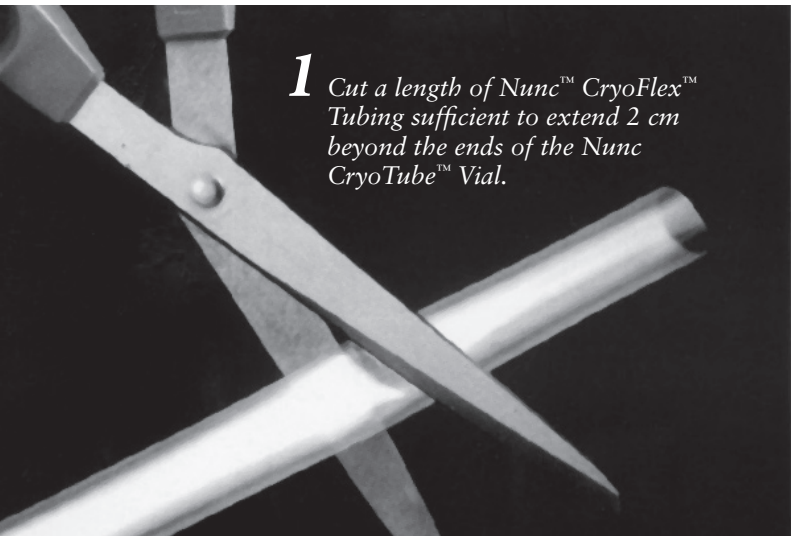
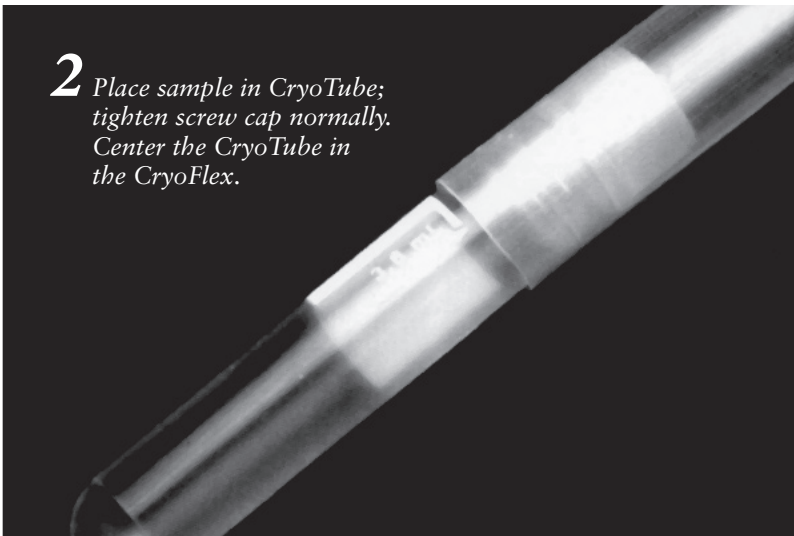


# Thermo Scientific Nunc CryoFlex Tubing

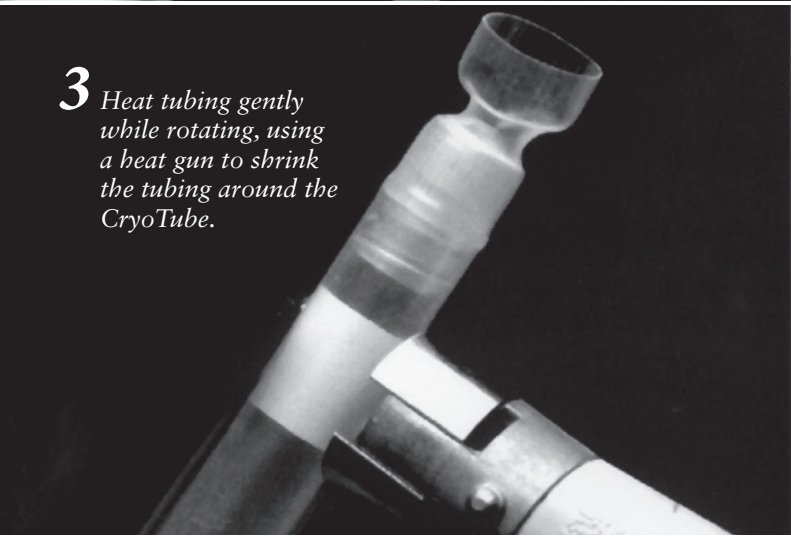
## – Instructions for Use



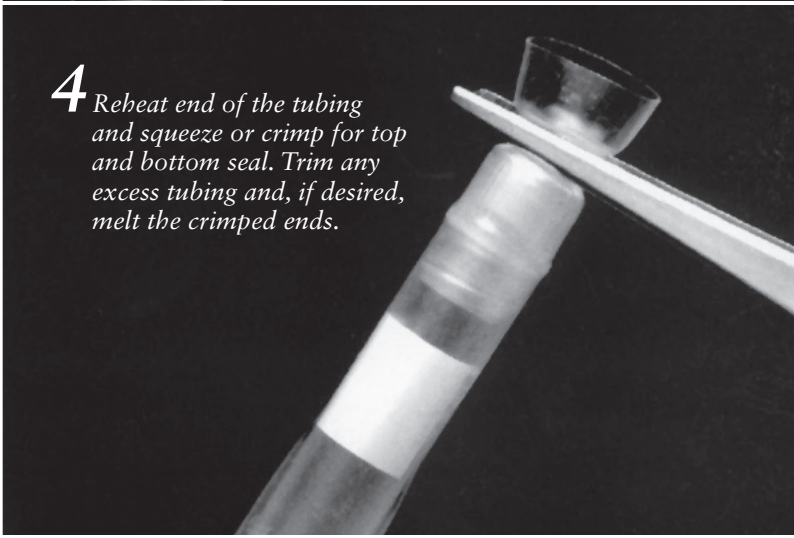
**1** Cut a length of Nunc™ CryoFlex™ Tubing sufficient to extend 2 cm beyond the ends of the Nunc CryoTube™ Vial.



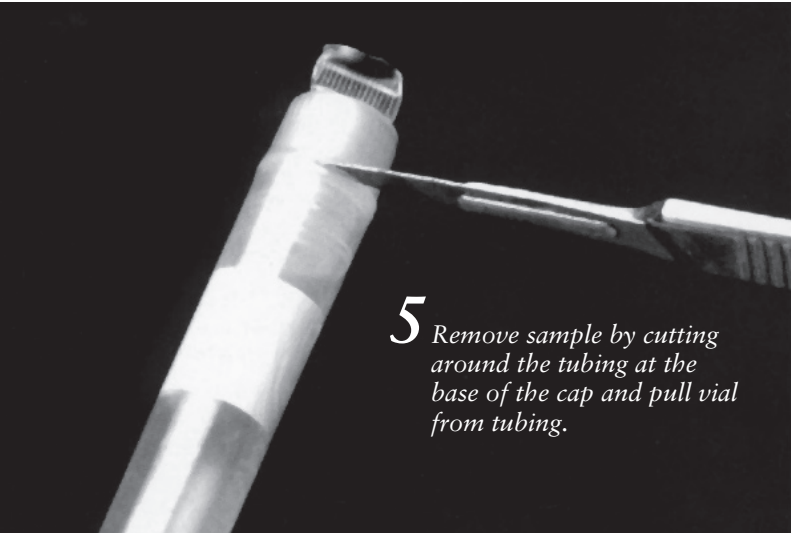
**2** Place sample in CryoTube; tighten screw cap normally. Center the CryoTube in the CryoFlex.



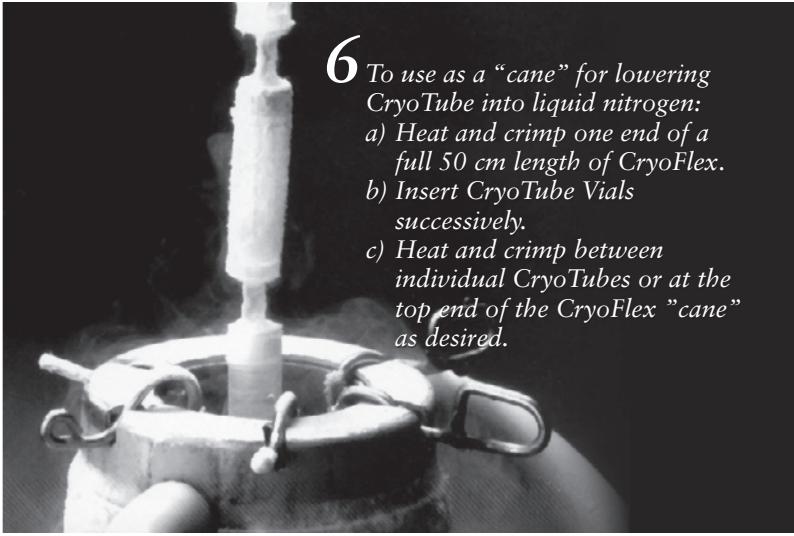
**3** Heat tubing gently while rotating, using a heat gun to shrink the tubing around the CryoTube.



**4** Reheat end of the tubing and squeeze or crimp for top and bottom seal. Trim any excess tubing and, if desired, melt the crimped ends.



**5** Remove sample by cutting around the tubing at the base of the cap and pull vial from tubing.

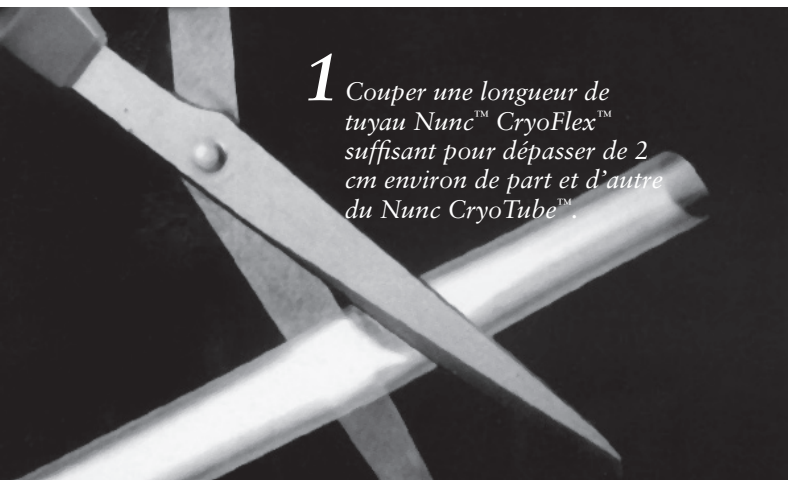


**6** To use as a "cane" for lowering CryoTube into liquid nitrogen:

- Heat and crimp one end of a full 50 cm length of CryoFlex.
- Insert CryoTube Vials successively.
- Heat and crimp between individual CryoTubes or at the top end of the CryoFlex "cane" as desired.

# Thermo Scientific Nunc CryoFlex

– Instructions pour l'usage



**1** Couper une longueur de tuyau Nunc™ CryoFlex™ suffisant pour dépasser de 2 cm environ de part et d'autre du Nunc CryoTube™.



**2** Mettre l'échantillon dans le CryoTube; visser normalement le capuchon. Placer le CryoTube à l'intérieur de la longueur de CryoFlex.



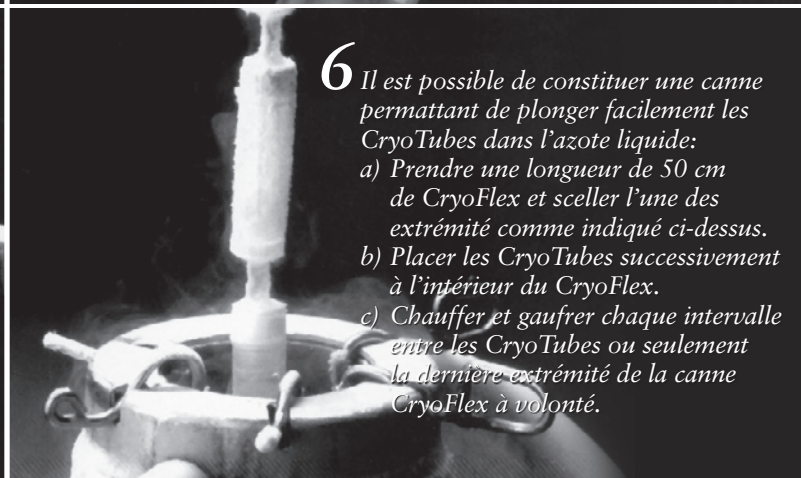
**3** Chauffer le tuyau doucement à l'aide d'un pistolet chauffant, en vue de rétrécir le tuyau autour de CryoTube.



**4** Chauffer à nouveau les extrémités du CryoFlex et presser ou gaufrer pour sceller définitivement les deux extrémités gaufrées.



**5** Pour retirer l'échantillon, découper le CryoFlex autour du joint silicone du CryoTube et dévisser le capuchon de ce dernier.



**6** Il est possible de constituer une canne permettant de plonger facilement les CryoTubes dans l'azote liquide:  
a) Prendre une longueur de 50 cm de CryoFlex et sceller l'une des extrémités comme indiqué ci-dessus.  
b) Placer les CryoTubes successivement à l'intérieur du CryoFlex.  
c) Chauffer et gaufrer chaque intervalle entre les CryoTubes ou seulement la dernière extrémité de la canne CryoFlex à volonté.

© 2010 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.

**Asia:** China Toll-free: 800-810-5118 or 400-650-5118; India: +91 22 6716 2200, India Toll-free: 1 800 22 8374; Japan: +81 3 3816 3355

Other Asian countries: 65 68729717

**Europe:** Austria: +43 1 801 40 0; Belgium: +32 53 73 42 41; Denmark: +45 4631 2000; France: +33 2 2803 2180;

Germany: +49 6184 90 6940, Germany Toll-free: 08001-536 376; Italy: +39 02 02 95059 or 434-254-375;

Netherlands: +31 76 571 4440; Nordic/Baltic countries: +358 9 329 100; Russia/CIS: +7 (812) 703 42 15;

Spain/Portugal: +34 93 223 09 18; Switzerland: +41 44 454 12 12; UK/Ireland: +44 870 609 9203

**North America:** USA/Canada +1 585 586 8800; USA Toll-free: 800 625 4327

**South America:** USA sales support: +1 585 899 7198

**Countries not listed:** +49 6184 90 6940 or +33 2 2803 2180

MNLSPCRYOFL70246 0310  
70246, Ver. 4.1

**Thermo**  
SCIENTIFIC