

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 helps with running sample sequences error-free by connecting barcoded sample containers with lab data.

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 has been updated to the required 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 CDS WorkStation Plus

  - ▶ See Advanced Sample Linking Installation and Configuration Guide (advanced-sample-linking-v1.1-ws+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.

## Components of the Advanced Sample Linking Solution

### Sample Linking software M8270AA

#### Bench Applications Portal

Sample Linking software  
Acquisition Adapter  
Control Center  
Workflow Dashboard

#### Foundational software

OpenLab CDS Workstation Plus v2.8  
update 5 or later

### Sample Linking hardware G4752A

#### Sample Linking Interface G4752-64000

InfinityLab Assist Interface  
Display stand  
PoE Cable  
Quick Start Guide

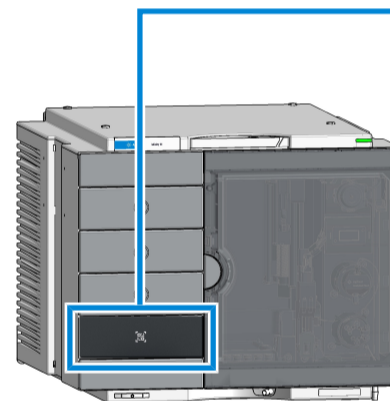


InfinityLab Assist Interface

Handheld scanner includes stand and cables

Display stand

### InfinityLab Sample ID Reader G4756A



Sample ID Reader  
Scanner module for  
Multisampler

Barcoded vials

40-Vial-Rack Sample ID

Palette Sample ID  
(Sample tray)

## Overview of the Advanced Sample Linking Solution Workflow

