

Thermo Scientific™ Partisol™ 2000iD

Dichotomous Air Samplers

Maintenance and troubleshooting guide

Operator	
Site location	
Instrument serial number	
Shelter temperature	
Software version	

Date		
Time		

		Unit displayed value	Calibration date	Audit date	Measured results	Calibration offset/ intercept	Calibration Span
Instrument diagnostics	Ambient temperature						
	Ambient pressure						
	Filter temperature						
	Filter 2 temperature						
	Filter pressure						
	Filter 2 pressure						
	Filter compartment temperature						
	Electronics compartment temperature						
	Ambient RH						
	Flow						
	Flow 2						
	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

