

Thermo Scientific™ Particulate Matter Continuous Emissions Monitoring System

INSTRUMENT PM AND VERIFICATION SHEET

Probe Controller serial #		Instrument Firmware Version	
Probe Monitor Serial #		Instrument location	

Recommended preventive maintenance intervals				
Activity	PM interval	Last activity date	Current activity date	Next activity date
Sample volume check	Daily			
Upscale drift (System check schedule)	Daily			
Zero drift check (System check schedule)	Daily			
Zero nephelometer	Daily			
Sample flow check with flow standard	Quarterly			
Bypass filter replacement	Quarterly			
Fan filter cleaning or replacement	Quarterly			
In-line filter monitor path	Quarterly			
In-line filter reference path	Quarterly			
TEOM filter replacement (may vary with higher concentrations)	Quarterly			
Replace Orings (dilution module, connection points in sample path)	Every 6 months			
External sample pump rebuild or replacement	Annually			
Sample line cleaning	Annually			
Nozzle & Dilution module cleaning	Annually			
Heated block exchange (Factory service only)	Annually			
Nephelometer assembly exchange (Factory service only)	Annually			

Recommended instrument verification interval (unless regional requirements dictate otherwise)				
Activity	Verification frequency	Last audit date	Current audit date	Next calibration date
Schedule Mass Reference	Every 24 hours @ peak load			
System check	Every 24 hours			
Blowback schedule	Every 24 hours/heavy loads every 8 hours			
Sample volume check	Every 24 hours			
Multi point span check	Every 24 hours to every 7 days			
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) 520-2800
orders.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
A Thermo Fisher Scientific Brand