

Thermo Scientific™ Model 49i Ozone 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 uptime of your Thermo Scientific Instrument during its life cycle.

Consumables Parts Kit # 112457-00			Extended Maintenance Parts Kit # 112848-00		
Part Numbers	Description	Quantity	Part Numbers	Description	Quantity
4124	Capillary, 15 Mil Short (Purple)	2	100907-00	Fan, 24VDC, I-Series	1
14697	Scrubber Assembly	1	101023-00	Pressure Transducer	1
108002-00	Pump Repair Kit	1	101681-00	Power Supply W/O Cables	1
4510	Fuse 3 Amp ,S/B, 110V	2	102055-00	Flow Transducer	1
11290	Receptacle Fuse Drawer, 110V (Add)	1	102441-01	Sample / Reference Solenoid Valve Assembly	1
			111553-01	Pump Assembly	1
			8540	Photometer Lamp	1
			8542	Absorption Tube Cell	2
			8592	Detector Assembly	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