

Thermo Scientific™ TEOM™ 1405

Ambient Particulate Monitor

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	
Instrument diagnostics	Ambient temperature					
	Ambient pressure					
					Measured	Difference
	Total flow					
	Bypass flow					
	Main flow					
	Leak check					
	Bypass flow					
	Zero					
	Actual					
	Difference					
	Main flow					
	Zero					
	Actual					
	Difference					
	Mass verification					
	Initial KO					
	Audit KO					
	% difference					
	F ₀ (initial frequency)					
F ₁ (final frequency)						
Pump pressure						

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

