

# Thermo Scientific™ 5014i Beta

## Continuous Particulate Monitor

Instrument PM and verification sheet

|                                 |  |
|---------------------------------|--|
| <b>Instrument serial number</b> |  |
| <b>Instrument location</b>      |  |

### Recommended preventative maintenance intervals

| Activity          | PM interval | Last activity date | Current activity date | Next activity date |
|-------------------|-------------|--------------------|-----------------------|--------------------|
| Clean PM10 inlet: | Monthly     |                    |                       |                    |
| Clean VSCC inlet: | Monthly     |                    |                       |                    |

### Recommended instrument verification interval (unless regional requirements dictate otherwise)

|                            |   |
|----------------------------|---|
| <b>Equipment required:</b> | 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 $\pm 5$ 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 .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 RH:               | Audit monthly/Calibrate annually |                 |                    |                       |
| Ambient pressure:         | Audit monthly/Calibrate annually |                 |                    |                       |
| Vacuum flow:              | Calibrate annually               |                 |                    |                       |
| Vacuum pressure:          | Calibrate annually               |                 |                    |                       |
| Flow check:               | Audit monthly/Calibrate annually |                 |                    |                       |
| Leak check:               | Audit monthly                    |                 |                    |                       |
| Integrity of filter spot: | Audit monthly                    |                 |                    |                       |
| Detector calibration:     | Annually                         |                 |                    |                       |
| Mass calibration:         | 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.epm@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**  
Takkelijsters 1  
Breda Netherlands 4801EB  
+31 765795641  
info.aq.breda@thermofisher.com

**Thermo**  
SCIENTIFIC

A Thermo Fisher Scientific Brand