

Thermo Scientific™ TEOM™ 1405DF

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			
					Base	Reference	Difference base	Difference reference
Instrument diagnostics	Ambient temperature							
	Ambient pressure							
	Total flow							
	Bypass flow							
	PM2.5 flow							
	Coarse flow							
	Leak check	Base	Reference					
	Bypass flow							
	Zero							
	Actual							
	Difference							
	PM2.5 flow							
	Zero							
	Actual							
	Difference							
	Coarse flow							
	Zero							
	Actual							
	Difference							
	Mass verification	PM2.5	PM coarse					
	Initial KO							
	Audit KO							
	% difference							
	F ₀ (initial frequency)							
	F ₁ (final frequency)							
Dryer dew point								
Chiller temperature								
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

Thermo
SCIENTIFIC
 Part of Thermo Fisher Scientific