

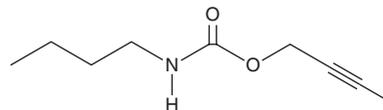
# PRODUCT INFORMATION



## 3-Iodo-2-propynyl Butylcarbamate

Item No. 36028

**CAS Registry No.:** 55406-53-6  
**Formal Name:** N-butyl-carbamic acid, 3-iodo-2-propyn-1-yl ester  
**Synonyms:** Iodopropynyl Butylcarbamate, IPBC  
**MF:** C<sub>8</sub>H<sub>12</sub>INO<sub>2</sub>  
**FW:** 281.1  
**Purity:** ≥95%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

3-Iodo-2-propynyl Butylcarbamate (IPBC) is supplied as a solid. A stock solution may be made by dissolving the IPBC in the solvent of choice, which should be purged with an inert gas. IPBC is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of IPBC in these solvents is approximately 2, 15, and 10 mg/ml, respectively.

### Description

IPBC is a carbamate fungicide and an algicide.<sup>1,2</sup> It inhibits the growth of *A. niger* when applied to linear low-density polyethylene at concentrations ranging from 1,000 to 50,000 ppm.<sup>1</sup> IPBC is active against dairy cow isolates of *P. zopfii*, *P. blaschkeae*, and *Prototheca* parasitic algae (MIC<sub>50S</sub> = 4, 1, and 4 mg/L, respectively).<sup>2</sup> Formulations containing IPBC have been used as antifungal agents in cosmetic and industrial applications.

### References

1. Gitchaiwat, A., Kositchaiyong, A., Sombatsompop, K., *et al.* Assessment and characterization of antifungal and antialgal performances for biocide-enhanced linear low-density polyethylene. *J. Appl. Polym. Sci.* **128(1)**, 371-379 (2013).
2. Jagielski, T., Bakuła, Z., Di Mauro, S., *et al.* A comparative study of the *in vitro* activity of iodopropynyl butylcarbamate and amphotericin B against *Prototheca* spp. isolates from European dairy herds. *J. Dairy Sci.* **100(9)**, 7435-7445 (2017).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/19/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897

[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM