

INSTRUMENT PM AND VERIFICATION SHEET

Thermo Scientific™ TEOM™ 1405DF

Ambient particulate monitor

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 $\pm 0.5^\circ\text{C}$ 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 .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			

USA

27 Forge Parkway
Franklin, MA 02038
Ph: (508) 520-0430
Fax: (508) 520-2800
orders.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
Ph: +86 10 84193588
info.eid.china@thermofisher.com

Europe

Ion Path, Road Three,
Winsford, Cheshire CW73GA UK
Ph: +44 1606 548700
Fax: +44 1606 548711
sales.epm.uk@thermofisher.com

For more information, please consult your **technical support representative**

ThermoFisher
SCIENTIFIC