

Advanced Sample Linking is an Agilent solution that simplifies basic workflows and helps to avoid error-prone steps.

The user-centric Sample Linking software is the core of the solution. It is accessed and used within the Bench Applications portal. It links the barcodes for samples and target vials, and communicates with Sample Scheduler for OpenLab to submit sequences.

A handheld scanner connected to a display (InfinityLab Assist Interface) serves as an interface for scanning barcoded containers and vials.

- Preparatory Steps
1. Prepare your system.

a. Ensure that your OpenLab CDS software, and Sample Scheduler for OpenLab have been updated to the current version.

b. Ensure the Sample ID Reader upgrade kit is available and installed to your Multisampler.

▶ See installation video or Agilent InfinityLab LC Series Multisampler Sample ID Upgrade Kit Installation Note (G4764-Sample-ID-Reader-Upgrade-TecPu-en-D0032639), provided on the USB media.
2. Install the Sample Linking software.

This prepares the software on the OpenLab Server

▶ See Advanced Sample Linking Installation and Configuration Guide (advanced-sample-linking-v1.0-installation-en.pdf).
3. Set up the Sample Linking hardware.

▶ Described in the Quick Start Guide provided with the kit (G4752-90000) and in the installation guide.
4. Log into the Bench Applications portal.

