

Thermo Scientific™ Partisol™ 2000iD

Dichotomous Air Samplers

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
Clean PM10 inlet:	Monthly			
Clean virtual impactor:	Monthly			
Replace in-line filters:	Every six months			
Clean filter cassettes/magazines:	Inspect regularly, clean as required			
Replace V-seals:	Annually			
Clean rain hood:	Every month			
Clean air screen/replace as necessary:	Every month			
Rebuild vacuum pump:	Inspect every six months, rebuild annually as a minimum			

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) transfer standard for relative humidity calibrated for $\pm 2.5\%$ 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 quarterly			
Filter temperature:	Audit monthly/Calibrate quarterly			
Filter compartment temperature:	Audit monthly/Calibrate annually			
Electronics compartment temperature:	Audit monthly/Calibration annually			
Ambient pressure:	Audit monthly/Calibrate quarterly			
Filter pressure:	Audit monthly/Calibrate quarterly			
Ambient RH:	Audit monthly/Calibrate annually			
Flow check:	Audit monthly/Calibrate quarterly			
Flow check 2:	Audit monthly/Calibrate quarterly			
External leak check:	Every five days of inlet usage			
Internal leak check:	Monthly			

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