

Thermo Scientific News

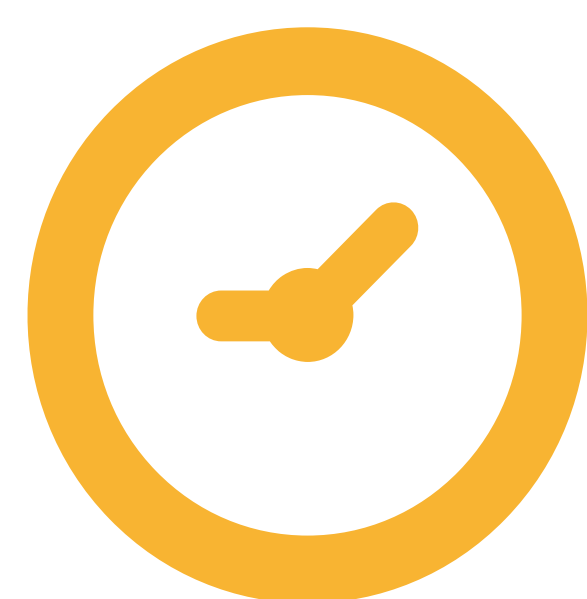
The Latest Evolution of the iQ Series!

May 7, 2019

Environmental and Process Monitoring

Issue 100

SERVICE TIPS that will SAVE you TIME!

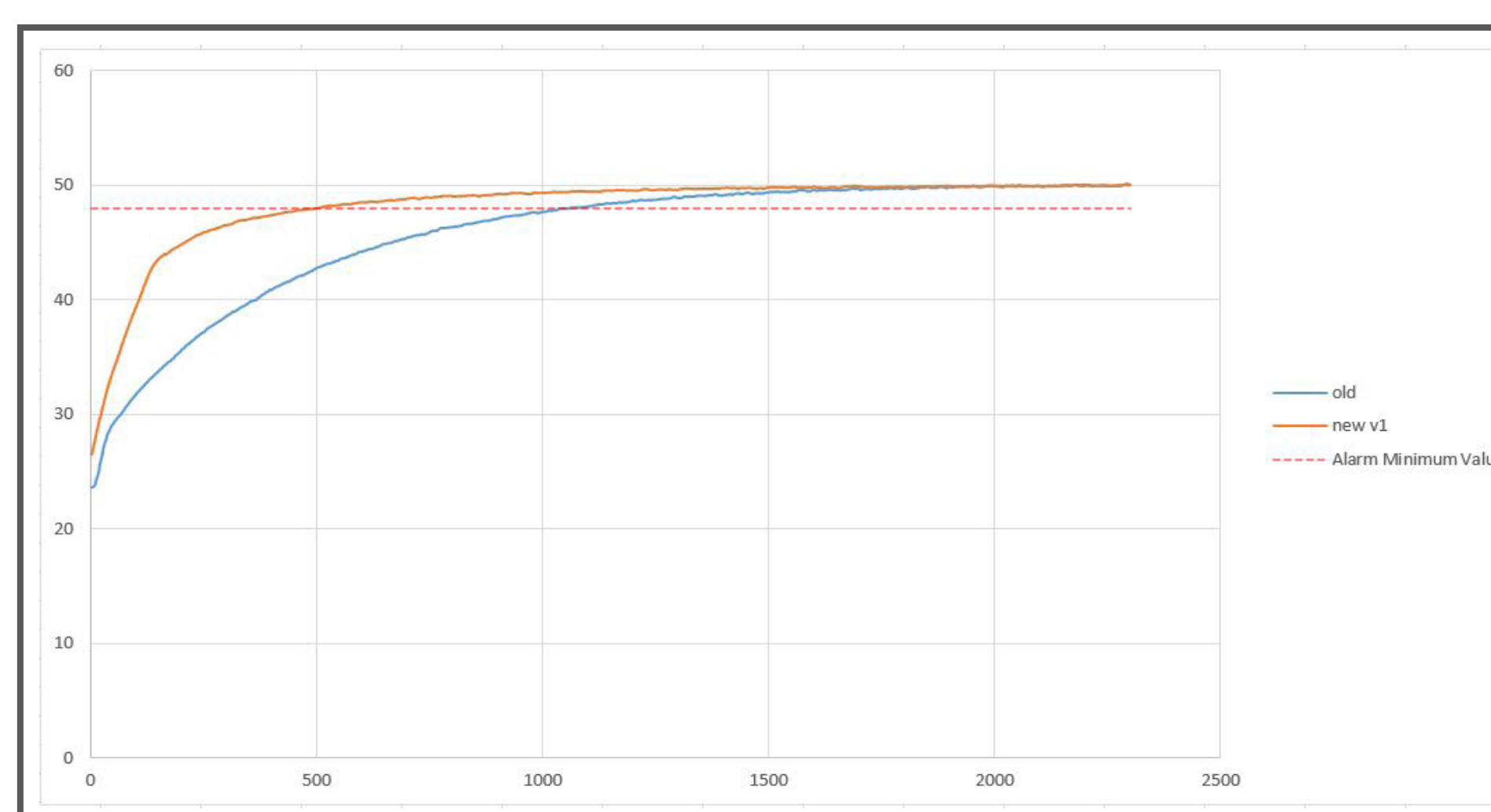
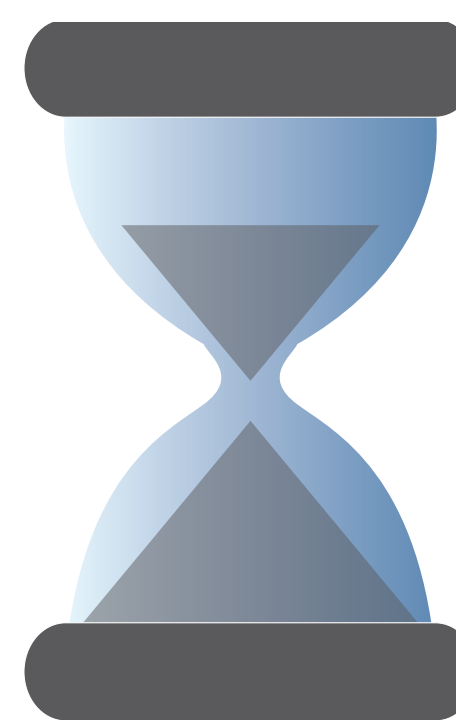


TIPS table

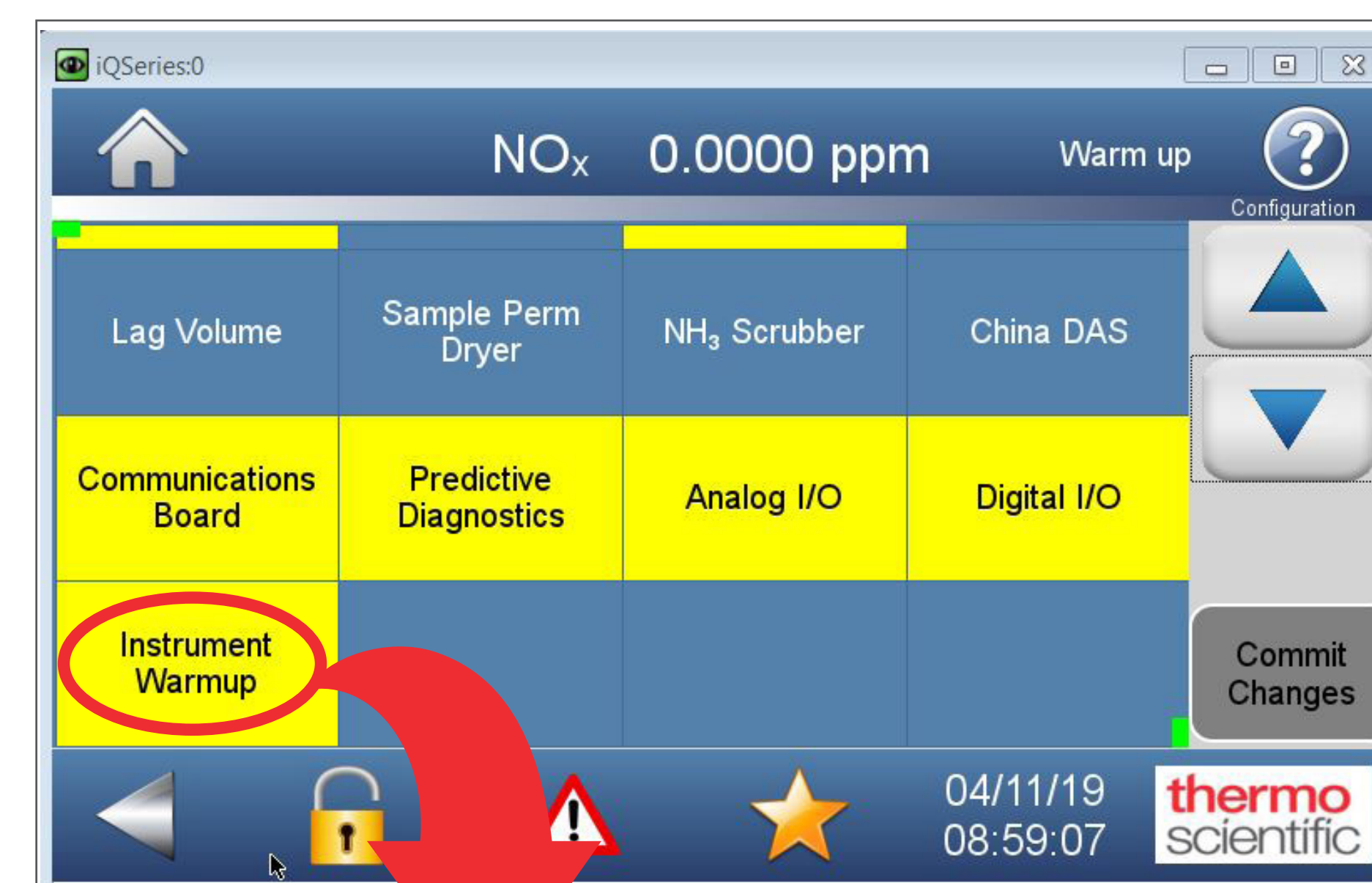
Product	Pump rebuild video	Step POL board	Leak check	Purge housing plumbing	Hydro-carbon kicker	PMT replacement	Reaction chamber cleaning
42iQ	✓	✓	✓			✓	✓
43iQ	✓	✓	✓	✓			
48iQ	✓	✓	✓		✓		
49iQ	✓	✓	✓				
146iQ	✓	✓	✓				
410iQ	✓	✓	✓				

Are you tired of waiting for the Warm-up?

Our cold-start Warm-up times have been reduced by an average of 10 minutes per instrument model!



Don't Like it? Disable it!



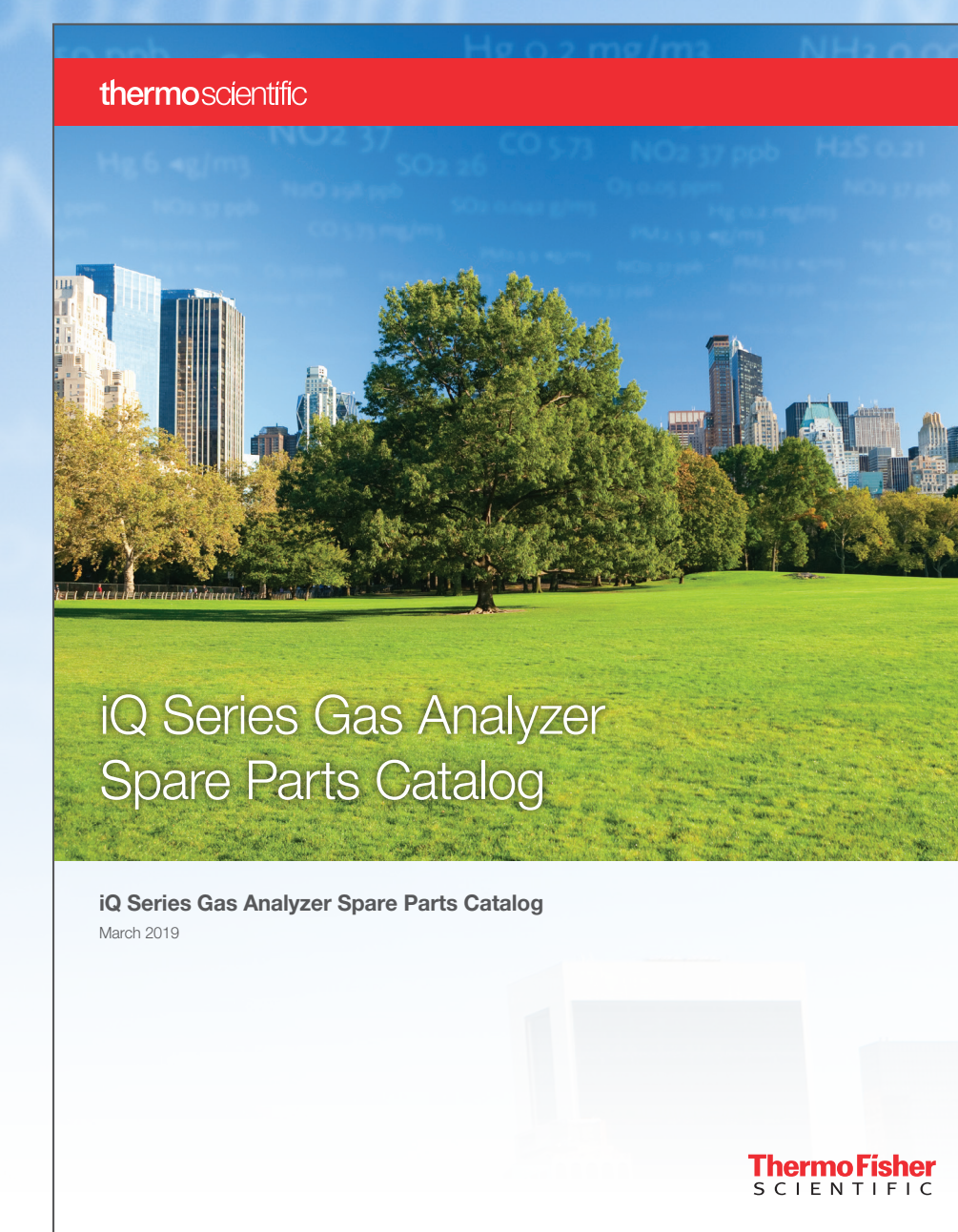
You can now disable the Warm-up feature in the instrument's Configuration menu. Remember, concentration readings are not accurate until the instrument stabilizes and is **Alarm FREE**.

Who can use an Alert Assistant?



FREE UP your time!

New Digital Parts Catalog Now Available!



42iQ Optical Bench DMC

Item	Description	P/N	Qty	Component
1a	ASSY. PMT COOLER DMC, 42iQ	117272-01	1	Spare Part
1b	ASSY. PMT COOLER DMC, 42iQLS	117272-04	1	Spare Part
2	ASSY. BASE, PHOTOMULTIPLIER	101324-00	1	Spare Part
3	O-RING, 1.739 ID, 1.879 OD, 070 CSD, VITON	4806	1	Spare Part
4	O-RING, 1.424 ID, 1.633 OD, 103 CSD, VITON	4892	1	Spare Part
6	FEM RUN TEE	9452	1	Spare Part
7	FTG. ELBOW, 1/4" OD, 1/8"NPT, SS	4430	1	Spare Part
8	CAP, 1/8" ID, 25 OD	4786	1	Spare Part
9	FITTING, UNION, CAJON, SS, 1/4"	9211	2	Spare Part
10	O-RING, 4/14 ID, 072 CSD, VITON	5212	2	Spare Part
11	FITTING, ELBOW, 1/8" OD, 1/8 NPT, SS (not shown)	4403	2	Spare Part
12	TUBING, 1/8, SS, RG TO OZONE CAP OUTLET	117402-00	1	Spare Part
13	PCB ASSY. DMC PMT COOLER	116902-00	1	Spare Part
14	REACTION CHAMBER, 42iQ	117127-00	1	Spare Part
17	TUBE, PHOTOMULTIPLIER, MODEL 42iQ	100606-00	1	Spare Part
18	TUBE, PHOTOMULTIPLIER, MODEL 42, TL, Y, 17	114050-00	1	Spare Part
20	NIPPLE, RECOVER	9453	1	Spare Part
21	ASSY. HEATER, 42iQ DMC	117012-00	1	Spare Part
5	SCREW, PAN HEAD, PHILLIPS, W/EXT WASHER, FT, #8-32, 3/4, SS	21-005034-0012	3	Not Sellable
15	SCREW, PAN HEAD, PHILLIPS, W/EXT WASHER, FT, #8-32, 3/8, SS	21-005087-0008	1	Not Sellable
16	TAPE, SEALANT, PIPE THREAD, 1/4 W	5766	A/R	Not Sellable
19	CAP, 1/8" ID	4785	2	Not Sellable

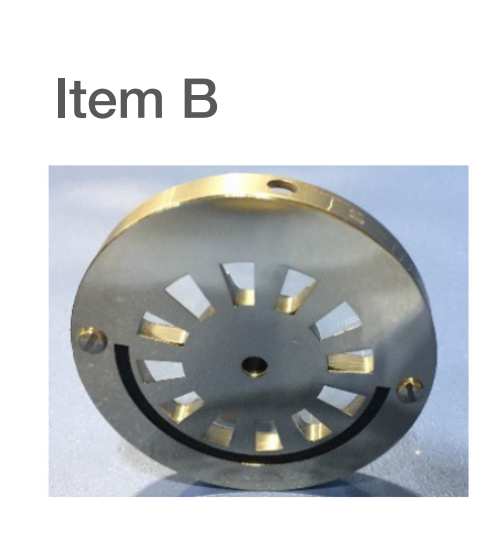
Component	Description
Consumable	Required every year
Spare part	Available for purchase
Optional	Not standard component
Not sellable	Not available for purchase

43iQ Optical Bench DMC

Item	Description	P/N	Qty	Component
4	ASSY. BENCH, OPTICAL, 43iQ, STANDARD	116556-00	1	Spare Part
9	PCB DMC PMT OPTICAL BENCH	118485-00	1	Spare Part
10	ASSY. HEATER, 43iQ	115948-01	1	Spare Part
11	LED ASSEMBLY	8050	1	Spare Part
12	ASSY. KICKER, HEATED, 43iQ	117035-00	1	Spare Part
13	TRIGGER PACK ASSEMBLY	117779-00	1	Spare Part
1	ASSY. BRACKET, DMC, 43iQ	116601-00	1	Not Sellable
2	SCREW, CAP, SOCKET HEAD, PARTIAL THREAD, #10-32, 2.75, SS	21-000352-0044	4	Not Sellable
3	CABLE ASSY., DETECTOR, UV	116607-01	1	Not Sellable
6	SCREW, PAN HEAD, PHILLIPS, W/EXT WASHER, FT, #8-32, 1/4, SS	21-005034-0004	1	Not Sellable
7	TIE, CABLE, NYLON, 4"	5385	2	Not Sellable

48iQ Motor Assembly

Item	Description	P/N	Qty	Component
10	IR SOURCE	7381	1	Consumable
1	WHEEL HOUSING	7544	1	Spare Part
3	ASSY. OPTICAL SWITCH	101714-00	1	Spare Part
6	MOTOR, STEP, NEMA 23	115552-00	1	Spare Part
7	ASSY. SOURCE, IR, 48iQ/19iQ	115976-00	1	Spare Part
8	WHEEL FILTER, TESTED, CO	7359	1	Spare Part
2	SCREW, MACHINE, FLAT HEAD, PHILLIPS, FULL THD, #8-32, 3/8, SS	21-004277-0008	2	Not Sellable
4	TAPE, SEALANT, PIPE THREAD, 1/4 WIDE, 0035 THICK	5766	A/R	Not Sellable
5	O-RING, 1/8"NPT, BARB, NYLON, 01070	116788-0100	1	Not Sellable
9	SCREW, PAN HEAD, PHILLIPS, W/EXT WASHER, #4-40, 3/8, SS	102175-00	2	Not Sellable
10	PLATE, MOTOR, 48iQ/19iQ	116268-00	1	Not Sellable
11	SCREW, MACHINE, PAN HEAD, PHILLIPS, FULL THD, #8-32, 1/4, SS	41453	1	Not Sellable
12	SCREW, CAP, SOCKET HEAD, FULL THREAD, #10-32, 1.00, SS	21-000352-0018	2	Not Sellable
13	SCREW, CAP, SOCKET HEAD, FULL THREAD, #8-32, 1/2, SS	21-000268-0008	2	Not Sellable
14	PLATE, DEFLECTOR, AIR	116768-00	1	Not Sellable
15	SCREW, PAN HEAD, PHILLIPS, W/EXT WASHER, FT, #8-32, 1/4, SS	104488-00	2	Not Sellable



ThermoFisher
SCIENTIFIC
The world leader in serving science