

REFERENCES:

- 1• Pedersen, C.J. *Aldrichimica Acta* **1971**, *4*, 1.
- 2• Gokel, G.W.; Durst, H.D. *ibid.* **1976**, *9*, 3.
- 3• Valentine, J.S.; Curtis, A.B. *J. Am. Chem. Soc.* **1975**, *97*, 224.
- 4• Christensen, J.J. *Science* **1971**, *459*, 174.
- 5• Pedersen, C.J. *Fed. Proc.* **1965**, *27*, 1305.
- 6• Liotta, C.L.; Harris, H.P. *J. Am. Chem. Soc.* **1974**, *96*, 2250.
- 7• Cook, F.L.; Bowers, C.W.; Liotta, C.L. *J. Org. Chem.* **1974**, *39*, 3416.
- 8• Zubrick, J.W.; Dunbar, B.I.; Durst, H.D. *Tetrahedron Lett.* **1975**, 71.
- 9• Landini, D.; Maia, A.M.; Montanari, F.; Pinsi, F.M. *Gazz Chim Ital.* **1975**, *105*, 863.
- 10• Landini, D.; Montanari, F. *Chem. Commun.* **1974**, 879.
- 11• Herriott, A.W.; Picker, D. *J. Am. Chem. Soc.* **1975**, *97*, 2345.
- 12• Makosza, M.; Ludwikow, M. *Angew. Chem.* **1974**, *86*, 744.
- 13• Moss, R.A.; Pilkiewica, F.G. *J. Am. Chem. Soc.* **1974**, *96*, 5632.
- 14• Fedorynski, M. *Synthesis* **1977**, 784.
- 15• Matsuda, T.; Koida, K. *Bull. Chem. Soc. Jpn.* **1973**, *46*, 2259.
- 16• Svoboda, M.; Hapala, J.; Zarada, J. *Tetrahedron Lett.* **1972**, 265.
- 17• Bartsch, R.A.; Wiegers, K.E. *ibid.* **1972**, 3819.
- 18• Sam, D.J.; Simmons, H.E. *J. Am. Chem. Soc.* **1974**, *96*, 2252.
- 19• Durst, H.D. *Tetrahedron Lett.* **1967**, 2421.
- 20• Knöchel, A.; Oehler, J.; Rudolph, G. *Angew. Chem.* **1975**, *87*, 3167.
- 21• Komarynsky, M.A.; Weissman, S.I. *J. Am. Chem. Soc.* **1975**, *97*, 1589.
- 22• Nelson, G.W.; von Zelewsky, A. *ibid.* **1975**, *97*, 6279.
- 23• Sam, D.J.; Simmons, H.E. *ibid.* **1972**, *94*, 4024.
- 24• Hunter, D.H.; Lee, W.; Sim, S.K. *Chem. Commun.* **1974**, 1018.
- 25• Sepp, D.T.; Scherer, K.V.; Weber, W.P. *Tetrahedron Lett.* **1974**, 2983.
- 26• *Chem. Eng. News*, March 2, 1970, p 26.
- 27• Roeske, R.W.; Gesellchen, P.D. *Tetrahedron Lett.* **1976**, 3369.
- 28• Leong, B.K.J. *Chem. Eng. News*, January 27, 1975, p 5.

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.

