

Thermo Scientific™ Model 49i-PS Ozone Primary Standard Analyzer

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

Consumables Parts Kit # 115896-00		
Part Numbers	Description	Quantity
4510	Fuse, 3.0 Amp, S/B, 110V	2
108002-00	Pump Repair Kit	1
4124	Capillary, 0.015-Inch ID (Purple)	2
4111	Capillary, 0.028-Inch ID (Grey)	1
7336	Capillary, 0.018-Inch ID (Purple/Blue)	1
4800	Capillary O Ring (10 Per Pack)	1

Extended Maintenance Parts Kit # 115887-00		
Part Numbers	Description	Quantity
100907-00	Fan, 24VDC, I-Series	1
101023-00	Pressure Transducer	1
101681-00	Power Supply W/O Cables	1
102055-00	Flow Transducer	1
102441-02	Sample / Reference Solenoid Valve Assembly	1
111553-01	Pump Assembly	1
8540	Photometer Lamp	1
8542	Absorption Tube Cell	2
8592	Detector Assembly	1
8645	Ozonator Lamp	1

© 2015 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.aqj@thermofisher.com