

Procedure for Transferring Dodecafluorocyclohexane from the Ampule

PRODUCT No.
13,393-0

revised 4 / 96

2 pages

Dodecafluorocyclohexane (C₆F₁₂) is a colorless solid which sublimes **very rapidly at room temperature**. Appreciable loss occurs within minutes of breaking the ampule.

The following procedure is recommended for opening the ampule and transferring material without significant loss.

- 1• Cool the ampule in a dry-ice bath for at least 15 minutes.
- 2• Use gloves and a towel to break the ampule safely; have an appropriate spatula on hand to transfer the material.
- 3• Remove the cooled ampule from the dry-ice bath wearing gloves. Place the scored end of the ampule up. Put the towel over the ampule (to contain any glass fragments) and break the ampule.
- 4• Quickly transfer the material using the spatula.

Aldrich Chemical Company, Inc.

1001 West Saint Paul Ave., Milwaukee, WI 53233

Telephone 414-273-3850

Fax 414-273-4979

Internet aldrich@sial.com

800-231-8327

800-962-9591

Aldrich warrants that its products conform to the information contained in this and other Aldrich publications. Purchaser must determine the suitability of the product for its particular use. See reverse side of invoice or packing slip for additional terms and conditions of sale.

