

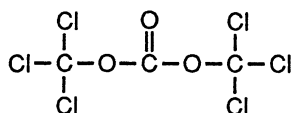
Triphosgene

PRODUCT No.
33,075-2

revised 4 / 96

6 pages

PROPERTIES



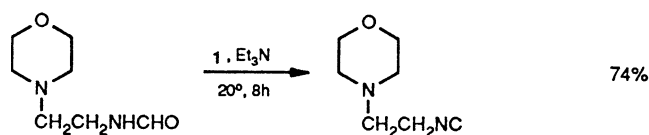
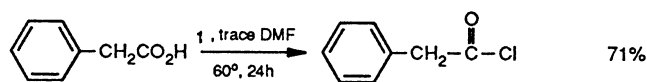
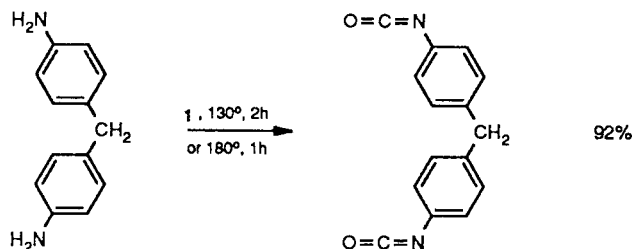
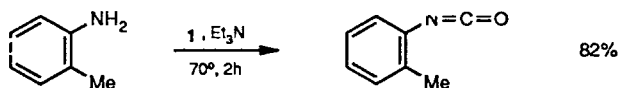
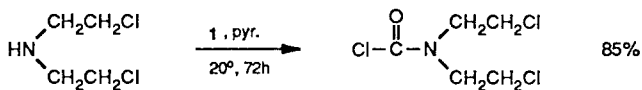
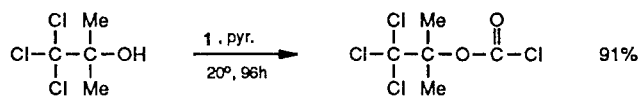
Appearance white crystalline solid
F.W. 296.75
b.p. 203-206°V
m.p. 81-83°C
Solubility alcohol, benzene, ether, hexane,
tetrahydrofuran; decomposes in hot
water, reacts slowly with cold water

APPLICATION

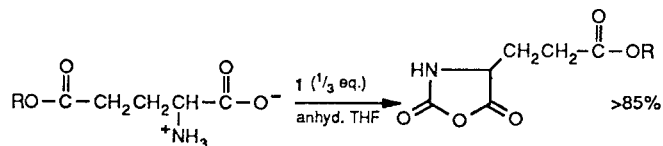
Phosgene has long been of tremendous importance to organic chemists for a wide variety of synthetic applications (e.g., chloroformylation, carbonylation, chlorination, and dehydration). However, because of its highly toxic nature, phosgene gas is difficult to handle safely in the laboratory.

Triphosgene [bis(trichloromethyl) carbonate, **1**] is a stable, crystalline solid which has proved to be a useful substitute for phosgene.¹ It is safer and more convenient to handle, transport and store. Exact amounts may be weighed easily and used to perform desired chemical transformations. Reaction requires only one-third equivalent of triphosgene in most cases.¹⁻³

Below are examples of some reactions employing triphosgene.¹



Recently, triphosgene has been shown to react with several α -amino acids to give the corresponding N-carboxyanhydrides in good yields.³



TOXICITY, HANDLING AND STORAGE

Triphosgene is severely irritating to the eyes and skin. Avoid contact and inhalation. Use adequate personal protection. Handle only in a chemical fume hood. Keep tightly closed. Store in a cool, dry place.

REFERENCES

- 1• Eckert, H.; Forster, B. *Angew. Chem.* **1987**, *99*, 922 (*Angew. Chem., Int. Ed. Engl.* **1987**, *26*, 894).
- 2• Eckert, H. Ger. Offen. DE 3 440 141, 1986; *Chem. Abstr.* **1987**, *106*, 4294d.
- 3• Daly, W.H.; Poché D. *Tetrahedron Lett.* **1988**, *29*, in press.

For pricing and ordering information, call 800-558-9160



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.

