

Advancing Glycosciences. Together.

With the addition of ProZyme products and services, Agilent provides a single source offering for instruments and consumables, from sample to trusted answer. In addition to other critical quality attributes, our expertise now covers the complete glycan analysis workflow, so you can easily get the reliable, reproducible results you need.

Learn more: www.agilent.com/chem/better-together



- Key**
- Agilent Milestones
 - ProZyme Milestones
 - Agilent and ProZyme Collaborative Project

