


Control Panel

CONTROL PANEL

 Use the Desktop icon to launch OpenLab Control Panel.

MANAGE USERS

ROLE: SYSTEM ADMIN

- Assign users to groups.
- Assign roles to groups.
- Use default roles or assign privileges to roles (use role **Dissolution user** for Instrument Control and Data Analysis).
- Edit security policy.
- Manage licenses.

PROJECTS

ROLE: PROJECT ADMIN

Create projects to organize your data:

- Activate method audit trails.
- Apply project-specific options.
- Add custom parameters for samples and compounds.

★ Create shortcut for Data Analysis with a project.

 Launch Data Analysis for a project.

INSTRUMENTS

ROLE: INSTRUMENT ADMIN

Create and configure instruments
Monitor instruments

★ Create shortcut for acquisition with an instrument.

 Launch instrument control.

 Define Accessory Serial Numbers.

Dissolution Instrument Control

CREATE DISSOLUTION METHOD

Navigate to Method Layout

- Click **Method** (Method type/Apparatus type).
- Click **Create a new Method**.
- Define Method Type Manual (708-DS only) or 708-DS/850-DS.
- Define Apparatus Type.
- Define Sampling and instrument parameters for runs using this instrument configuration.


RUN A DISSOLUTION METHOD

- Enter sample parameters.
- Define Dissolution method.
- Define Dosage Introduction.
- Start **Run Options**.
- Number of active vessels.
- Click **Start Run**.

MANUAL CONTROL

Right click the instrument in the instrument status page to access manual control, cleaning cycle and to reset instrument.

STATUS

 View Activity Log, Method Run, Online Signals, Status and Timers..

Data Analysis

REVIEW DATA

- To load data, double click node in the **Data Selection** view.
- Upon double clicking a **Result Set**, the **DWM Method** and **DWM Run List** will be displayed.

PROCESSING METHOD

- Select the Dissolution Processing method to process the dissolution run data.
- This method can be edited via the edit processing method defined in Online Help.

REPORT AND PRINT

Dissolution specific parameters can be found under the DWM section of Reporting Items, these can be added to any custom report.



- Import default templates into your project** (start customizing templates from here):
Data Selection > Import Default Templates



- Print results manually:**
In Data Analysis, edit the processing method **Reports** section, and link the method. Click **Print all** in the ribbon.
Print processing method:
Go to **Data Processing > File** ribbon tab > **Print Method**



- Print Individual PDF:**
Go to **Reporting > Preview > Save as PDF**
Sequence summary report: submenu **Save result set summary report**.