

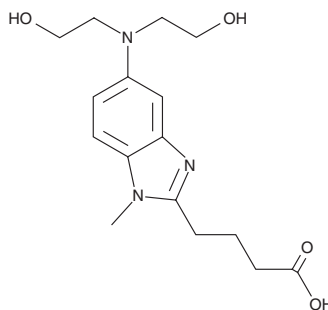
PRODUCT INFORMATION



Dihydroxy Bendamustine

Item No. 46211

CAS Registry No.: 109882-30-6
Formal Name: 5-[bis(2-hydroxyethyl)amino]-1-methyl-1H-benzimidazole-2-butanoic acid
Synonym: HP2
MF: C₁₆H₂₃N₃O₄
FW: 321.4
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

Dihydroxy bendamustine is supplied as a solid. A stock solution may be made by dissolving the dihydroxy bendamustine in the solvent of choice, which should be purged with an inert gas. Dihydroxy bendamustine is sparingly soluble (1-10 mg/ml) in DMSO.

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 dihydroxy bendamustine can be prepared by directly dissolving the solid in aqueous buffers. Dihydroxy bendamustine is sparingly soluble in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Dihydroxy bendamustine is an inactive metabolite of the DNA alkylating agent bendamustine (Item No. 23693).^{1,2}

References

1. Iwamoto, K., Uehara, Y., Inoue, Y., *et al.* Inhibition of STAT3 by anticancer drug bendamustine. *PLoS One* **12**(1), e0170709 (2017).
2. Dubbelman, A.C., Tibben, M., Rosing, H., *et al.* Development and validation of LC-MS/MS assays for the quantification of bendamustine and its metabolites in human plasma and urine. *J. Chromatogr. B Analyt. Technol. Biomed. Life Sci.* **893-894**, 92-100 (2012).

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, 08/27/2026

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