

# Thermo Scientific™ Partisol™ 2000i

## Air Sampler

### Maintenance and troubleshooting guide

<b>Operator</b>	
<b>Site location</b>	
<b>Instrument serial number</b>	
<b>Shelter temperature</b>	
<b>Software version</b>	

<b>Date</b>		
<b>Time</b>		

		Unit displayed value	Calibration date	Audit date	Measured results	Calibration offset/ intercept	Calibration Span	
<b>Instrument diagnostics</b>	Ambient temperature							
	Ambient pressure							
	Filter temperature							
	Filter pressure							
	Filter compartment temperature							
	Electronics compartment temperature							
	Ambient RH							
	Flow							
	Maximum pump pressure	atm						
	Internal leak check	mm Hg						
	External leak check	mm Hg						

Status code		Comments/notes
Code	Description	

For more information, please consult your technical support representative.

©2015 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.

**USA**  
 27 Forge Parkway  
 Franklin, MA 02038  
 Ph: (866) 282-0430  
 Fax: (508) 520-1460  
 customerservice.aqi@thermofisher.com

**India**  
 C/327, TTC Industrial Area  
 MIDC Pawane  
 New Mumbai 400 705, India  
 Ph: +91 22 4157 8800  
 india@thermofisher.com

**China**  
 +Units 702-715, 7th Floor  
 Tower West, Yonghe  
 Beijing, China 100007  
 +86 10 84193588  
 info.eid.china@thermofisher.com

**Europe**  
 Takkebijsters 1  
 Breda Netherlands 4801EB  
 +31 765795641  
 info.aq.breda@thermofisher.com

**Thermo**  
 SCIENTIFIC  
 A Thermo Fisher Scientific Brand