

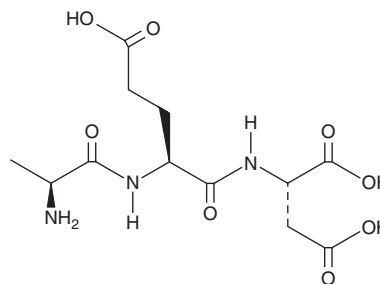
PRODUCT INFORMATION



H-Ala-Glu-Asp-OH

Item No. 45082

CAS Registry No.: 85806-95-7
Formal Name: L-alanyl-L- α -glutamyl-L-aspartic acid
Synonyms: AED, T-31
Peptide Sequence: H-AED-OH
MF: C₁₂H₁₉N₃O₈
FW: 333.3
Purity: \geq 98%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

H-Ala-Glu-Asp-OH (AED) is supplied as a solid. A stock solution may be made by dissolving the AED in the solvent of choice, which should be purged with an inert gas. AED 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 AED can be prepared by directly dissolving the solid in aqueous buffers. AED is soluble (\geq 10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

AED is a synthetic tripeptide with geroprotective effects *in vitro*.^{1,2} It increases proliferation and reduces apoptosis in young and aged rat skin fibroblast cultures but reduces matrix metalloproteinase-9 (MMP-9) levels only in young cultures.¹ AED (20 ng/ml) increases sirtuin 1 (SIRT1), SIRT6, and collagen I levels in young and aged human skin fibroblast cultures.² It also prevents mortality in a mouse model of cisplatin-induced acute kidney injury (AKI) when administered at a dose of 3 μ g/kg.³

References

1. Lin'kova, N.S., Drobintseva, A.O., Orlova, O.A., *et al.* Peptide regulation of skin fibroblast functions during their aging *in vitro*. *Bull. Exp. Biol. Med.* **161**(1), 175-178 (2016).
2. Fridman, N.V., Linkova, N.S., Kozhevnikova, E.O., *et al.* Comparison of the effects of KE and AED peptides on functional activity of human skin fibroblasts during their replicative aging. *Bull. Exp. Biol. Med.* **170**(1), 154-157 (2020).
3. Shchudrova, T.S., Zamorskii, I.I., and Korotun, O.P. Evaluation of a renoprotective potential of organospecific peptides under the conditions of acute kidney injury of different etiology. *J. Sport Health Sci.* **8**(1), 221-227 (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, 05/05/2026

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