

Thermo Scientific™ Particulate Matter Continuous Emissions Monitoring System (CEMS)

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.

POS Parts Kit # 116606-00		
Part Numbers	Description	Quantity
57-009727-0020	Box of 20 TEOM Filters (PDM 3600)	1
55-012013	Filter Assembly Jumbo HEPA	1
113552-00	Pump Re-build Kit (bypass)	1
59-012051	O-Ring Kit	1
55-012005	Filter Assembly Medium In-line	1
113547-00	Sample Check Valve Assembly	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