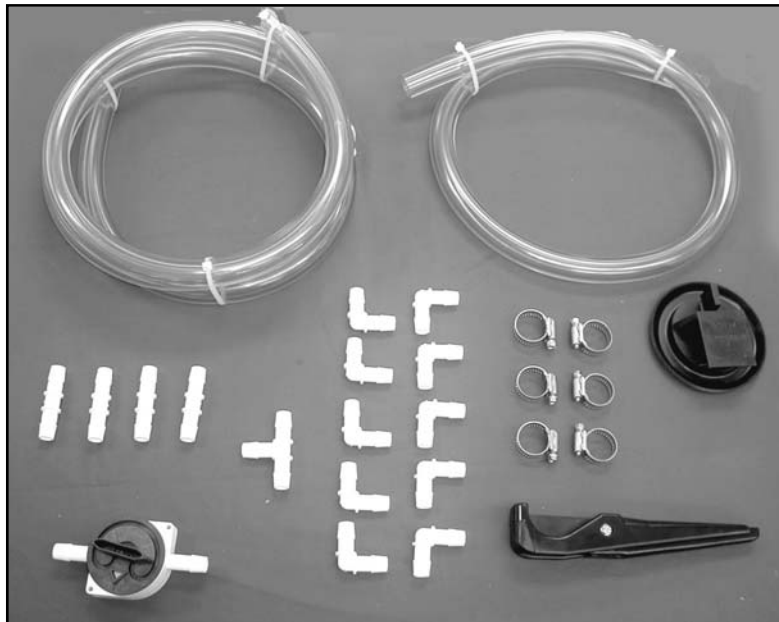
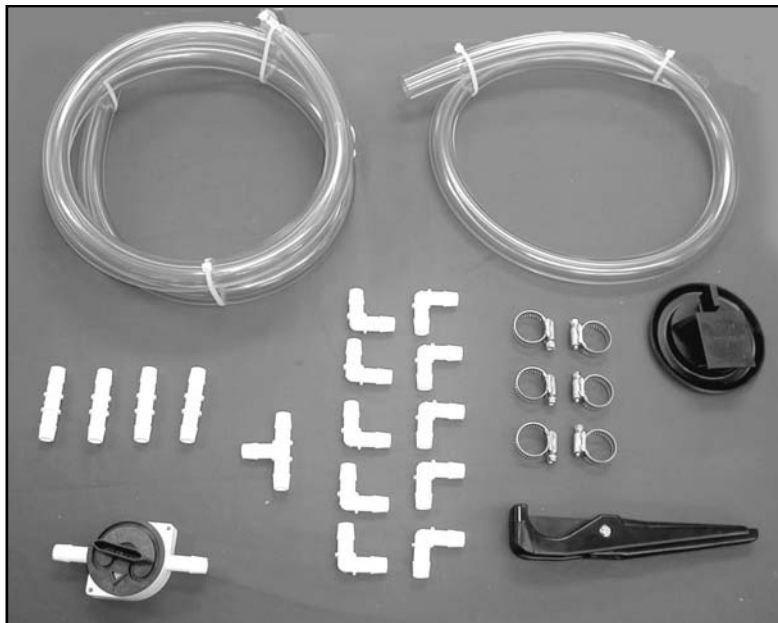

UTP-TYG

INSTRUCTION MANUAL





UTP-TYG

Universal tubing packages for SC110, SC110A, SC210A, SPD101B and SPD111V

Note: GCF400's sold separately

Quantity	Part Number & description
1	M30-0007-03 8' Vacuum tubing
1	M30-0007-02 4' Vacuum tubing
6	M30-0013-04 Hose clamps
10	M60-0044-11 Elbow fittings
4	M60-0044-14 Straight fittings
1	M60-0044-16 Tee fitting
1	TC100 Tubing cutter
1	FC400 Flask cover
1	BV130 Bleeder valve
1	Instructions

BASIC COMPONENT BENCH-TOP SET-UP

Section 1



SC110A, RVT400, and VLP Pump

OR



SPD111V OR SPD121P, RVT400, and VLP Pump



Step 1

Pour 750 ml of CryoCool® (SCC1) into RVT unit. (RVT400, RVT4104).



Step 2

Insert GCF400 into refrigeration chamber of RVT400 or RVT4104 (not pictured).



Step 3

Install white foam insulating ring by pushing down into black 'O' Ring of refrigeration unit.



Step 4

Install black flask cover (FC400) by pushing cap down onto rim of glass jar.



Universal Tubing Packages Available:

- 1) UTP-TYG - For SC110, SC110A, SPD101B and SPD111V
- 2) CAFS400 - For SC110, SC110A, SPD101B and SPD111V
- 3) UTP-TEF0.05
- 4) UTP-TEF0.75

Step 5

Installation of Tubing Packages

1. Using appropriate vacuum tubing (4' or 8' clear tygon, 4' corrugated or 3' TEFLON®) install tubing onto fitting

(Note: Do Not Twist Tubing on, apply small amount of vacuum grease to fitting, then push on) see Pic 1A



Picture 1

Step 5A

Picture 1

Note: Do Not Twist tubing onto fitting. Apply small amount of vacuum grease to the nipple end of fitting or apply small amount of CryoCool® to nipple end. Then “push” vacuum tubing onto fitting!



Step 6

Measure approximate length of tubing to be cut. (There is no standard length to be given, each installation can be different due to location of equipment).



Step 7

Insert 90 ° (M60-0044-11) elbow into vacuum tubing (See step 5A for proper installation of tubings to fitting).



Step 8

Insert remaining PVC vacuum tubing into 90 ° elbow (M60-0044-11).



Step 9

Measure approximate length of tubing to be cut (Normal cuts should be made even with top of refrigeration unit).



Step 10

Insert 90 ° elbow (M60-0044-11, see Step 5A).



Step 11 & 12

Inset remaining vacuum tubing onto 90 ° elbow (See Step 5A).



Step 11 & 12

Measure distance to flask cap even with hole in back of cap and then cut tubing (See Step 5A).

Step 12



Step 13

Insert 90 ° elbow (M60-0044-11) into vacuum tubing (See Step 5A).



Step 14 & 15

Push 90 ° elbow into empty hole in back of flask cover.



Step 15

Tubing & Fitting now inserted into flask cover.



Step 16

Insert 90 ° (M60-0044-11) elbow into remaining empty hole of flask cover.



Step 17

Insert remaining vacuum tubing onto 90 ° elbow (See Step 5A). Instructions for proper method of installing vacuum tubing onto fitting.

This PVC tubing should reach to vacuum inlet of vacuum pump.



Installation of manual bleeder valve. (BV130)

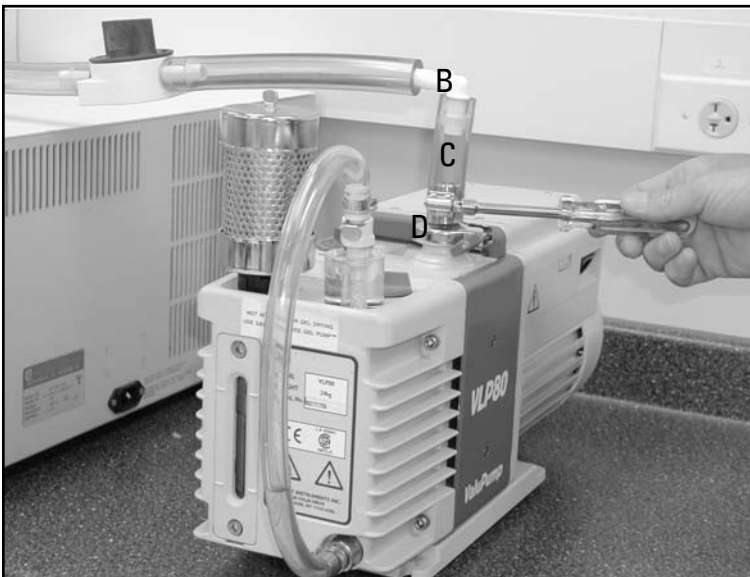
Step 18

Measure approximate distance needed to install BV130 valve. Push both ends of valve into vacuum tubing (See Step 5A).



Step 18

Complete installation of BV130.



Step 19

Cut tubing A to approximate length to reach vacuum inlet D of VLP pump, then install 90 ° elbow B. Cut vacuum tubing C to level height of tube A and install onto vacuum inlet D of VLP & 90 ° elbow. NOTE: Clamp provided must be installed on tube C where it meets vacuum inlet D of VLP pump.



Finished installation of component system using UYP-TTG tubing package.

NOTE: This same system could have been installed using a CAFS400 kit.

ADDENDUM



A. How to Hook up DTK120R/SCT120



- B.
1. Disconnect existing vacuum line from BV130
 2. using supplied "L" vacuum line, connect "L" line to bottom fitting on SCT120 trap. Connect remaining line to fitting of BV130 (See Figure 5A instructions for proper method of installing vacuum tubing onto fitting).



C.
A connection now must be made from the top fitting on the SCT120 back to the vacuum pump.

1. You will need 2 elbow fittings (Part # M60-0044-11)
2. 2 pieces 1/2 ID vacuum tubing approximately 3" & 11".

All tubing should be cut to fit.



Finished installation of
DTK120R/SCT120



VN100DDA Installation

Place VN100DDA on top of RVT unit.
(RVT4104 not shown)



Step 1

1. Remove vacuum tubing containing BV130 from back of flask cover.
2. Push on vacuum line containing BV130 into hole marked vacuum pump (See Figure 1 Step1).



Figure 1

Step 1

Existing vacuum line & BV130.



Figure 2

Step 2

A. Parts needed:

- 3– M60-0044-11 (90 °) Elbows
- 1– 5” Vacuum tubing
- 1– 6” Vacuum tubing
- 1– 8” Vacuum tubing

Cut and assemble according to (Figure 2).

B. Proceed to Step 3 Figure 3 push on vacuum tubing (5”) into remaining hole marked.

Vapor Trap: Add flask cap (FC400) to 8” vacuum tubing and 90 ° elbow.

See Figure 1 step 1 for completed vapornet assembly.



Figure 3

Step 3

Push on 5” PVC tubing into hole marked “vapor trap”.

See figure 5A instructions for proper method of installing vacuum tubing onto fittings.



Figure 4

Step 4
Final assembly of VN100DDA onto
SpeedVac® System!

450 Fortune Boulevard, Milford, MA 01757 USA
Tel: (508) 482-7000 • 1-866-984-3766 • Fax: (508) 634-2199
www.thermo.com

Thermo
ELECTRON CORPORATION

UTP-TYG-3000 Rev. A
© 2003 Thermo Savant

A Thermo Electron business