

PRODUCT INFORMATION



UAMC 0039 (hydrochloride)

Item No. 15340

CAS Registry No.: 697797-51-6

Formal Name: (2S)-2-amino-4-[[[(4-chlorophenyl)methyl]amino]-1-(1-piperidiny)]-1-butanone, dihydrochloride

MF: $C_{16}H_{24}ClN_3O \cdot 2HCl$

FW: 382.8

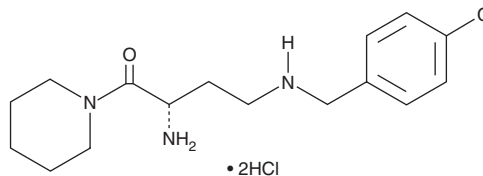
Purity: $\geq 98\%$

UV/Vis.: λ_{max} : 220 nm

Supplied as: A crystalline solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



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

Laboratory Procedures

UAMC 0039 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the UAMC 0039 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. UAMC 0039 (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of UAMC 0039 (hydrochloride) in these solvents is approximately 5, 14, and 2 mg/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 UAMC 0039 (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of UAMC 0039 (hydrochloride) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

UAMC 0039 is a potent inhibitor of dipeptidyl peptidase 2 (DPP-2; $IC_{50} = 0.48$ nM).¹ It also inhibits DPP-4 ($IC_{50} = 165$ μ M). DPP-2 is thought to be involved in degradation of collagen, substance P, and some proline-containing neuropeptides.¹⁻³

References

1. Senten, K., Van der Veken, P., De Meester, I., et al. γ -amino-substituted analogues of 1-[(S)-2,4-diaminobutanoyl]piperidine as highly potent and selective dipeptidyl peptidase II inhibitors. *J. Med. Chem.* **47**(11), 2906-2916 (2004).
2. Andersen, K.J., and McDonald, J.K. Lysosomal heterogeneity of dipeptidyl peptidase II active on collagen-related peptides. *Ren. Physiol. Biochem.* **12**(1), 32-40 (1989).
3. Mentlein, R., and Struckhoff, G. Purification of two dipeptidyl aminopeptidases II from rat brain and their action on proline-containing neuropeptides. *J. Neurochem.* **52**(4), 1284-1293 (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, 11/08/2022

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