

# Agilent AdvanceBio Gly-X InstantPC

## Quick N-glycan Sample Preparation

Agilent AdvanceBio Gly-X InstantPC reagents and consumables let you prepare samples for analysis including N-glycan release, labeling and cleanup, in as little as 1 hour using this simple step by step process.



You will need these additional items to prepare your samples:

### Lab supplies and equipment

Pipettes and pipette tips  
Eppendorf tubes  
Graduated cylinder (250 mL or larger)  
Glass storage vessel (250 mL or larger)  
Thermocycler

### Additional reagents

Formic acid  
Acetonitrile



**1. Prepare N-glycanase working solution.** Pipette 1.2  $\mu$ L of N-glycanase to one Eppendorf tube per sample. **Make sure to change the pipette tip between transfer steps.**  
Duration: 1 min      Total time: 1 min



**2. Add 2  $\mu$ L of Digestion Buffer to each tube. Make sure to change the pipette tip between transfer steps.** Return the N-glycanase and Digestion Buffer to 4 °C.  
Duration: 1 min      Total time: 2 mins



**3. Prepare InstantPC Dye solution.** Add 150  $\mu$ L of InstantPC Dye Solvent (green cap) to the InstantPC Dye (orange cap). Vortex until dissolved. The dye is a dry solid that goes into solution after solvent is added. Return the InstantPC Dye Solvent to -20 °C.  
Duration: 5 mins      Total time: 7 mins



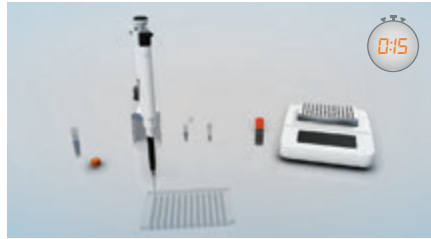
**4. Prepare Load/Wash solution.** Add 6 mL of formic acid to a 6 mL glass graduated cylinder.  
Duration: 3 mins      Total time: 10 mins



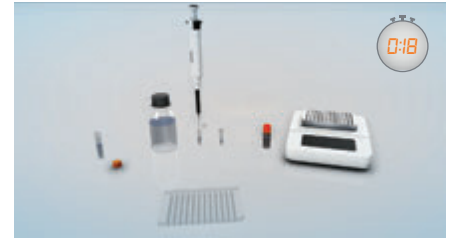
**5. Bring the volume up to 240 mL with acetonitrile.**  
Duration: 3 mins      Total time: 13 mins



**6.** Transfer the solution in the graduated cylinder to a glass storage vessel. Cap tightly and swirl to mix.  
Duration: 1 min      Total time: 14 mins



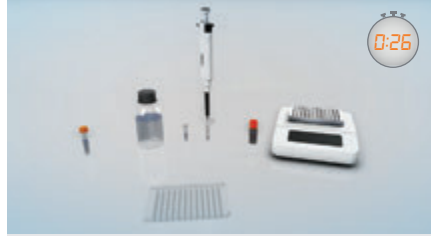
**7. Gly-X deglycosylation.** Pipette 2  $\mu$ L of Gly-X Denaturant into wells of a Gly-X Deglycosylation Plate for each sample you are preparing.  
Duration: 1 min      Total time: 15 mins



**8.** Add 20  $\mu$ L of glycoprotein sample to each well. Mix well using the pipette. **Make sure to change the pipette tip between transfer steps.**  
Duration: 3 mins      Total time: 18 mins



**9.** Tap plate on the benchtop to collect samples at bottom of wells. Put the Gly-X Deglycosylation Plate on heater and incubate uncovered at 90  $^{\circ}$ C for 3 min.  
Duration: 5 mins      Total time: 23 mins



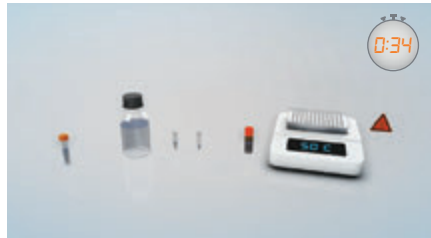
**10.** Remove plate and let sit on the bench at room temp for 2 min. Add 2  $\mu$ L N-Glycanase working solution to each sample well on the plate. Mix well using the pipette.  
Duration: 3 mins      Total time: 26 mins



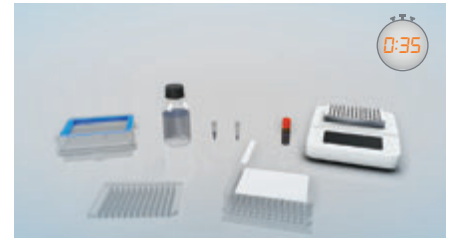
**11.** Tap plate on the benchtop to collect samples at bottom of wells. Put the Gly-X Deglycosylation Plate on heater and incubate uncovered at 50  $^{\circ}$ C for 3 min.  
Duration: 3 mins      Total time: 29 mins



**12. Instant PC labeling.** Add 5  $\mu$ L InstantPC Dye solution to each sample well on the plate. Mix well using the pipette.  
Duration: 3 mins      total time: 32 mins



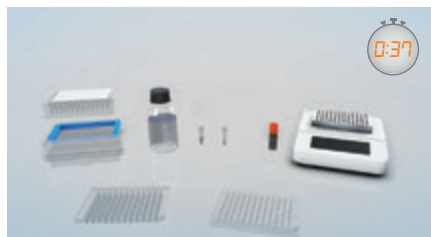
**13.** Tap plate on benchtop to collect samples at bottom of wells. Put Gly-X Deglycosylation Plate on heater and incubate uncovered at 50  $^{\circ}$ C for 1 min.  
Duration: 2 mins      Total time: 34 mins



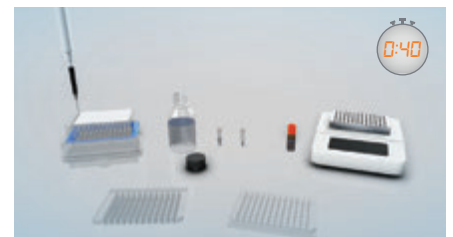
**14. InstantPC cleanup.** Carefully remove the white caps on the Gly-X Cleanup Plate.  
Duration: 1 min      Total time: 35 mins



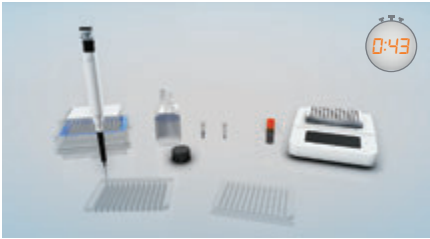
**15.** Install the waste tray in the vacuum manifold.  
Duration: 1 min      Total time: 36 mins



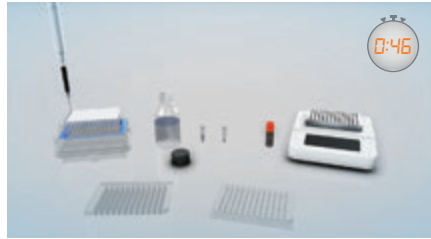
**16.** Put the Gly-X Cleanup Plate on top of the vacuum manifold.  
Duration: 1 min      Total time: 37 mins



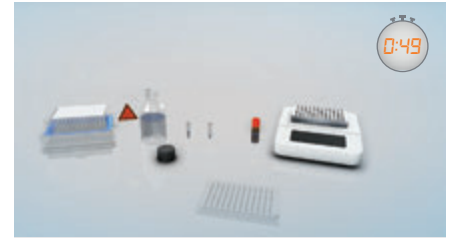
**17.** Add 400  $\mu$ L of Load/Wash solution into wells of the Gly-X Cleanup Plate for every sample in your Gly-X Deglycosylation Plate.  
Duration: 3 mins      Total time: 40 mins



**18.** Add 150  $\mu\text{L}$  of Load/Wash solution to the first sample well on the Gly-X Deglycosylation Plate. Mix well using the pipette.  
Duration: 3 mins      Total time: 43 mins



**19.** Transfer the entire sample ( $\sim 172 \mu\text{L}$ ) from the Gly-X Deglycosylation Plate into the corresponding well of the Gly-X Cleanup Plate. Mix well using the pipette. Repeat steps 18 and 19 until each sample has been transferred to the Gly-X Cleanup Plate.  
Duration: 3 mins      Total time: 46 mins



**20.** Apply vacuum ( $<5 \text{ in Hg}$ ) using a pump to draw the liquid through the well into the waste plate.  
Duration: 3 mins      Total time: 49 mins



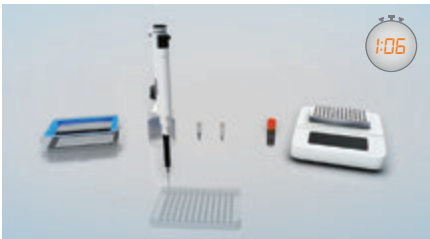
**21.** Add 600  $\mu\text{L}$  of Load/Wash Solution to each well. Apply 2 in Hg vacuum again, let the solution pass through. Repeat this step two more times for a total of three washes.  
Duration: 10 mins      Total time: 59 mins



**22.** Set the Gly-X Cleanup Plate aside. Install the Gly-X vacuum manifold spacer (black) and the Collection Plate (PCR plate).  
Duration: 1 min      Total time: 60 mins



**23.** Place the Gly-X Cleanup Plate back on the vacuum manifold. Add 100  $\mu\text{L}$  Gly-X InstantPC Eluent to each well of the Cleanup Plate containing sample. Apply  $<2 \text{ in Hg}$  vacuum.  
Duration: 3 mins      Total time: 63 mins



**24.** Take out the collection plate (PCR plate). Mix each sample prior to analysis. Cover plate with foil.  
Duration: 3 mins      Total time: 66 mins



**25.** Put the collection plate inside the Multi Sampler and run analysis.  
Duration: 1 min      Total time: 67 mins

## Complete Suite of Glycan Sample Preparation and Analysis Consumables

Everything you need is now available from Agilent:

- Fast, easy-to-use sample prep kits with a variety of N-glycan labels
- Labeled N-glycan standards and libraries
- Unlabeled N-glycan standards and libraries
- Endo- and exoglycosidases for structural characterization of N-glycans
- HILIC columns and other LC supplies for every level of analysis from intact glycoprotein to released glycans or monosaccharides

To learn more, visit [www.agilent.com/chem/glycananalysis](http://www.agilent.com/chem/glycananalysis)

[www.agilent.com/chem](http://www.agilent.com/chem)

DE44343.6400462963

**For Research Use Only. Not for use in diagnostic procedures**

This information is subject to change without notice.

© Agilent Technologies, Inc. 2021  
Printed in the USA, June 2, 2021  
5994-3482EN

