

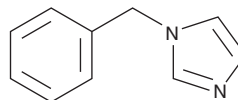
# PRODUCT INFORMATION



## 1-Benzylimidazole

Item No. 70510

**CAS Registry No.:** 4238-71-5  
**Formal Name:** 1-(phenylmethyl)-1H-imidazole  
**MF:** C<sub>10</sub>H<sub>10</sub>N<sub>2</sub>  
**FW:** 158.2  
**Purity:** ≥95%  
**Supplied as:** A crystalline 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

1-Benzylimidazole is supplied as a crystalline solid. A stock solution may be made by dissolving the 1-benzylimidazole in the solvent of choice, which should be purged with an inert gas. 1-Benzylimidazole is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 1-benzylimidazole in these solvents is approximately 500, 560, and 530 µg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of 1-benzylimidazole can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 1-benzylimidazole in PBS, pH 7.2, is approximately 6.5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

1-Benzylimidazole is a selective inhibitor of thromboxane synthase. It reduces TXB<sub>2</sub> levels and increases blood flow in ischemia-reperfusion injury of rat brain.<sup>1</sup>

### Reference

1. Pettigrew, L.C., Grotta, J.C., Rhoades, H.M., *et al.* Effect of thromboxane synthase inhibition on eicosanoid levels and blood flow in ischemic rat brain. *Stroke* **20(5)**, 627-632 (1989).

#### 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, 10/19/2022

#### 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