

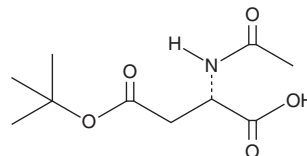
PRODUCT INFORMATION



Ac-Asp(OtBu)-OH

Item No. 30539

CAS Registry No.: 117833-18-8
Formal Name: N-acetyl-L-aspartic acid, 4-(1,1-dimethylethyl) ester
MF: C₁₀H₁₇NO₅
FW: 231.2
Purity: ≥95%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Ac-Asp(OtBu)-OH is supplied as a crystalline solid. A stock solution may be made by dissolving the Ac-Asp(OtBu)-OH in the solvent of choice, which should be purged with an inert gas. Ac-Asp(OtBu)-OH is soluble in the organic solvent DMSO at a concentration of approximately 1 mg/ml.

Ac-Asp(OtBu)-OH is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, Ac-Asp(OtBu)-OH should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Ac-Asp(OtBu)-OH has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Ac-Asp(OtBu)-OH is an amino acid-containing building block.¹⁻³ It has been used in the synthesis of caspase-3 and -7 peptide substrates and inhibitors, as well as peptide inhibitors of HIV integrase.

References

1. Peterson, Q.P., Goode, D.R., West, D.C., *et al.* Preparation of the caspase-3/7 substrate Ac-DEVD-pNA by solution-phase peptide synthesis. *Nat. Protoc.* **5**(2), 294-302 (2010).
2. Goode, D.R., Sharma, A.K., and Hergenrother, P.J. Using peptidic inhibitors to systematically probe the S1' site of caspase-3 and caspase-7. *Org. Lett.* **7**(16), 3529-3532 (2005).
3. Gordon, C.P., Dalton, N., Vandergraaff, N., *et al.* A structure-based design approach to advance the allyltyrosine-based series of HIV integrase inhibitors. *Tetrahedron* **74**(12), 1253-1268 (2018).

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, 05/04/2020

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