

Thermo Scientific™ TEOM™ 1405

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			
Clean PM10 inlet:	With every TEOM filter exchange			
Clean VSCC inlet:	With every TEOM filter exchange			
Replace in-line filters:	Every six months			
Clean air inlet-system:	Annually			
Rebuild vacuum pump:	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 $\pm 0.5^{\circ}\text{C}$ 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 0.9-19 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/Calibrate annually			
Ambient pressure:	Audit monthly/Calibrate annually			
Bypass flow:	Audit monthly/Calibrate annually			
Main flow:	Audit monthly/Calibrate 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.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

A Thermo Fisher Scientific Brand