

PRODUCT INFORMATION



Bromelain

Item No. 17337

CAS Registry No.: 37189-34-7

Supplied as: A powder

Storage: -20°C

Stability: 4 years

Specific Activity: 0.5 DMC-U/mg

Unit Definition: 1 DMC-U catalyzes the cleavage of 1 μ mole peptide bond from dimethyl casein per minute at 25°C, pH 7.0 expressed in terms of newly formed terminal amino groups (determined with TNBS).

Laboratory Procedures

Bromelain is supplied as a powder. Aqueous solutions of bromelain can be prepared by directly dissolving the powder in aqueous buffers. The solubility of bromelain in 50 mM NaH_2PO_4 (pH 7.0) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Bromelain is a mixture of proteolytic enzymes and non-protease components (phosphatases, glucosidases, peroxidases, cellulases, glycoproteins, and carbohydrates) derived from the stem and fruit of pineapple plants.^{1,2} Bromelain has been studied for a number of potentially beneficial effects, including anti-inflammatory, anti-thrombotic, fibrinolytic, and anti-cancer functions.^{1,2}

References

1. Chobotova, K., Vernallis, A.B., and Majid, F.A.A. Bromelain's activity and potential as an anti-cancer agent: Current evidence and perspectives. *Cancer Lett.* **290(2)**, 148-156 (2010).
2. Yuan, G., Wahlqvist, M.L., He, G., *et al.* Natural products and anti-inflammatory activity. *Asia Pac. J. Clin. Nutr.* **15(2)**, 143-152 (2006).

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, 10/28/2022

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