

SPECIFICATIONS OF THE PYRIS DSC 9



Introduction

The PerkinElmer Pyris™ DSC 9 marks the next generation in high temperature differential scanning calorimetry with the ability to reach temperatures of up to 750 °C. The DSC 9 provides a high-performance solution to a variety of solutions across a wide range of markets.

| Technical Specification | | Comments |
|--------------------------|---|---|
| Dynamic Range | +/- 175 mW | |
| Temperature Performance | | |
| Range | -100 to 750 °C | New furnace design allows for higher temperatures than achievable using most heat flux systems |
| Accuracy | +/- 0.05 °C | |
| Precision | < 0.02% ¹ | |
| Data Points/Sec | 10 | |
| Controlled Heating Rate | 0.1 to 100 °C/min | |
| Controlled Cooling Rate | 0.1 to 100 °C/min | |
| Calorimetric Performance | | |
| Calorimetric Accuracy | +/- 0.2% | |
| Calorimetric Precision | < 0.2% ¹ | |
| Indium Response Ratio | >19 mW/°C | Indium melting peak height/peak width. 7mg Indium, 1 °C/min. No mathematical correction for thermal resistance applied. |
| Indium Melting Time | 3.3 seconds | Time from the onset of the melting peak to the peak maximum |
| Purge gas | Nitrogen, Helium (Inert) Compressed Air, Oxygen (Oxidative) Argon (Inert) | |
| Weight | 12 kg | |
| Physical dimensions | 37x 34 x 32 cm (w x d x h) | |
| Instrument Control | Pyris Software control or stand-alone operation using the integrated touch screen | |

Data presented in this specification sheet were collected under ideal laboratory conditions and are used to demonstrate the specifications achievable by this system rather than specifications guaranteed in every analytical setting.

¹ Precision calculated as relative standard deviation.

PerkinElmer U.S. LLC
710 Bridgeport Ave.
Shelton, CT 06484-4794 USA
(+1) 855-726-9377
www.perkinelmer.com



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright ©2024, PerkinElmer U.S. LLC. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer U.S. LLC. All other trademarks are the property of their respective owners.