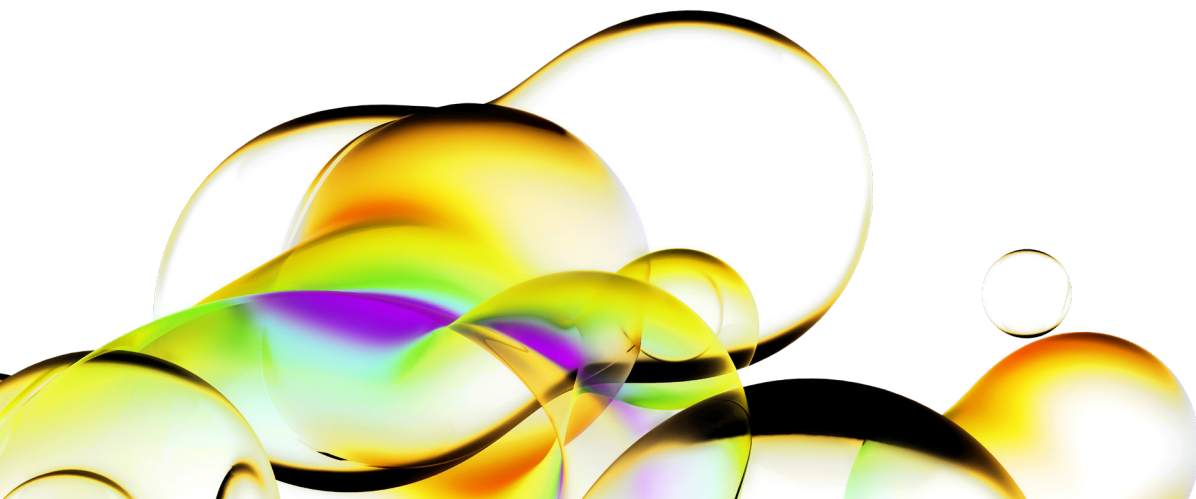




# Upgrade your LabChip instrument with Microsoft Windows 11.



A Windows 11 upgrade is now available for the LabChip GXII Touch protein Characterization System and LabChip GX Touch Nucleic Acid Analyzer.

Microsoft® Windows® 11 is the next evolution of Windows, which has been reimagined to boost productivity and enhance collaboration. The Windows 11 operating system is designed to enhance the performance, stability and security of your valuable IT investments.

It offers several improvements, enabling:

- **Emphasis on productivity:** Windows 11 provides a simple, powerful user experience that helps you focus and be more effective.
- **Streamlined user interface:** Layouts and personalized widgets to make it easier to stay organized.
- **Upgraded security features:** Windows 11 is the most secure version of Windows yet, ensuring your data and privacy is protected.



## We'll handle the Microsoft Windows 11 upgrade for your LabChip instrument

When you're ready to upgrade, we can help. A Revvity service engineer will come to your lab to install Microsoft Windows 11 on your instrument and configure the LabChip software saving your precious time for scientific work.

Included in the Microsoft Windows 11 software upgrade installation:

- Calibration of your instrument
- Transfer and testing of your current protocols

Because our service engineers are trained instrument maintenance experts, disruption to your scientific work will be minimized.

Allowing Revvity to upgrade your LabChip system allows for you to focus on value-added activities while we handle the upgrade process.

**To learn more, contact your Revvity representative today.**