



COMPLEX CHARACTERIZATION MADE SIMPLE

The solutions and expertise you need to optimize your polymers

The polymer industry is made up of a diverse set of players, all facing similar challenges. Better characterization of materials and derivatives in the polymer value chain is vital to understand their thermal properties and suitability for use in a variety of applications. Faster product development, ensuring quality control, and reducing costs to stay ahead of the competition demands enhanced material characterization capabilities from a trusted partner. Our lineup of advanced solutions for thermogravimetric analysis (TGA) are designed to help you in:

- Identification, quantification, and characterization of polymers
- Ensuring quality control for raw materials and finished products
- New product development and improvement
- High-throughput and confident analyses
- Deeper insights with advanced hyphenation capabilities (FT-IR, GC/MS/MS)

Still not convinced? Here's why TGA is right for you.

Are you looking for complete sample atmosphere control, easy, unattended operation, and the ability to add hyphenation systems for maximum application flexibility? Look no further than the TGA 8000™ system with Pyris™ software.



TGA 8000

- Industry-leading atmosphere control
- 48-position autosampler for unattended use
- Wide temperature range (-20 °C to 1200 °C)
- Rapid heating (0.1 °C to 500 °C/min) and cooling for high throughput
- Fast, simple, seamless coupling for a complete hyphenation solution

Pyris Software

- Intuitive, easy-to-use interface that minimizes risk of errors
- Analysis of blends with exceptional separation of overlapping events and quantitative compositional analysis
- Allows real-time calculation during sample run
- Simplified data visualization at the touch of a screen
- Convenient monitoring capabilities from your iPhone® and iPad®

What consumables do I need for my polymer analyses?

You invest great efforts into your polymer testing – and we do the same with our consumables and accessories, tested and validated to fit your characterization needs. That's why we developed a full range of quality consumables such as sample pans, sample holders, and kits that work seamlessly with your Pyris software to tackle your daily challenges in the lab. Browse our consumables portfolio for products that offer reliable performance, reduce operating costs, and maximize uptime of your instruments. Like our trusted instruments, our consumables offer the best performance over and over.

Choose your TGA consumables and accessories here:

▶ [Thermal Analysis Consumables](#)

Trust the Consumables Engineered for Your Instruments

OneSource® Asset Management Services offers a single point of contact and custom service plans built to suit your needs. Some services include:

- **Improved Service Levels:** Proactive monitoring and management of service providers
- **Reduced Administration:** One contract and PO to manage, one contact to request service
- **Global Coverage:** 1,600 extensively trained and experienced team members located worldwide



Polymer Applications for Your Analysis

Characterization
of Polymers
Using TGA
Application Note

Polymer
Applications
Compendium

Plasticizer
Characterization
by TG-IR
Application Note

TGA 8000
Reference Guide