

Thermo Scientific™ 1405-DF TEOM™, Continuous Dichotomous Ambient Air Monitor

Recommended Spare Parts Kits

To keep your system running and properly collecting data, it's important to have spare parts on hand to avoid costly downtime. We developed Spare Parts Kits to ensure maximum uptime of your Thermo Scientific Instrument during its life cycle.

| Consumables Parts Kit # 116270-00 | | |
|-----------------------------------|---------------------------------|----------|
| Part Numbers | Description | Quantity |
| 59-010993 | 1405DF Consumables PKG | 1 |
| 22-009863 | TEOM V-Seal | 2 |
| 56-009872-EXCH | Dryer Refurb Program | 2 |
| 22-000485-1036 | PM10 Inlet Lg O-Ring | 1 |
| 22-002853-3026 | PM10 Inlet Sm O-Ring | 2 |
| 109773-00 | Virtual Impactor O-Ring Kit | 1 |
| 15-011185-00DF | Compact flash card with program | 1 |
| 111754-00 | Pump Re-build Kit | 1 |

| Extended Maintenance Parts Kit # 116271-00 | | |
|--|-----------------------------|----------|
| Part Numbers | Description | Quantity |
| 103674-00 | Fuse in corcom | 2 |
| 55-010022 | Mass Flow Controller Dichot | 1 |
| 56-011934 | Switching Valve Dichot | 1 |
| 110836-00 | Vacuum Pump | 1 |
| 50-011081 | Interface PCB | 1 |
| | | |
| | | |

Please have PO made out to THERMO ENVIRONMENTAL INSTRUMENTS, INC.

© 2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo 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.

Contact Customer Service to Order Today

USA
27 Forge Parkway
Franklin, MA 02038
Ph: (866) 282-0430
Fax: (508) 520-2800
orders.aqi@thermofisher.com