

TriPlus RSH: Spare Parts Guide

This guide contains part numbers for all of the replaceable components in the TriPlus RSH. Refer to the *TriPlus RSH Hardware Manual* for information about installing these components in your instrument. Throughout the TriPlus RSH documentation, any component with a part number can be ordered from us. Components without a part number are not available.

IMPORTANT Only components with a part number are available for purchase. The part you are looking for may be shown in an illustration, but if it does not list a part number, it is not available.

You can purchase the following column components for the TriPlus RSH autosampler. Be sure to reference the component's part number when placing an order with your local Sales/Service Representative. See the *TriPlus RSH Hardware Manual* for information about installing these components or [Table 1](#) for a comprehensive list of available component.

Table 1. Parts That Can Be Reordered (Sheet 1 of 3)

Component	Thermo Scientific Parts Number
TriPlus RSH - TRACE Ultra Mounting Kit	1R77010-1001
TriPlus RSH - TRACE Ultra Injector Adapter Kit	1R77010-1002
TriPlus RSH - FOCUS GC Mounting Kit	1R77010-1003
TriPlus RSH - FOCUS GC Single Injector Adapter Kit	1R77010-1004
TriPlus RSH - TRACE 1300/1310 Mounting Kit	1R77010-1005
TriPlus RSH - TRACE 1300/1310 Injector Adapter Kit	1R77010-1006
TriPlus RSH - Liquid Tool: D7, 57 mm	1R77010-1007
TriPlus RSH - Liquid Tool: D7, 85mm	1R77010-1008
TriPlus RSH - Liquid Tool: D8, 57mm	1R77010-1009
TriPlus RSH - Liquid Tool: D8, 85 mm	1R77010-1010
TriPlus RSH - Liquid Tool: D18, 57 mm	1R77010-1011
TriPlus RSH - SPME Syringe Fiber Tool Kit for USA only.	1R77010-1099
TriPlus RSH - SPME Syringe Fiber Tool Kit for World Wide Distribution except USA.	1R77010-1100
TriPlus RSH - HS Tool: 1000 µL	1R77010-1012
TriPlus RSH - HS Tool: 2500 µL	1R77010-1013
TriPlus RSH - HS Tool: 5000 µL	1R77010-1014
TriPlus RSH - Automatic Tool Change Station	1R77010-1019
TriPlus RSH - Tool Releasing Station	1R77010-1020

Table 1. Parts That Can Be Reordered (Sheet 2 of 3)

Component	Thermo Scientific Parts Number
TriPlus RSH - Standard Tray Holder	1R77010-1021
TriPlus RSH - VT15, 15-vial Tray: 10/20 mL	1R77010-1022
TriPlus RSH - VT54, 54-vial Tray: 2 mL	1R77010-1023
TriPlus RSH - VT70, 70-vial Tray: 1mL	1R77010-1024
TriPlus RSH - R60, 60-vial Tray: 10/20mL	1R77010-1025
TriPlus RSH - Liquid Cooled Tray Holder Kit	1R77010-1026
TriPlus RSH - Liquid Cooled 32-Vial Tray: 10/20 mL	1R77010-1027
TriPlus RSH - Temperature Controlled Drawer: Single	1R77010-1028
TriPlus RSH - Standard Wash Station: 5x10 mL Vials	1R77010-1029
TriPlus RSH - Large Wash Station: 2x100 mL Vials	1R77010-1030
TriPlus RSH - Solvent Station: 3x100 mL Vials	1R77010-1031
TriPlus RSH - Fast Washing Module for Two Solvents Kit	1R77010-1098
TriPlus RSH - Agitator/Incubator	1R77010-1032
TriPlus RSH - Vortexer	1R77010-1033
TriPlus RSH - 1D Barcode Reader	1R77010-1034
TriPlus RSH - SPME Fiber Conditioning Station	1R77010-1035
TriPlus RSH - Multiple HS Extraction Module	1R77010-1036
TriPlus RSH - Actuator for TRACE GC Ultra OCI	1R77010-1037
TriPlus RSH - Hand-held Terminal	1R77010-1038
TriPlus RSH - Headspace Upgrade Kit	1R77010-1039
TriPlus RSH - Preventive Maintenance Kit	1R77010-1040
TriPlus RSH - Gas Tube Replacement Kit	1R77010-1041
TriPlus RSH - Z-Head for Exchange	1R77010-1042
TriPlus RSH - Z-Axis Motor Assy Kit	1R77010-1045
TriPlus RSH - Needle Guide Motor Assy Kit	1R77010-1046
TriPlus RSH - Syringe Plunger Motor Assy Kit	1R77010-1047
TriPlus RSH - Z-Head PCB Kit	1R77010-1053
TriPlus RSH - Agitator PCB Kit	1R77010-1054
TriPlus RSH - ATC Station PCB Kit	1R77010-1055
TriPlus RSH - Temperature Controller Drawer PCB Kit	1R77010-1056
TriPlus RSH - Main Power Supply	1R77010-1057
TriPlus RSH - 100 mL Solvent Vial for Washing	1R77010-1060
TriPlus RSH - Pe Snap Caps and Seals (Set of 10)	1R77010-1061

Table 1. Parts That Can Be Reordered (Sheet 3 of 3)

Component	Thermo Scientific Parts Number
TriPlus RSH - 10 mL Solvent Vial for Washing (Set of 10)	1R77010-1062
TriPlus RSH - TRACE GC Ultra Mounting Left Leg	1R77010-1063
TriPlus RSH - TRACE GC Ultra Mounting Right Leg	1R77010-1064
TriPlus RSH - FOCUS GC Mounting Leg	1R77010-1065
TriPlus RSH - TRACE 1300/1310 Mounting Leg	1R77010-1066
TriPlus RSH - Agitator Adapter for 2 mL Vial (Set of 6)	1R77010-1067
TriPlus RSH - Agitator Adapter for 10mL Vial (Set of 6)	1R77010-1068
TriPlus RSH - Standard Washing Station Adapter for 2 mL Vial	1R77010-1069
TriPlus RSH - Foil Cutter for Microtiter and Deep Well Plates	1R77010-1070
TriPlus RSH - Main PCB Kit	1R77010-1073
The kit includes: TMO Key GC1, TMO CPU, Right Side Cover including LED Light Pipe	
TriPlus RSH - XY Adapter Replacement PCB Kit	1R77010-1074
TriPlus RSH - Y-axis Front Cover Kit	1R77010-1075
TriPlus RSH - Vortexer Replacement PCB Kit	1R77010-1076
TriPlus RSH - Headspace 1000 µL O-ring Kit	1R77010-1077
TriPlus RSH - Headspace 2500 µL O-ring Kit	1R77010-1078
TriPlus RSH - Headspace 5000 µL O-ring Kit	1R77010-1079
TriPlus RSH - Cable BUS, 250 mm	1R77010-1080
TriPlus RSH - Cable BUS, 850 mm	1R77010-1081
TriPlus RSH - DC Cable, 2 m	1R77010-1082
TriPlus RSH - Y-Interface Cable for two Thermo Scientific GC's	1R77010-1083
TriPlus RSH - Z-Cover Kit	1R77010-1084
TriPlus RSH - Z-Cover Window Kit	1R77010-1085
TriPlus RSH - Z-Spindle Nut Kit	1R77010-1086
TriPlus RSH - Z-Motor Brake	1R77010-1087
TriPlus RSH - Z-Ribbon Cable Kit	1R77010-1088
TriPlus RSH - Pressure Regulator	1R77010-1089
TriPlus RSH - Connector TP RSH to Pressure Regulator	1R77010-1090
TriPlus RSH - Tube Connector from Pressure Regulator to 1/8-inch Tubing	1R77010-1091
TriPlus RSH - Purge Valve Kit	1R77010-1092
TriPlus RSH - 10 µL syringe for PTV, FN 57 mm, gauge 26, cone tip	365D0291
TriPlus RSH - 10 µL syringe for PTV, FN 57 mm, gauge 26, bevel tip	365D0391