

Thermo Scientific™ Model 51i-HT Total Hydrocarbon Analyzer High Temperature 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 # 115898-00		
Part Numbers	Description	Quantity
4070	Fan Filter	1
11290	Fuse Drawer	1
12190	Filter Inlet Element, 0.5 micron	1
12188	Aluminum Gasket (For P/N 12190)	1
11032	O-Ring, FID Flame Sensor (2/Pk)	1
105198-00	Fuse, 250 VAC, 6.25 Amp	2
105567-00	Ignitor/Cable Assembly	1
108091-00	Pump Repair Kit-(51i-HT)	1
103581-00	Flame Sensor Assembly	1
104139-00	Sample Frit, FID	1

Extended Maintenance Parts Kit # 115891-00		
Part Numbers	Description	Quantity
103803-00	Fan 4", 24v	1
112286-00	Sample Pump (HT)	1
101681-00	Power supply 24V	1
104151-00	Exhaust Frit Bulkhead Fitting	1
105554-01	Bypass Orifice Bulkhead (51iHT)	1
103950-00	Air Frit, FID	1
103687-00	Pogo Detector Assembly	1
103762-00	Solenoid Valve Assembly	1
104872-00	Pressure Regulator	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