

UNLEASH DISSOLUTION CAPABILITIES

AUTOSAMPLING WITH THE 850-DS: SAVING TIME AND REDUCING ERRORS

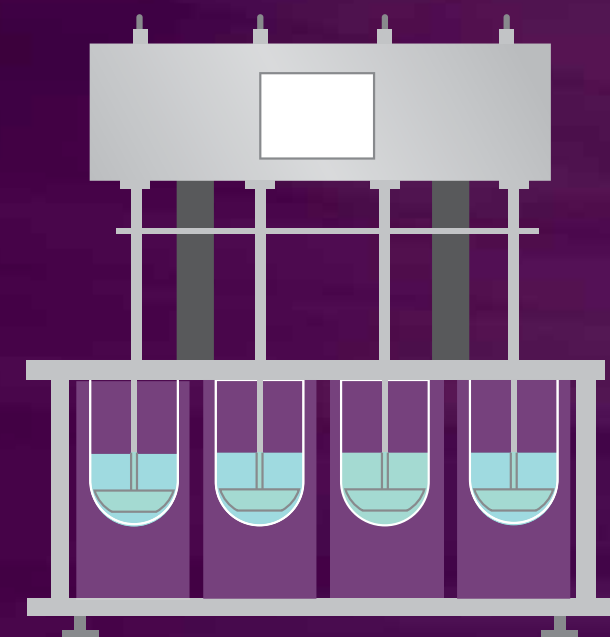
Whether you want to improve productivity or simply eliminate errors associated with manual sampling, the 850-DS Dissolution Sampling Station is the answer. The 850-DS controls your dissolution tests from start to finish. Enter and store your methods, initiate dissolution testing, monitor your results and perform a post-run cleaning, all from this easy-to-use autosampler. Learn more at www.agilent.com/lifesciences/850-DS.

The Measure of Confidence

01

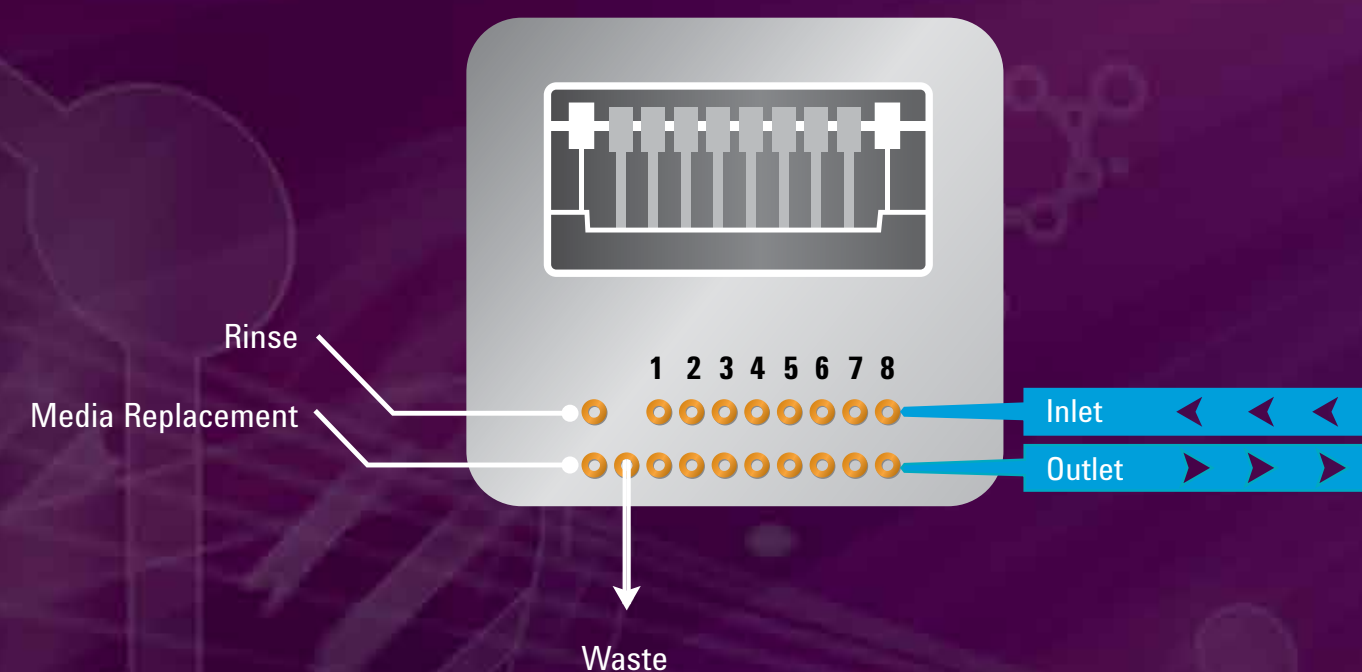
DISSOLUTION APPARATUS

Control USP Apparatus 1, 2, 3, 5, 6 and 7, including all Agilent, Varian and VanKel dissolution apparatus



Internal rotary piston pump for accurate, fast and precise liquid handling

- Media replacement to maintain vessel volume
- Handles common media types, including surfactants (up to 5% SLS)
- Custom cleaning cycle performed immediately following each test
- Sampling timepoints possible as often as every 2 minutes



02

850-DS DISSOLUTION SAMPLING STATION

Setup

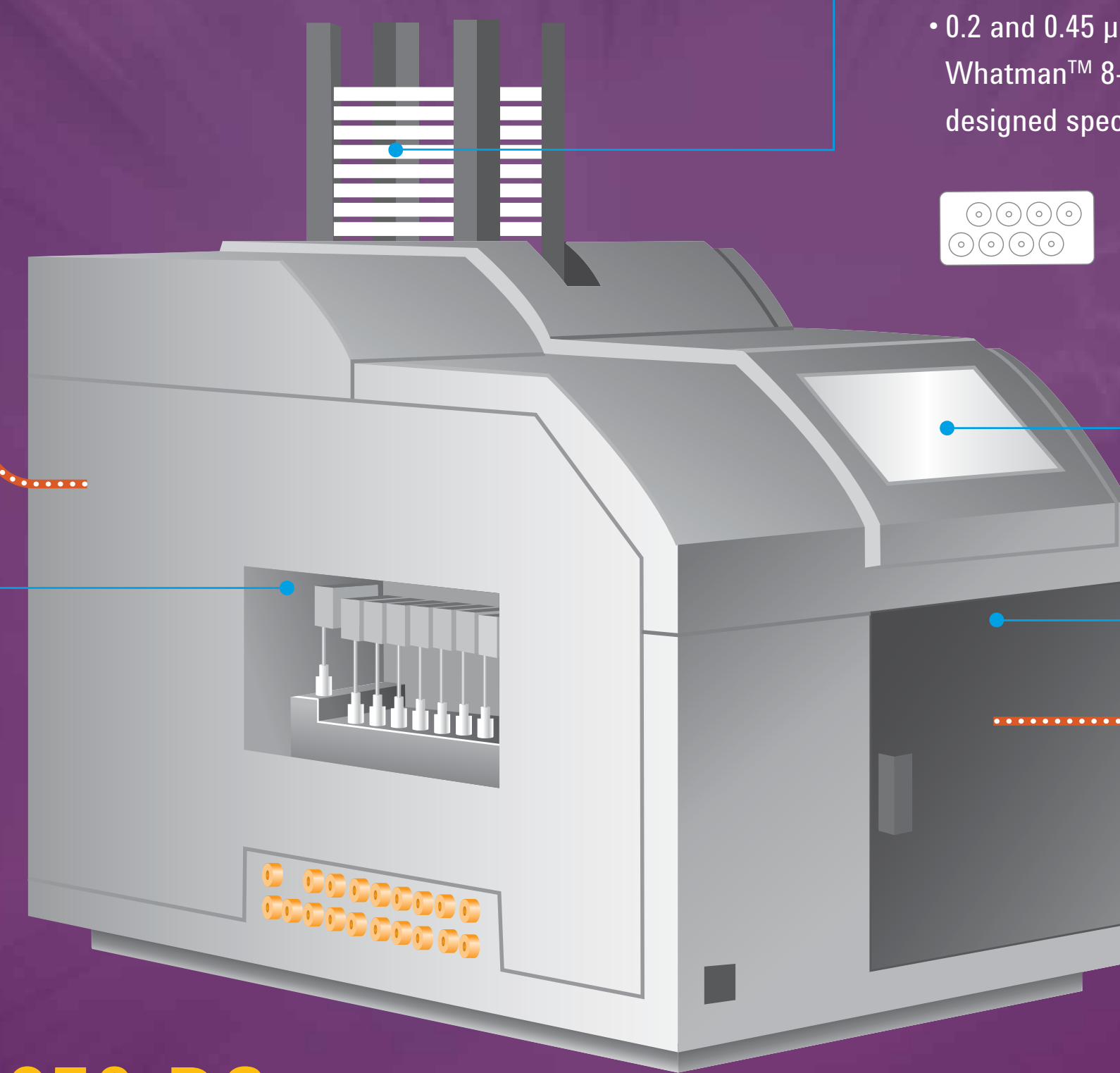
- Create and store methods for USP Apparatus 1, 2, 3, 5, 6 and 7
- Controlled access to secure dissolution methods and instrument settings
- Transfer methods to other 850-DS stations using SD card

Operation

- Repeatable, unattended sampling at the USP sampling location
- Eliminates analyst variability for repeatable results
- Link two 850-DS stations to two dissolution apparatus for synchronized starting and sampling

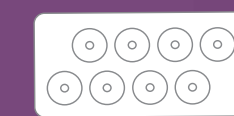
All-in-one Solution

- Complete 21 CFR Part 11 compliance with Dissolution Workstation Software
- Online UV-Vis analysis using the Cary 60 Spectrophotometer
- Automated cleaning cycle for long-term, reliable operation and a quick turnaround



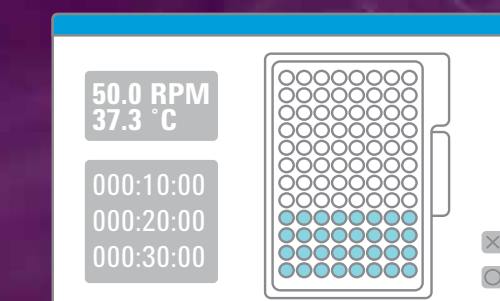
Filtration for a variety of methods

- Filtration can be easily enabled or disabled for any method
- Cannula tip and Full Flow filters ranging from 2 to 70 μm
- 0.2 and 0.45 μm membranes using the optional Whatman™ 8-channel filter plate from GE Healthcare, designed specifically for the 850-DS



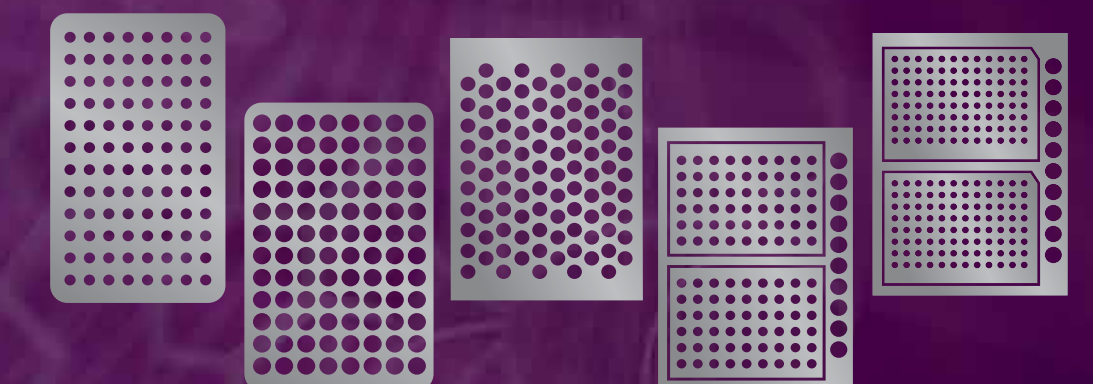
Intuitive user interface

- Data export options
- Synchronized start feature initiates two dissolution tests in parallel
- Supports 8 languages



Multiple sample tray options

- Supports test tubes, HPLC vials or multiple 96-well plates
- Agilent HPLC trays eliminate vial transfer, and can be placed directly in an Agilent LC



03

CARY 60 UV-VIS SPECTROPHOTOMETER

Samples can be analyzed online by the Cary 60, archived or collected for HPLC analysis

