

Thermo Scientific™ TEOM™ 1405F

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						
	Main flow						
	Leak check	Base	Reference				
	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)						
Dryer dew point							
Main 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