

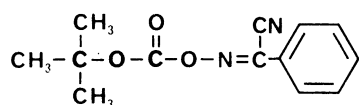
[2-(*tert*-butoxycarbonyloxyimino)-2-phenylacetoneitrile]

revised 4 / 96

4 pages

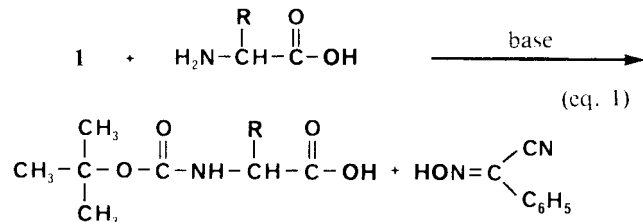
PROPERTIES

Molecular formula	C ₁₃ H ₁₄ N ₂ O ₃
Molecular weight	246.27
Melting point	87-89°C
Solubility	Very soluble in ethyl acetate, ether, benzene, chloroform, dioxane and acetone. Soluble in methanol, 2-propanol and <i>tert</i> -butanol. Insoluble in petroleum ether and water.



APPLICATION

BOC-ON, **1**, reacts simple and rapidly with amino acids to provide protected derivatives in excellent yields (eq.1). *t*-Butoxy-carbonylation of amino acids is carried out at room temperature using a 10% excess of BOX-ON and a 50% excess of triethylamine in either 50% aqueous dioxane or 50% aqueous acetone.



This reagent offers a distinct advantage over *t*-BOC azide which can require reaction temperatures of 50-60°C (*t*-BOC azide is thermally unstable and decomposes with apparent detonation at temperatures above 80°C).¹ The oxime by-product can be easily and completely removed from the reaction mixture by extraction with ether, ethyl acetate or benzene. Deprotection of *t*-BOC amino acids is accomplished under mildly acidic conditions.

The bibliography on page 2 lists references to the use of BOC-ON since its introduction in 1975 by Itoh *et al.*² A number of patents to its use in pharmaceutical and agricultural applications have been issued but are not included in this bibliography.

GENERAL PROCEDURE FOR THE INTRODUCTION OF THE *t*-BOX GROUP INTO AMINO ACIDS³

To a stirred solution of L-tryptophan (2.04g, 10mmol) and triethylamine (2.10ml, 15mmol) (Note 1) is added BOC-ON (2.71g, 11mmol) in a mixture of water (6ml) and dioxane (6ml) at room temperature. The mixture is homogeneous within 1 hour and stirring is continued for 2 hours (Note 2). After addition of water (15ml) and ethyl acetate (20ml), the aqueous layer is separated, washed with ethyl acetate (20ml) (Note 3), acidified with 5%

citric acid solution (Note 4) and extracted with ethyl acetate. The dry, concentrated extract gave *t*-butoxycarbonyl-L-tryptophan, 3.00g (98.6%), mp 137-138°C (dec.).

Examples of other *t*-BOC-amino acids prepared by this procedure are shown in Table 1 (Note 5).

TABLE 1

<i>t</i> -BOX-Amino acids	Reaction time (h)	Oxime extraction solvent	Yield (%)	m.p. (°C)
-Gly-OH	2	EtOAc	86.9	86-88
-Leu-OH•½H ₂ O	3	EtOAc	72.0	78-84
-Met-OH•DCHA*	3	EtOAc	82.1	137-139
-Thr-OH•DCHA	3	EtOAc	99.7	151-153
-Phe-OH•DCHA	5	Ether	98.2	221-223

* = dicyclohexylamine

STABILITY AND STORAGE

BOC-ON should be stored in a brown bottle in the freezer (-20°C) to prolong its shelf life. After several weeks at room temperature, BOC-ON undergoes gradual decomposition with evolution of carbon dioxide.

TOXICITY AND HANDLING

Contact with the product may cause irritation. To the best of our knowledge, the chemical, physical and toxicological properties have not been thoroughly investigated. The usual safe laboratory precautions and procedures should be observed when handling this material.

NOTES

- 1• Sodium hydroxide, sodium bicarbonate and other inorganic bases (10 to 15mmol) may be used in place of triethylamine as long as the amino acids are solubilized in the system. In such cases, excess base over 10mmol should be neutralized before the extraction of 2-hydroxyimino-2-phenylacetoneitrile to prevent contamination of the *t*-BOC-amino acids.
- 2• When using a solvent other than dioxane, it is recommended that the solvent be removed before extraction to ensure high yield and purity of the *t*-BOX-amino acid derivative.
- 3• Other suitable solvents for the extraction of 2-hydroxyimino-2-phenylacetoneitrile are benzene and ether. For amino acids with lipophilic side chains, the use of ether is recommended in order to obtain contaminant-free products in high yields.
- 4• Alternatively, cold 1N hydrochloric acid may be used.
- 5• Dicyclohexylammonium salts are crystallized by the addition of dicyclohexylamine (1.8g per 10mmol amino acid) to an ether solution of *t*-butoxycarbonylamino acids.³

REFERENCES

- 1• *Org. Syn.* **1977**, *57*, 122.
- 2• Itoh, M.; Hagiware, D.; Kamiya, T. *Tetrahedron Lett.* **1975**, 4393.
- 3• *Idem Bull. Chem. Soc. Jpn.* **1977**, **50**, 718.

BIBLIOGRAPHY

- 1• Acton, N.; Komoriya, A. Synthesis of Pseudopeptides. *Org. Prep. Proced. Int.* **1982**, *14*, 381.
- 2• Aimoto, S.; Richards, F. Synthesis of Carriers of differing Stokes Radius with Activated Acyl groups for Use as Reagents in Labeling Membrane Proteins. *J. Biol. Chem.* **1981**, *256*, 5134.
- 3• Andok, S.; Aoyagi, H.; Waki, M.; Kato, T.; Izumiya, N. Studies of Peptide Antibiotics. XLIII. Synthesis of Gramicidin S Analogs Containing D-Serine or Dehydroalanine in Place of D-Phenylalanine and Asymmetric Hydrogenation of the Dehydroalanine Residue. *Int. J. Pept. Protein Res.* **1983**, *21*, 313.
- 4• Anonymous. *t*-Butyl Azidoformate. *Org. Syn.* **1977**, *57*, 122.
- 5• Armstrong, D.W.; Seguin, R.; Saburi, M.; Fendler, J.H. Synthesis of Amino Acyl Adenylates Using the *tert*-Butoxycarbonyl Protecting Group. *J. Mol. Evol.* **1979**, *13*, 103.
- 6• Baldwin, J.E.; Otsuka, M.; Wallace, P.M. Synthesis of a Naturally Occurring Inhibitor of Glutamine Synthetase, Tabtoxinine- β -lactam. *Chem. Commun.* **1985**, 1549.
- 7• Bartlett, P.A.; Meadows, J.D.; Brown, E.G.; Morimoto, A.; Jernstedt, K.K. Carbonate Extension. A Versatile Procedure for Functionalization of Acyclic Homoallylic Alcohols with Moderate Stereocontrol. *J. Org. Chem.* **1982**, *47*, 4013.
- 8• Bartlett, P.A.; Meadows, J.D.; Ottow, E. Enantiodivergent Syntheses of (+)- and (-)- Nonactic Acid and the Total Synthesis of Nonactin. *J. Am. Chem. Soc.* **1984**, *47*, 4013.
- 9• Beamer, R.L.; Griffith, O.W.; Gass, J.D.; Anderson, M.E.; Meister, A. Interaction of L- and D-3-Amino-1-chloro-2-pentanone with γ -Glutamylcysteine Synthetase. *J. Biol. Chem.* **1980**, 689.
- 10• Bergeron, R.J.; Stolorow, J.J.; Porter, C.W. Reagents for the Selective Secondary N-Acylation of Linear Triamines. *Synthesis* **1982**, 689.
- 11• Bodanszky, M.; Bednarek, M.A. Derivatives of S-9-Fluorenylmethyl-L-cysteine. *Int. J. Pept. Protein Res.* **1982**, *20*, 434.
- 12• Bodanszky, M.; Martinez, J.; Priestley, G.P.; Gardner, J.D.; Mutt, V. Cholecystokinin (Pancreozymin). 4) Synthesis and Properties of a Biologically Active Analog of the C-Terminal Heptapeptide with ϵ -Hydroxynorleucine Sulfate Replacing Tyrosine Sulfate. *J. Med. Chem.* **1978**, *21*, 1030.
- 13• Cachia, P.J.; Sykes, B.D.; Hodges, R.S. Calcium-dependant Inhibitory Region of Troponin: A protein Nuclear Magnetic Resonance Study of the Interaction between Troponin C and the Synthetic Peptide N α -Acetyl-[Fphe¹⁰⁶]TnI-(104-115) Amide. *Biochemistry* **1983**, *22*, 4145.
- 14• Fisher, G.H.; Berryer, P.; Ryan, J.W.; Chauhan, V.; Stammer, C.H. Dehydrophenylalanyl Analogs of Bradykinin: Synthesis and Biological Activities. *Arch. Biochem. Biophys.* **1981**, *211*, 269.
- 15• Galaray, R.E. Inhibition of Angiotensin-Converting Enzyme by Phosphoramidates and Polyphosphates. *Biochemistry* **1982**, *21*, 5777.
- 16• Gigot, D.; Penninckx, M. New Compounds: Peptide Derivatives of the Antitumor Agent N-Phosphonoacetyl-L-aspartic Acid. *J. Pharm. Sci.* **1984**, *73*, 275.
- 17• Glass, D.; Dembure, P.; Priest, J.; Elsas, L., II. A [³H]Lysine-containing Synthetic Peptide Substrate for Human Protocollagen Lysyl Hydroxylase. *Biochim. Biophys. Acta* **1985**, *840*, 143.
- 18• Heimer, E.P.; Wang, C.T.; Lambros, T.J.; Felix, A.M. Direct Synthesis of N α -Benzoyloxycarbonyl-N ϵ -*tert*-butyloxycarbonyl-L-lysine from L-Lysine. *Org. Prep. Proced. Int.* **1983**, *15*, 379.
- 19• Horiuchi, Y.; Ikeda, D.; Gomi, S.; Kondo, S.; Umezawa, H. Isolation and Synthesis of 2"-N-Formimidoylistamycins A and B, New Istamycin Components. *Carbohydr. Res.* **1982**, *109*, 25.
- 20• Hunt, A.; Dorman, D.; Debono, M.; Molloy, R.M. Structure of Antibiotic A41030AA. *J. Org. Chem.* **1985**, *50*, 2031.
- 21• Inman, J.K.; DuBois, G.C.; Appella, E. Synthesis of Mesoporous Polystyrene Derivatives for Supports in Sequential Degradation of Peptides. *INSERM Symp.* **1977**, *5* (Solid Phase Methods Protein Sequence Anal.), 81.
- 22• Itoh, M.; Hagiware, D.; Kamiya, T. A New Reagent for *tert*-Butoxycarbonylation: 2-*tert*-Butoxycarbonyloxymino-2-phenylacetonitrile. *Org. Syn.* **1980**, *59*, 95.
- 23• Itoh, M.; Hagiware, D.; Kamiya, T. Peptides. VI. Some Oxime Carbonates as Novel *tert*-Butoxycarbonylating Reagents. *Bull. Chem. Soc. Jpn.* **1977**, *50*, 718.
- 24• Itoh, M.; Hagiware, D.; Kamiya, T. *Tetrahedron Lett.* **1975**, 4393.
- 25• Johnson, R.L. Renin Inhibitors. Substitution of the Leucyl residues of Leu-Leu-Val-Phe-OCH₃ with 3-Amino-2-hydroxy-5-methylhexanoic Acid. *J. Med. Chem.* **1982**, *25*, 605.
- 26• Komai, L.; Matsuyama, K. Synthesis and Thermal Decomposition of Di-*tert*-butylperoxydicarbonate. *Bull. Chem. Soc. Jpn.* **1985**, *58*, 2207.
- 27• Lee, T.T.; Williams, R.E.; Fox, C.F. Photoaffinity Inactivation of the Enkephalin Receptor. *J. Biol. Chem.* **1979**, *254*, 11787.
- 28• Matsuda, K.; Yasuda, J.; Tsutsumi, H.; Takaya, T. Studies on Antiviral Activity on New Kanamycin A Derivatives Having Higher Acyl Group at N-1 Position. *J. Antibiot.* **1985**, *38*, 1050.
- 29• Matsuda, K.; Yasuda, N.; Tsutsumi, H.; Takaya, T. Studies on Antiviral Agents. III. Synthesis and *in vitro* Antiviral Activity of 1-N-Higher-acyl-3"-N-functionalized Acylkanamycin A Derivatives. *J. Antibiot.* **1985**, *38*, 1719.

BIBLIOGRAPHY (CONTINUED)

- 30• Mitra, A.K.; Ostashevsky, I.; Brewer, C.F. Synthesis and Fluorine-19 Spectra of Tetra-L-alanine Residues. *Int. J. Pept. Protein Res.* **1983**, *22*, 495.
- 31• Nakajima, K.; Tanada, T.; Neya, M.; Okawa, K. Total Synthesis of Actinomycin D and Its Analog. *Pept. Chem.* **1981**, *19*, 143.
- 32• Nicholson, A.W.; Cooperman, B.S. Photoaffinity Labeling of *Escherichia coli* Ribosomes with an Aryl Azide Analog of Puromycin. *FEBS Lett.* **1978**, *90*, 203.
- 33• Nomoto, S.; Shiba, T. Synthesis of N β -Methyl-L-arginine, a Component Amino Acid in a New Antibiotic, LL-BM547 β . *Chem. Lett.* **1978**, 589.
- 34• Nutt, R.F.; Jouillie, M. Four-Component Condensation: a New Versatile Method for the Synthesis of Substituted Prolyl Peptides. *J. Am. Chem. Soc.* **1982**, *104*, 5852.
- 35• Nutt, R.F.; Veber, D.F.; Curley, P.E.; Saperstein, R.; Hirshchmann, R. Somatostatin Analogs Which Define the Role of the Lysine-9 Amino Group. *Int. J. Pept. Protein Res.* **1983**, *21*, 66.
- 36• Ohfuné, Y.; Tomita, M.; Nomoto, K. Total Synthesis of 2'-Deoxymugineic Acid, the Metal Chelator Excreted from Wheat Root. *J. Am. Chem. Soc.* **1981**, *103*, 2409.
- 37• Otsuka, M.; Kittaka, A.; Iimorei, T.; Yamashita, H.; Kobayashi, S.; Ohno, M.; Synthetic Studies on an Antitumor Antibiotic, Bleomycin. XII. Preparation of an L-2,3-Diaminopropionic Acid Derivative as a Synthetic Intermediate. *Chem. Pharm. Bull.* **1985**, *33*, 509.
- 38• Platen, M.; Steckhan, E. Oxidative Deblocking of the 4-Methoxybenzyl Thioether Protecting Group: Application to the Directed Synthesis of Polycystinyl Peptides. *Liebigs Ann. Chem.* **1984**, 1563.
- 39• Prochazka, Z.; Jost, K. Amino Acids and Peptides. CLXIII. L-Cystathionine Derivatives for the Synthesis of Peptides. *Collect. Czech. Chem. Commun.* **1980**, *45*, 1982.
- 40• Rao, S.P.; Dunn, B.M. Preparation of Photoaffinity Labels of Pepsin and *p*-Nitro-, *p*-Azido- and *p*-Diazophenyl Ligands and a study of the Effects of Irradiation of Pepsin. *Biochim. Biophys. Acta* **1982**, *706*, 86.
- 41• Rosowsky, A.; Freisheim, J.; Moran, R.; Solan, V.; Bader, H.; Wright, J.; Radike-Smith, M. Methotrexate Analogs. 26. Inhibition of Dihydrofolate Reductase and Folylpolylglutamate Synthetase Activity and *in vitro* Tumor Cell Growth by Methotrexate and Aminopterin Analogs Containing a Basic Amino Acid Side Chain. *J. Med. Chem.* **1986**, *29*, 655.
- 42• Rosowsky, A.; Wright, J.E. N ω -Alkoxyacylation of α,ω -Diamino Acids with 2-(Trimethylsilyl)ethyl 4-Nitro-phenyl Carbonate. *J. Org. Chem.* **1983**, *48*, 1539.
- 43• Shashoua, V.E.; Jacob, J.N.; Ridge, R.; Campbell, A.; Baldessarini, R. γ -Aminobutyric Acid Esters. I. Synthesis, Brain Uptake and Pharmacological studies of Aliphatic and Steroid Esters of γ -Aminobutyric Acid. *J. Med. Chem.* **1984**, *27*, 659.
- 44• Shoelson, S.; Fickova, M.; Haneda, M.; Nahum, A.; Musso, G.; Kaiser, E.T.; Rubenstein, A.H.; Tager, H. Identification of a Mutant Human Insulin Predicted to Contain a Serine-for-Phenylalanine Substitution. *Proc. Nat. Acad. Sci. USA* **1983**, *80*, 7390.
- 45• Staros, J.V.; Knowles, J.R. Photoaffinity Inhibition of Dipeptide Transport in *Escherichia coli*. *Biochemistry* **1978**, *17*, 3321.
- 46• Stewart, F.H.C. Synthesis of L-Kynurenine Peptides Conducted without Masking the Side-chain Amino Group. *Aust. J. Chem.* **1980**, *33*, 633.
- 47• Talma, A.; Jouin, P.; DeVries, J.; Troostwijk, C.; Bunine, G.; Waninge, J.; Visscher, J.; Kellogg, R. Reductions of Activated Carbonyl Compounds with Chiral-Bridged 1,4-Dihydropyridines. An Investigation of Scope and Structural Effects. *J. Am. Chem. Soc.* **1985**, *107*, 3981.
- 48• Takada, S.; Ishizuka, N.; Sasatani, T.; Makisumi, Y.; Jyoyama, H.; Hatakeyama, H.; Asanuma, F.; Hirose, K. Studies on Fused Indoles. II. Structural Modifications and Analgesic Activity of 4-Aminomethyltetrahydrothiopyrano-[2,3-*b*]indoles. *Chem. Pharm. Bull.* **1984**, *32*, 877.
- 49• Tsunematsu, H.; Imamura, T.; Makisumi, S. Kinetics of Hydrolysis of N α -Benzoyl-*p*-guanidino-L-phenylalanine *p*-Nitroanilide by Trypsin. *J. Biochem. (Tokyo)* **1983**, *94*, 123.
- 50• Van den Braken-Van Leersum, A.M.; Maat, L. Synthesis of [L-*m*-Tyr¹]- and [D-*m*-Tyr¹]-leucine-enkephalin *via* the REMA Method. *J.R. Neth. Chem. Soc.* **1984**, *103*, 110.
- 51• Venn, R.F.; Barnard, E.A. A Potent Peptide Affinity Reagent for the Opiate Receptor. *J. Bio. Chem.* **1981**, *256*, 1529.
- 52• Waki, M.; Kitajima, Y.; Isumiya, N. A Facile Synthesis of N²-Protected L-2,3-Diaminopropanoic Acid. *Synthesis* **1981**, 266.
- 53• Wakimasu, M.; Kitada, C.; Fujino, M. 4-Methoxy-2,3,6-tri-methylbenzenesulfonyl (Mtr): A New Amino and Imidazole Protecting Group in Peptide Synthesis. *Chem. Pharm. Bull.* **1982**, *30*, 2766.
- 54• Zemlicka, J.; Bhuta, A.; Bhuta, P. Analogs of 2'(3')-O-L Phenylalanyladenosine as Substrates and Inhibitors of Ribosomal Peptidyltransferase. *J. Med. Chem.* **1983**, *26*, 167.

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.

