

# The Agilent ICP Workflow Automation Solution

Less work. More flow.

## Reactive Dilution

If a sample result exceeds the calibration range or has an internal standard recovery issue, the ADS 2 can automatically perform an appropriate dilution and rerun the sample measurement.

## Prescriptive Dilution

Set a dilution factor for automatic sample preparation.

For instance, to dilute a sample 10 times, simply enter "10" for autodilution in the sample list.

## Optimized for Agilent ICP-OES/ICP-MS

The ADS 2 is optimized for Agilent ICPs, designed to work as an integrated system.

All settings are included in the method, so there's only one software application to learn.

Tight integration of the autodilutor with the instrument allows advanced features that can only be achieved when software and hardware are designed as one.

## Reduced Cost-Per-Analysis

The ADS 2 reduces operator time, labware usage (such as gloves, sample vials, and pipette tips), and waste disposal costs compared to manual dilutions.

## Automatic Standard Preparation

The ADS 2 will automatically calibrate your ICP.

Autocalibration assistant is used to define the calibration range. Calibration standards are then prepared by accurate auto-dilutions of a stock standard placed into the autosampler rack to generate multipoint calibration curves.

## Smart and Easy

Software tools streamline method development, data analysis, reporting, and troubleshooting when using the ADS 2.

Interactive flow path diagrams provide real-time system information about the ADS 2.

The Help and Learning Center includes detailed information on how to use and easily maintain the system.

## Faster Turnaround Times

When not actively diluting a solution, the ADS 2 has been optimized to minimize the time added to analysis.

The flow of solution is directed through the dilution flow path of the ADS 2 only when a dilution is triggered.



Learn more:  
<https://aglt.co/ro9k>

DE-005148

This information is subject to change without notice.

© Agilent Technologies, Inc. 2025  
Published in the USA, April 2, 2025  
5994-8243EN