

# Thermo Scientific™ Model 146i Multi-Gas Calibrator

## Spare Parts Recommendations

Part Number	Description	Annual Maintenance	Age of Instrument or Fleet (Years)					
			1	2	3	4	5	
8606	PUMP REPAIR KIT, KNF (Original Style)	X	X					
108002-00	Pump Repair Kit (new technology pump)	X	X					
101426-00	Pump, KNF 110VAC with Plate and Fittings (Optional)			X				
103206-00	Solenoid Valve, Span Gas	X		X				
100880-00	Measurement Interface Board (146i)				X			
100533-00	Mother Board				X			
101023-00	Pressure Transducer Assembly				X			
103125-00	Solenoid Valve, External/internal				X			
103200-00	Solenoid Valve, Perm 2				X			
103207-00	Solenoid Valve, Reaction Chamber				X			
103209-00	Detector Assembly (Option)				X			
103334-00	Solenoid Valve, Perm 1				X			
103075-00	Back Pressure Regulator							X
100907-00	Fan, 24vdc						X	
101681-00	Power Supply, 24vdc						X	
102351-00	Photometer Lamp Heater Assembly						X	
100480-00	Keypad Membrane Assembly							X
100539-00	Digital Output Board							X
100542-00	I/O Expansion Board (optional)							X
100895-00	Ozonator Powersupply Board							X
101055-00	Assembly, Power Receptical							X
101491-17	Processor Board, Model 146i							X
102496-00	Front Panel Display							X
102554-00	Photometer Board							X
110608-00	Ozonator Assembly							X
102604-00	Pressure Regulator							X
8630	Fan Filter / Guard Assembly	X						
4121	Capillary 10 Mil Black (If GPT Option is Installed)	X						
8540	Photometer Lamp (Optional)	X						
8645	Ozonator Lamp (If GPT Option is Installed)	X						
4510	Fuse, 3 AMP S/B (pk/5)	X						

© 2014 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher 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.

**USA**  
27 Forge Parkway  
Franklin, MA 02038  
Ph: (866) 282-0430  
Fax: (508) 520-1460  
customerservice.aq@thermofisher.com

**India**  
C/327, TTC Industrial Area  
MIDC Pawane  
New Mumbai 400 705, India  
Ph: +91 22 4157 8800  
india@thermofisher.com

**China**  
+Units 702-715, 7th Floor  
Tower West, Yonghe  
Beijing, China 100007  
+86 10 84193588  
info.eid.china@thermofisher.com

**Europe**  
Takkebijsters 1  
Breda Netherlands 4801EB  
+31 765795641  
info.aq.breda@thermofisher.com