

Thermo Scientific™ Model 15C Analyzer

Unique C-Series Spare Parts

Part Number	Description
8331	PCB A/D (48C)
8935	PCB TEMP. CONTROLLER (48C)
9952	PCB PUSH BUTTON (B2)
9963	MOTHER TO DISPLAY CABLE

© 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.aqj@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

Thermo Scientific™ Model 17C Analyzer

Unique C-Series Spare Parts

Part Number	Description
10223	PCB TEMPERATURE CONTROL (CONVERTER MODULE 17C)
9841	PCB ANALOG/DIGITAL
9952	PCB PUSH BUTTON (B2)
9963	MOTHER TO DISPLAY CABLE
9981	BASE SOCKET ("C" SERIES)

© 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

Thermo Scientific™ Model 41C Analyzer

Unique C-Series Spare Parts

Part Number	Description	41C	41C-HL
8331	PCB A/D (48C)	X	X
8935	PCB TEMP. CONTROLLER (48C)	X	X
9952	PCB PUSH BUTTON (B2)	X	X
9963	MOTHER TO DISPLAY CABLE	X	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

Thermo Scientific™ Model 42C Analyzer

Unique C-Series Spare Parts

Part Number	Description	42C	42C-HL	42C-LS	42C-TL	42C-Y	42C-D
10196	SCRUBBER ASSEMBLY, NH3 (42CLS)			X			
10284	ASSY S/S/CONVERTER (42CHL-220)	X	X	X			X
10765	PCB TEMP. CONTROLLER (MODEL 42CHL)		X	X			
10766	PCB INPUT (42CHL)		X	X			
10768	PCB POWER SUPPLY ASSY				X		
12522	FLOW SWITCH ASSEMBLY (42CHL) (STANDARD ONLY)		X				
64P7157-1	ASSY, 120V COOLER PARTS (42C, 42C-D)	X					X
8119-02	VALVE SOLENOID (146C)	X					
9841	PCB ANALOG/DIGITAL	X	X	X	X	X	X
9952	PCB PUSH BUTTON (B2)	X	X		X	X	X
9963	MOTHER TO DISPLAY CABLE	X					
9981	BASE SOCKET ("C" SERIES)	X	X	X	X		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

Thermo Scientific™ Model 43C Analyzer

Unique C-Series Spare Parts

Part Number	Description	43C	43C-HL	43C-TLE
10996	PCB FLASHER SUPPLY			X
57P7107-4	43C POWER SUPPLY, ASSY	X		
57P713-1	OPTICAL BENCH ASSY, 120V	X		
8119-02	VALVE SOLENOID (146C)	X		
8943	PCB A / D	X	X	X
8951	PCB INPUT	X	X	X
9681	PCB FLASHER SUPPLY (43B-10)	X	X	
9829	PCB MOTHER	X	X	X
9952	PCB PUSH BUTTON (B2)	X	X	X
9963	MOTHER TO DISPLAY CABLE			X
99976	INPUT BD CABLE, 43C	X	X	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

Thermo Scientific™ Model 46C Analyzer

Unique C-Series Spare Parts

Part Number	Description	46C	46C-HL
8331	PCB A/D (48C)	X	X
8935	PCB TEMP. CONTROLLER (48C)	X	X
9952	PCB PUSH BUTTON (B2)	X	X
9963	MOTHER TO DISPLAY CABLE	X	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

Thermo Scientific™ Model 48C Analyzer

Unique C-Series Spare Parts

Part Number	Description	48C	48C-HL	48C-TL	48C-TLE
8331	PCB A/D (48C)	X	X	X	X
8935	PCB TEMP. CONTROLLER (48C)	X	X	X	X
9952	PCB PUSH BUTTON (B2)	X	X	X	X
9963	MOTHER TO DISPLAY CABLE		X		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

Thermo Scientific™ Model 49C Analyzer

Unique C-Series Spare Parts

Part Number	Description	49C	49C-PS
10758	PCB LAMP POWER SUPPLY (49C)	X	X
10760	PCB OZONATOR	X	X
10761	PCB ANALOG/DIGITAL (49C)	X	X
13405	SOFTWARE, C-LINK (49C)	X	
45P752	ASSY OPTICAL BENCH 49C	X	X
8314	THERMISTOR ASSEMBLY (49)	X	X
9847	PCB POWER SUPPLY (49C)	X	X
9952	PCB PUSH BUTTON (B2)	X	X
9963	MOTHER TO DISPLAY CABLE	X	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

Thermo Scientific™ Model 51C Analyzer

Unique C-Series Spare Parts

Part Number	Description	51C-HT	51C-LT
112326-01	PUMP ASSY (51CHT) 110V	X	
14951	SOLENOID VALVE 2-WAY (55C)	X	X
16106	PCB ASSY, MOTHER BOARD	X	X
16107	PCB ASSY, A/D	X	X
81P729	THERMOCOUPLE ASSEMBLY	X	X
81P739	SIGNAL POGO ASSY	X	X
81P751	BIAS POGO ASSY	X	X
81P758	ZERO SPAN FRIT ASSEMBLY	X	X
81P759	DETECTOR REBUILD KIT (51C)	X	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

Thermo Scientific™ Model 55C Analyzer

Unique C-Series Spare Parts

Part Number	Description
10774-01	ELECTROMETER BD, LOW RANGE
10774-02	ELECTROMETER BD, MID RANGE
10774-06	ELECTROMETER BD, CHRYSLER SPECIAL RANGE
109315-00	PUMP VACUUM 115/60
109316-00	PUMP VACUUM 220/50
13258	PCB POWER SUPPLY (110V)
14695	FERRULES 1/8" (COLUMN)
14951	SOLENOID VALVE 2-WAY (55C)
14952	SOLENOID VALVE 3-WAY (55C)
78P775	SIGNAL POST ASSEMBLY
9952	PCB PUSH BUTTON (B2)
9963	MOTHER TO DISPLAY CABLE

© 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

Thermo Scientific™ Model 146C Analyzer

Unique C-Series Spare Parts

Part Number	Description
10760	PCB OZONATOR
10761	PCB ANALOG/DIGITAL (49C)
14291	SPARE PARTS 146C POWER SUPPLY
8119-02	VALVE SOLENOID (146C)
9952	PCB PUSH BUTTON (B2)
9963	MOTHER TO DISPLAY CABLE

© 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

Thermo Scientific™ Model 450C Analyzer

Unique C-Series Spare Parts

Part Number	Description	450C	450C-TL
450P702-1	CONVERTER ASSEMBLY (450C + 450CTL) 110V	X	X
450P702-2	CONVERTER ASSEMBLY (450C + 450CTL) 220V	X	X
9963	MOTHER TO DISPLAY CABLE	X	X
99976	INPUT BD CABLE, 43C	X	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