



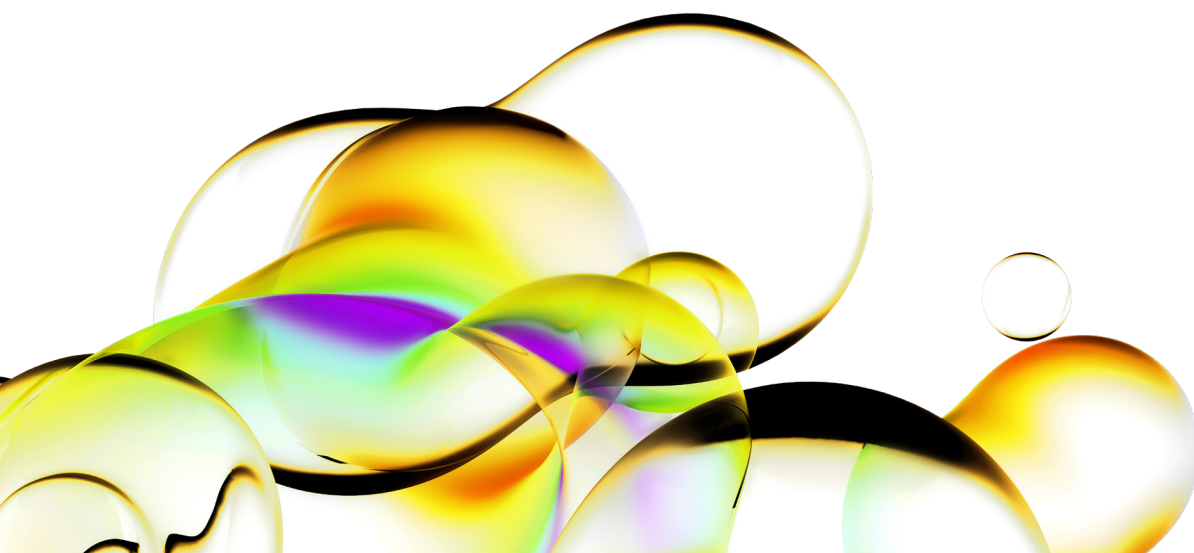
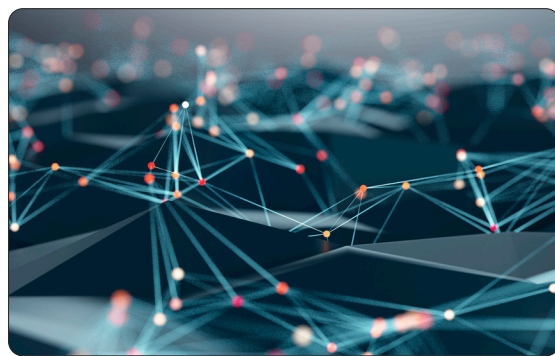
Empower your workflow with seamless integration.

LabChip Empower driver v1.0 for the LabChip systems

Take full control of your nucleic acid and protein assays on the LabChip GX Touch™ and LabChip GXII Touch™ systems with Waters Empower™ software. Seamlessly acquire and analyze your data with a single software with the new LabChip™ Empower™ v1.0 driver.

Key features:

- **Seamless Empower integration:** Manage LabChip assays from the Empower™ software's intuitive interface. Set up, start, and stop assays with ease, and monitor real-time assay status, chip info, and plate details—all from the Empower platform.
- **Optimized data management:** Automatically save raw data to the Empower Database (DB), eliminating manual data handling and ensuring secure, centralized data storage.
- **Advanced assay control:** Use the Method Editor to set up and run assays with precise control over assay type, plate, and sample selection.
- **Regulatory compliance:** Empower driver supports full audit trails, user access controls, and secure data storage, making it ideal for QA/QC labs in GxP-regulated environments.



Why upgrade to the LabChip Empower driver v1.0?

- **Improved efficiency:** By running LabChip systems directly from the Empower software, labs can streamline workflows, reduce hands-on time, and boost productivity.
- **Data integrity:** Ensure your data is securely stored and readily available for review, analysis, and audit through the Empower software's robust compliance features.
- **Future-proof:** LabChip Empower driver v1.0 is fully compatible with the Empower software, ensuring your system is ready for future software updates and advancements.

Technical requirements

- **Empower version:** 3.6.0 or above
- **Operating system:** Windows 11 (64-bit)
- **Software:** LabChip GX Touch SW 1.16 or above
- **LabChip instruments:**
Compatible with all LabChip Touch instruments

Upgrade and trade-in offers*

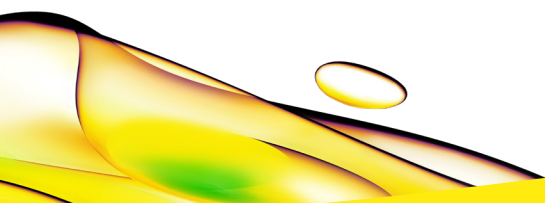
Upgrade package:

This package includes hardware and software upgrades to ensure full compatibility with the LabChip Empower driver v1.0.

Exclusive trade-in program:

Receive up to **\$20,000 USD** in trade-in credit when you upgrade older LabChip GX Touch systems. Plus, enjoy a **2-year extended service plan** at no extra cost when you trade in.

* Terms and conditions apply



revvity