

# Thermo Scientific™ PDM3700 Personal Dust Monitor

## Spare Parts Recommendations

| Part Number                   | Description   | Recommended to have on-site for regular maintenance |
|-------------------------------|---|---|
| <b>PDM Parts</b>              |   |   |
| 56-010897-2300                | Battery assembly, PDM3700   |   |
| 56-009961                     | Flow path seal kit, PDM3700 (3 O-rings, 1 cyclone gasket, 1 inlet plate gasket) | X   |
| 30-009596                     | Cyclone grit pot  | X   |
| <b>Charger Parts</b>          |   |   |
| 55-010924-0001                | Interface assembly, charger, PDM3700  |   |
| 07-000587                     | Cable, RS232 m/f 10 ft  |   |
| 04-003268                     | Fuse, 2A time delay   | X   |
| 07-000593                     | Cord, power, detachable (for charger input power)                               |   |
| 57-010925                     | Charger assembly, PDM3700   |   |
| 07-010239                     | Cable, USB to serial, 6 ft  |   |
| <b>Accessories</b>            |   |   |
| 59-010234-0005                | Overlay, Display peel-away (pkg of 5)   | X   |
| 57-009727-0020                | Filter, PDM3700 box of 20   | X   |
| 57-009727-1000                | Filter, PDM3700 box of 1000   | X   |
| 34-009958                     | Rugged carry case   |   |
| <b>Tools</b>                  |   |   |
| 55-009885                     | Filter tool assembly, PDM3700   | X   |
| 30-009908                     | Tool, hex key 7/64" ball end  |   |
| 59-010020                     | KO audit kit  |   |
| 59-010019                     | Flow calibration kit (includes orifice set)                                     |   |
| 59-010237                     | Leak check kit (suction cup, adapter, inlet plug)                               |   |
| <b>Samplet Inlet Assembly</b> |   |   |
| 55-012085                     | Inlet/tubing assembly   | X   |
| 37-012084                     | Inlet   |   |
| 38-012094                     | Inlet mounting bracket  |   |
| 35-012093                     | Sealing screw   |   |
| 22-000485-1012                | Inlet o-ring  | X   |
| 21-003721-0003                | Inlet screws  | X   |
| 55-012278                     | Inlet bracket clip attachment   |   |
| 32-009973-0001                | Tubing, pre-cut   | X   |
| 30-012091                     | Tubing clip   | X   |
| 30-012264                     | Inlet tool  |   |

For more information, visit [www.thermoscientific.com/endblacklung](http://www.thermoscientific.com/endblacklung)

© 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. This product is manufactured in a plant whose quality management system is ISO 9001 certified.

**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