

REIMAGINED. REDESIGNED.

AGILENT 8355 SULFUR CHEMILUMINESCENCE DETECTOR (SCD)–THE ONLY INTEGRATED SYSTEM FOR LOW-LEVEL SULFUR ANALYSIS

Simply put, we recognized that SCD technology must evolve to keep pace with growing demands for regulatory compliance and workflow efficiency.

We started by integrating the SCD with the Agilent 7890B—the world’s most reliable GC—to improve the overall user experience.. We then reengineered the burner assembly to maximize instrument uptime and simplify routine maintenance.

The result: the Agilent 8355 SCD—the only integrated system that delivers both selective and sensitive sulfur detection when time is critical and the right results are absolutely necessary.

Learn more about the *only* integrated system for low-level sulfur analysis. Visit www.agilent.com/chem/SCD

The Measure of Confidence



Meeting your needs for sensitivity, ASTM compliance, and ease of maintenance

Full integration

The Agilent 8355 SCD is now completely integrated into the Agilent 7890B GC with OpenLab CDS and MassHunter software—streamlining your lab’s analytical workflows and its ability to comply with regulatory requirements.

Worry-free maintenance, walk-up readiness

Maintaining the SCD burner, particularly replacing the inner ceramic tube, has never been easier. That’s because the simplified design of the 8355 SCD now allows you to replace the tube in just minutes, increasing workday productivity.

Robust and reliable

Electronic pneumatics control and digital electronics set a new benchmark for precision and repeatability—making the Agilent 8355 SCD the most sensitive and selective equimolar chemiluminescence detector that is linear in response.

Dual-plasma design

Eliminates quenching (signal suppression) caused by coelution of sulfur components and hydrocarbons.



Fully integrated Agilent 8355 SCD with Agilent 7890B GC and OpenLAB Chromatography Data System software
High selectivity and sensitivity for complex trace-sulfur methods



Work smarter with integrated GC, MSD, and software technologies



Standalone Agilent 8355 SCD
Compatible with legacy Agilent GC systems and other manufacturers’ GC systems

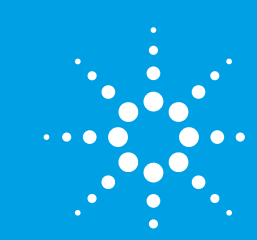


Standalone Agilent 8255 Nitrogen Chemiluminescence Detector For nitrogen detection



CP17989 Gas Clean Purification System (with Baseplate CP738407)
Offers fast, leak-free filter replacement

This information is subject to change without notice.
© Agilent Technologies, Inc., 2015
Printed in USA, September 24, 2015
5991-6294EN



Agilent Technologies