

# Thermo Scientific™ TEOM™ 1405DF

## Ambient particulate monitor

Operator	
Site location	
Instrument serial number	
Shelter temperature	
Software version	

Date			
Time			

		Unit displayed value	Calibration date	Audit date	Measured results			
Instrument diagnostics	Ambient temperature							
	Ambient pressure							
					Based	Reference	Difference base	Difference reference
	Total flow							
	Bypass flow							
	PM2.5 flow							
	Coarse flow							
	Leak check	Base	Reference					
	Bypass flow							
	Zero							
	Actual							
	Difference							
	PM2.5 flow							
	Zero							
	Actual							
	Difference							
	Coarse flow							
	Zero							
	Actual							
	Difference							
	Mass verification	Base	Reference					
	Initial K0							
	Audit K0							
	% difference							
	F <sub>0</sub> (initial frequency)							
	F <sub>1</sub> (final frequency)							
	Dryer dew point							
	Chiller temperature							
	Pump pressure							

Status code		Comments/notes
Code	Description	

### 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