will not significantly affect results.
6. If insoluble particulates are present, vortex the pre-heated samples from step 5 for 30 minutes.
7. Following incubation, briefly spin the vials to re-collect each sample at the bottom of the reaction vial. Allow the samples to cool to ambient temperature prior to proceeding with post-labeling analysis.
 - a. Extended periods of time between incubation and analysis may result in desialylation of labeled glycans and consequently should be avoided.
 - b. Sample clean up using GlycoProfile Glycan Clean-Up Cartridges (Product Code G 8169) is recommended to remove excess dye and labeling reagents. A purification protocol is included with the cartridges.

Analysis of 2-AB labeled glycans

The 2-AB reagent has an excitation range of 200-450 nm, with a peak excitation at 330 nm. The emission range for 2-AB is 300-750 nm, with a peak emission at 420 nm.

Once the glycans have been labeled, a variety of methods exist to analyze them. The most common techniques employ fluorescent detection after separation by HPLC or CE. These include separation by ion exchange (i.e. high-performance anion exchange), normal phase HPLC, and size exclusion chromatography.

Labeled glycans can also be detected using SDS-PAGE and mass spectrometry. The dye linkage is stable to both of these methods. Mass spectrometry can be performed with either an electrospray ionization (ESI) or by matrix assisted laser absorption ionization (MALDI) ion source. See References for more information.

Troubleshooting Guide

Estimating sample size	The glycan content of an unknown sample may be estimated. The average glycosylation of proteins is 2-5% with an average glycan molecular weight of 1,000 Da. A 1 mg sample of glycoprotein contains ~50 µg of glycans.
Poor dye incorporation	Improper temperature control during incubation. Ensure that samples remain at 65 °C throughout the entire incubation period by preheating heat blocks. Do not over or under heat. Dye incorporation declines sharply at incubation temperatures below 65 °C. Higher temperatures will not increase dye incorporation, but will contribute to desialylation.
	Incomplete solubilization. Glycans must be completely solubilized for maximum labeling efficiency. Ensure thorough mixing at each stage of the procedure. Vortexing for 30 minutes during the incubation period may be necessary (see Procedure, step 6).
Contamination	Sample may contain contaminants that interfere with dye incorporation. Ensure that glycans have been properly purified prior to labeling.
	Degraded Labeling Solution. The Labeling Solution must be freshly prepared immediately prior to use. The Labeling Solution begins to degrade after one hour.
Lower than expected concentration of glycans	Loss or absence of free reducing sugar. The aldehyde group on the free reducing sugars must be available for 2-AB conjugation. Glycans that have been conjugated at their reducing termini (glycolipids, previously labeled glycans, glycopeptides) leave no open site for conjugation to occur.
	Loss of glycans or incomplete binding of glycans to membrane during clean-up process. Refer to the technical bulletin provided with the GlycoProfile Glycan Clean-up Cartridges (Product Code G 8169) for troubleshooting guide and recommended solutions.
Desialylation of glycans	Exposure to acidic conditions or heating at temperatures greater than 65 °C may result in desialylation.
High background, non-glycan fluorescence	This may indicate the presence of an aldehyde-bearing contaminant that readily reacts with 2-AB dye. Ensure adequate clean-up procedure prior to labeling.

Related Products

Enzymatic Deglycosylation Kit (Product Code EDEGLY)
 PNGase F (Product Code P 7367)
 O-Glycosidase (Product Code G 1163)

References:

1. Bigge, J.C., *et al.*, Non-selective and efficient fluorescent labeling of glycans using 2-aminobenzamide and anthranilic acid. *Anal. Biochem.*, **230**, 229-238 (1995).
2. Merry, A.A., *et al.*, Recovery of intact 2-aminobenzamide-labeled O-glycans released from glycoproteins by hydrazinolysis. *Anal. Biochem.*, **304**, 91-99 (2002).
3. Delaney, J., and Vouros, P., Liquid chromatography ion trap mass spectrometric analysis of oligosaccharides using permethylated derivatives, *Rapid Commun. Mass Spectrom.*, **15(5)**, 325-334 (2001).
4. Rudd, P.M., *et al.*, Glycoproteins: rapid sequencing technology for N-linked and GPI Anchor glycans, *Biotechnol. Genet. Eng. Rev.*, **16**, 1-21(1999).
5. Anumula, K.R., and Dhume, S.T., High resolution and high sensitivity methods for oligosaccharide mapping and characterization by normal phase high performance liquid chromatography following derivitization with highly fluorescent anthranilic acid, *Glycobiology*, **8**, 685-694 (1998).
6. Hardy, M.R., *Glycan labeling with the fluorophores 2-aminobenzamide and anthranilic acid*, *Techniques in Glycobiology*, Dekker, Inc. (New York, NY: 1997).
7. Townsend, R., *et al.*, Multimode high-performance liquid chromatography of fluorescently labeled oligosaccharides from glycoproteins, *Anal. Biochem.*, **239**, 200-207 (1996).
8. Guile, G.R., *et al.*, A rapid and high-resolution high-performance liquid chromatographic method for separating glycans mixtures and analyzing oligosaccharide profiles, *Anal. Biochem.*, **240**, 210-226 (1996).

LICENSE TO USE

Purchase of this 2-AB Labeling Kit grants a Use Sublicense under Glyko's 2-AB patents. By accepting delivery of the GlycoProfile 2-AB Labeling Kit [Material(s)] and by subsequently using it in glycan analysis, Recipient agrees to be bound by the following terms and restrictions:

1. A Use Sublicense is granted Recipient for in-house use of Material(s) only.
2. The Material(s) will not be made available by Recipient to any outside parties in any form, separately or in combinations, for any monetary or other consideration or at no charge, except that the Materials may be made available to outside parties who agree to be bound by all the terms and restrictions of this Agreement for purposes of evaluation only.
3. Recipient will not make commercial use of the Material(s) unless it first secures a license agreement for such commercial use.
4. Recipient is solely responsible for qualification of the products for the Recipient's specific use.
5. The Material(s) will not be used *in vivo* in humans.

Labeling of glycans with 2-AB is covered under US Patent No 5,747,347 and its foreign equivalents .

JP,MAM 04/05-1

Sigma brand products are sold through Sigma-Aldrich, Inc.

Sigma-Aldrich, Inc. warrants that its products conform to the information contained in this and other Sigma-Aldrich publications. Purchaser must determine the suitability of the product(s) for their particular use. Additional terms and conditions may apply. Please see reverse side of the invoice or packing slip.