

MCPBA

(m-Chloroperoxybenzoic Acid)

PRODUCT No.
C6,270-0

revised 4 / 96

6 pages

PROPERTIES:

White powder

F.W. 172.57

m.p. 92-94° (dec.)

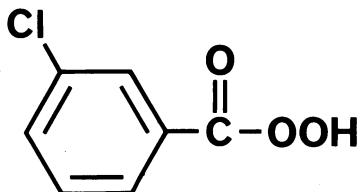
Contains up to 15% *m*-chlorobenzoic acid

Slight pungent odor

Bulk density 0.56g/cc.

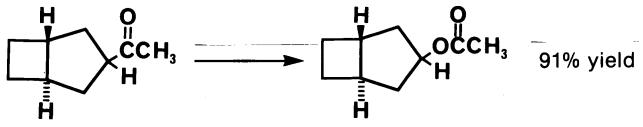
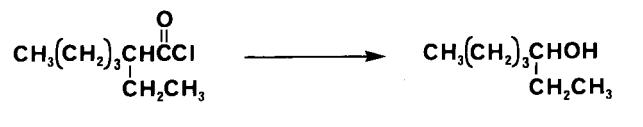
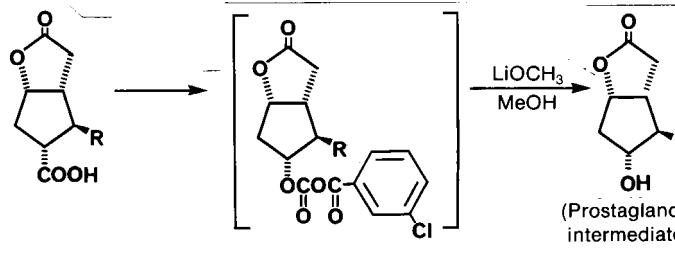
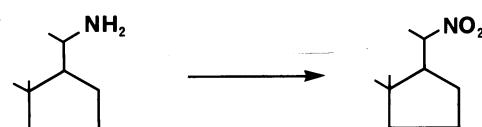
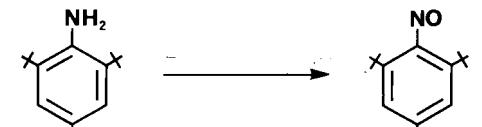
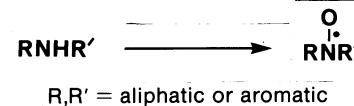
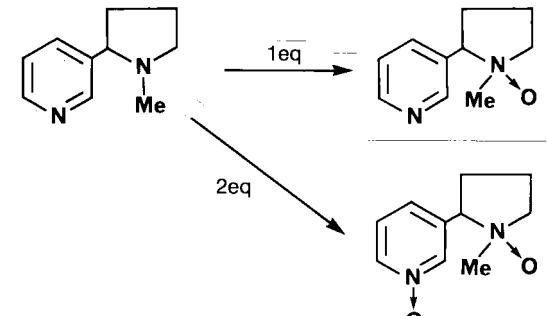
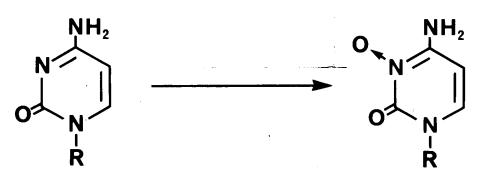
pH of a saturated aqueous solution at 25°C: 4.5

pK_a (in water at 25°C): 7.57



MCPBA (*m*-chloroperoxybenzoic acid), an oxidizing agent, is superior to hydrogen peroxide or other peracids in terms of reactivity, stereoselectivity, and in purity and yield of products. The scope of its reactivity is illustrated in the following table.

| Reactant | Product | Example | Ref. |
|--|--|---------|-------|
| Olefins | Expoxides ^a | | 1-7 |
| Cyclopropenes | α,β -Unsaturated aldehydes and/or ketones | | 8,9 |
| α,β -Unsaturated ketones and esters | Epoxides | | 10 |
| Disubstituted acetylenes | Oxirenes ^b | | 11,12 |
| Imines | Oxaziranes | | 13-16 |

| Reactant | Product | Example | Ref. |
|-------------------------------------|------------------------------|--|-------|
| Ketones (Baeyer-Villiger oxidation) | Esters |  | 17-21 |
| Acid Chlorides | Alcohols |  | 22 |
| Acids | Alcohols |  | 23 |
| Primary alkyl amines | Nitro alkanes ^c |  | 24 |
| Primary aromatic amines | Aromatic nitroso compounds |  | 25 |
| Secondary amines | Nitroxide radicals |  | 26 |
| Tertiary amines | <i>N</i> -oxides |  | 27 |
| Nucleic acid components | <i>N</i> -oxide ^d |  | 28-31 |

| Reactant | Product | Example | Ref. |
|--|--|--|-------|
| <i>N</i> -substituted aziridines | Olefin ^e | | 17-21 |
| 2-Pyridine-acetates | Corresponding glycolates ^f | | 34 |
| Sulfides | Sulfoxides or sulfones ^g | | 35-40 |
| Carbodiimides | Diaziridinones | | 41 |
| Ketals | Ortho esters | | 42 |
| Trimethylsilyl vinyl and allyl systems | Trimethylsilyl epoxides (latent precursors to carbonyl groups) | | 43-46 |
| Iminoethers | Esters and hydroxylamines | $\text{HC-O}^{\text{Me}} \quad \text{H}_3\text{O}^+$ $(\text{CH}_3)_3\text{C}-\text{N} \quad (\text{CH}_3)_3\text{C}-\text{NHOH}$ | 47 |
| α -Hydroxy ketones | Aldehydes and acids | | 48 |
| Mono-, di-, and trimethoxybenz aldehydes | Formate esters | | 49 |

| Reactant | Product | Example | Ref. |
|--------------------------------|-----------------------------------|---------|------|
| α -Diazoketones | α -Diketones | | 50 |
| β -Lactam acid chlorides | Aryl- β -lactam derivatives | | 51 |
| Secondary alcohols | Ketones | | 52 |
| Erythro thioether | Sulfone | | 53 |
| Terminal olefines | Primary alcohols | | 54 |
| Aromatics Hydrocarbons | Arene dioxides | | 55 |

- a In nonconjugated dienes the more substituted double bond is selectively epoxidized.
 b Oxirenes break down to ketones, carboxylic acids or esters depending on reaction conditions.
 c The yields of nitroalkanes decrease in the order: *tert*-alkyl > *sec*-alkyl > *n*-alkyl.
 d I.R. Subbaraman and co-workers²¹ reported that cytosine, adenine and their derivatives are oxidized to N-oxides while uracil, thymine, guanosine and their derivatives give ring-cleavage products. However, M.R. Harden *et al.*²² indicated that N(1)-oxides were obtained from adenine, cytosine and uracil derivatives while guanine derivatives yielded the N(3)-oxides.

- e N-substituted zairidines are presumably oxidized to the corresponding N-oxides. This reaction is successfully used in the sterospecific deamination of N-alkylaziridines to olefins.
 f the 3- and 4-pyridyl isomers gave the corresponding N-oxides in high yields.
 g Yields of sulfones or sulfoxides are excellent even in the presence of amino,²⁷ olefinic or acetylenic²⁸ moiety.

REFERENCES:

- Paquette, L.A.; Barrett, J.H. *Org. Syn.* **1969**, *49*, 62.
- Swern, D. *Chem. Rev.* **1949**, *45*, 1.
- Fieser, L.; Fieser, M. *Reagents for Org. Syn.* **1968**, *1*, 1136.
- Haywood-Farmer, J.; Friedlander, B.T.; Hammond, L.M. *J. Org. Chem.* **1973**, *38*, 3145.
- Anderson, W.K.; Veysoglu, T. *ibid.* **1973**, *38*, 2267.
- Ikegami, S.; *et al.* *Tetrahedron Lett.* **1980**, *21*, 3587.
- Kende, A.S.; Blacklock, T.J. *ibid.* **1980**, *21*, 3119.
- Ciabattoni, J.; Kocienski, P.J. *J. Am. Chem. Soc.* **1969**, *91*, 6534.
- Kocienski, P.J.; Ciabattoni, J. *J. Org. Chem.* **1974**, *39*, 388.
- Schwartz, N.N.; Blumbers, J.H. *ibid.* **1964**, *29*, 1976.
- Stille, J.K.; Whitehurst, D.D. *J. Am. Chem. Soc.* **1964**, *86*, 4871.
- Ciabattoni, J.; Campbell, R.A.; Renner, C.A. *ibid.* **1970**, *92*, 3826.
- Emmons, W.D. *ibid.* **1956**, *78*, 6208.
- Padwa, A. *ibid.* **1965**, *87*, 4365.
- Madan, V.; Clapp, L.B. *ibid.* **1969**, *91*, 6078.
- Madan, V.; Clapp, L.B. *ibid.* **1970**, *92*, 4902.
- Meinwald, J.; Tufariello, J.J.; Hurst, J.J. *J. Org. Chem.* **1964**, *29*, 2914.
- Palmer, B.W.; Fry, A. *J. Am. Chem. Soc.* **1970**, *92*, 2580.
- Syrkin, Y.K.; Moiseev, I.I. *Russ. Chem. Rev.* **1960**, *29*, 193.
- Marshall, J.A.; Ellison, R.H. *J. Org. Chem.* **1975**, *40*, 2070.
- Hudrik, P.R.; *et al.* *J. Am. Chem. Soc.* **1980**, *102*, 6894.
- Denney, D.B.; Sherman, N. *J. Org. Chem.* **1965**, *30*, 3760.

REFERENCES (CONTINUED):

- 23• Rosen, P.; et. al. 166th National Meeting of the American Chemical Society, Chicago, IL, August, 1973, MEDI Abstr. 61.
- 24• Robinson, C.H.; Milewich, L.; Hofer, P. *J. Org. Chem.* **1966**, *31*, 524.
- 25• Terabe, S.; Konaka, R. *J. Chem. Soc., Perkin Trans. 2* **1973**, 369.
- 26• Chapelet-Letourneux, G.; Lemaire, H.; Rassat, A. *Bull. Soc. Chim. France* **1965**, 3283.
- 27• Craig, J.C.; Purushothaman, K.K. *J. Org. Chem.* **1970**, *35*, 1721.
- 28• Delia, T.J.; Olsen, J.J.; Brown, G.B. *ibid.* **1965**, *30*, 2766.
- 29• Subbaraman, L.R.; Subbaraman, J.; Behrman, E.J. *Biochemistry* **1969**, *8*, 2059.
- 30• Harnden, M.R.; Brown, A.G.; Verg Hodge, R.A. *J. Chem. Soc., Perkin Trans. I* **1973**, 333.
- 31• Deady, L.W. *Synth. Commun.* **1977**, *7*, 509.
- 32• Padwa, A.; Hamilton, L. *J. Org. Chem.* **1966**, *31*, 1995.
- 33• Heine, H.W.; Myers, J.D.; Peltzer III, F.T. *Angew. Chem., Int. Ed. Engl.* **1970**, *9*, 374.
- 34• Skattebol, L.; Boulette, B. *J. Org. Chem.* **1969**, *34*, 4150.
- 35• Curci, R.; Giovine, A.; Modena, G. *Tetrahedron* **1966**, *22*, 1235.
- 36• Brown, D.J.; Ford, P.W. *J. Chem. Soc. (C)* **1969**, 2720.
- 37• Russell, G.A.; Ochrymowycz, L.A. *J. Org. Chem.* **1970**, *35*, 2106.
- 38• Trost, B.M.; Mao, M.R.T. *Tetrahedron Lett.* **1980**, *21*, 3523.
- 39• Ager, D.J. *ibid.* **1980**, *21*, 4759.
- 40• Heissler, D.; Richl, J.-J. *ibid.* **1980**, *21*, 4707.
- 41• Greene, F.D.; Bergmark, W.R.; Pazos, J.F. *J. Org. Chem.* **1970**, *35*, 2813.
- 42• Gaoni, Y. *J. Chem. Soc. (C)* **1968**, 2934.
- 43• Stork, G.; Jung, M.E. *J. Am. Chem. Soc.* **1974**, *96*, 3682.
- 44• Pennanen, S.I. *Tetrahedron Lett.* **1980**, *21*, 657.
- 45• Paquette, L.A.; et al. *J. Org. Chem.* **1980**, *45*, 3028.
- 46• Magnus, P.; Ehlinger, E. *J. Am. Chem. Soc.* **1980**, *102*, 5004.
- 47• Aue, D.H.; Thomas, D. *Tetrahedron Lett.* **1973**, 1807.
- 48• Greibrokk, T. *Acta Chem. Scand.* **1973**, *27*, 3365.
- 49• Sargent, M.V.; Godfrey, I.M. *J. Chem. Soc., Perkin Trans. 1* **1974**, 1353.
- 50• Curci, R.; DiFuria, F.; Ciabattoni, J.; Concannon, P.W. *J. Org. Chem.* **1974**, *39*, 3295.
- 51• Bose, A.K.; Kapur, J.C. *Tetrahedron Lett.* **1973**, 1811.
- 52• Cellia, J.A.; Kelley, J.A.; Kenahan, E.F. *J. Org. Chem.* **1975**, *40*, 1860.
- 53• Alcudia, F.; et al. *J. Chem. Soc., Perkin Trans. 2* **1979**, 564.
- 54• Kumada, M.; Tamao, K.; Kakui, T. *J. Am. Chem. Soc.* **1978**, *100*, 2268.
- 55• Griffin, G.W.; Ishikawa, K. *Angew. Chem., Int. Ed. Engl.* **1977**, *16*, 171.

SOLUBILITY, g / 100ml

| | |
|----------------------------|------|
| Benzene | 8.0 |
| Hexane | 1.4 |
| Methylene chloride | 11.2 |
| Chloroform | 9.8 |
| Carbon tetrachloride | 2.1 |
| 1,2-Dichloroethane | 10.6 |

| | |
|--------------------------|-------|
| Diethyl ether | 89.4 |
| Ethyl alcohol | 113.0 |
| tert-Butyl alcohol | 69.0 |
| Ethyl acetate | 51.0 |
| Water | 0.154 |

HANDLING

m-Chloroperoxybenzoic acid irritates the mucous membranes and respiratory tract, eyes and skin. Skin contact with MCPBA results in burns and blisters similar to those cause by hydrogen peroxide. the material should be used only in a chemical fume hood. Safety goggles, rubber gloves, and a dust mask should be worn.

EMERGENCY PROCEDURES

FIRE

Extinguish with "alcohol" foam, dry powder, or carbon dioxide. Water may be ineffective. Fire conditions may cause explosions. Wear a self-contained breathing apparatus.

SPILL

Wear safety goggles and rubber gloves. Treat with excess sodium bisulfite solution. Test with acidic starch-iodide paper to ensure complete decomposition of the peracid. wash down the drain.

SKIN CONTACT

Wash immediately with soap and water or sodium bicarbonate solution. Remove contaminated clothing.

EYE CONTACT

Flush immediately with large amounts of water and call a physician.

INGESTION

Take large quantities of milk or water immediately. Wash out mouth. Call a physician.

WASTE DISPOSAL

Treat with excess sodium bisulfite solution. Test with acidic starch-iodide paper to ensure complete decomposition. Dispose of properly. Observe all state, local and federal laws.

STORAGE AND STABILITY

Solid MCPBA shows less than 1% decomposition when stored at room temperature for 1 year. It has been determined that 95-100% metarial can be detonated by shock or sparks. 85% MCPBA is not shock-sensitive, but may be decomposed violently with heat. It should be stroed in a refrigerator in tightly closed containers. MCPBA is a flammable solid and contact with heat or oxidizable material should be avoided.

Questions? or for ordering information call 800-558-9160 (USA/Canada) or 414-273-3850

Aldrich Chemical Company, Inc.

1001 West Saint Paul Ave., Milwaukee, WI 53233

Telephone 414-273-3850

800-231-8327

Fax 414-273-4979

800-962-9591

Internet aldrach@sial.com

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.

