

# 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.

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

- Full integration
- Robust and reliable
- Worry-free maintenance, walk-up readiness
- Dual-plasma design

Learn more about the *only* integrated system for low-level sulfur analysis.  
Visit [www.agilent.com/chem/SCD](http://www.agilent.com/chem/SCD)



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



**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



**Work smarter** with  
integrated GC, MSD,  
and software technologies

The Measure of Confidence

 **Agilent Technologies**

This information is subject to change without notice.

© Agilent Technologies, Inc., 2015  
Printed in USA, September 24, 2015  
5991-6285EN