

Thermo Scientific™ TEOM™ 1405DF

Ambient Particulate Monitor

Instrument PM and verification sheet

Instrument serial number	
Instrument location	

Recommended preventative maintenance intervals				
Activity	PM interval	Last activity date	Current activity date	Next activity date
Replace TEOM filter(s):	Filter loading exceeds 75% or every 30 days			
Replace FDMS filter(s):	With every TEOM filter exchange			
Clean PM10 inlet:	With every TEOM filter exchange			
Clean virtual impactor:	With every TEOM filter exchange			
Replace in-line filters:	Every 6 months			
Clean coolers:	Annually			
Clean switching valve:	Annually			
Clean air inlet system:	Annually			
Rebuild vacuum pump:	Annually			
Dryer replacement:	Annually			
Replace mass transducer seals:	Annually			

Recommended instrument verification interval (unless regional requirements dictate otherwise)				
Equipment required:	A NIST-traceable (or equivalent) transfer standard for temperature calibrated for $\pm 0.5C$ precision			
	A NIST-traceable (or equivalent) transfer standard for pressure calibrated for ± 5 mm Hg precision			
	A NIST-traceable (or equivalent) flow transfer standard capable of measuring flow range of .9–19 lpm is required (depending on the flow transfer standard used, a flow inlet adapter may be required)			
Activity	Verification frequency	Last audit date	Current audit date	Next calibration date
Ambient temperature:	Audit monthly/calibration annually			
Ambient pressure:	Audit monthly/calibration annually			
Bypass flow:	Audit monthly/calibration annually			
PM2.5 flow:	Audit monthly/calibration annually			
Coarse flow:	Audit monthly/calibration annually			
Leak check:	Monthly (or if an audit dictates)			
Analog outputs:	Annually (or if an audit dictates)			
Mass verification:	Annually			

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.aq@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