

# Thermo Scientific™ 5014i Beta

## Continuous Particulate Monitor

### Maintenance and troubleshooting guide

<b>Operator</b>	
<b>Site location</b>	
<b>Instrument serial number</b>	
<b>Shelter temperature</b>	
<b>Software version</b>	

<b>Date</b>	
<b>Time</b>	

		Unit displayed value	Calibration date	Audit date	Measured results
<b>Instrument diagnostics</b>	Ambient temperature				
	Ambient pressure				
	Ambient RH				
	Flow				
		<b>Unrestricted</b>	<b>Restricted</b>		
	Leak Check				
	L Recs	Download last 48 hours minimum. Switch unit to collect 1-minute data for 24 hours minimum.			

Status/Alarm code		Comments/Notes
Code	Description	

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