

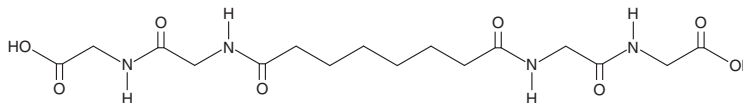
PRODUCT INFORMATION



CAY10679

Item No. 9001475

CAS Registry No.: 143673-93-2
Formal Name: 1,1'-(1,8-dioxo-1,8-octanediyl)
bis[glycyl-glycine
Synonym: N,N'-bis(glycyl-glycine)-hexane-
1,6-Dicarboxamide
MF: C₁₆H₂₆N₄O₈
FW: 402.4
Purity: ≥95%
Supplied as: A crystalline 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

CAY10679 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10679 in the solvent of choice, which should be purged with an inert gas. CAY10679 is soluble in the organic solvent DMSO (warmed) at a concentration of approximately 1 mg/ml.

Description

CAY10679 is an anionic oligopeptide-based bola-amphiphile that contains a hexane moiety flanked by a glycylglycine group at each end, which creates distinct hydrophilic and hydrophobic regions.¹ The microtube self-assembling properties of this amphiphilic peptide have been studied in aqueous acidic solution.² Bola-amphiphiles are of interest in the construction of nanotubes, nanovesicles, nanowires, nanocarriers for drug delivery, and other applications of bionanotechnology.¹

References

1. Santoso, S.S., Vauthey, S., and Zhang, S. Structures, function and applications of amphiphilic peptides. *Curr. Opin. Colloid In.* **7**, 262-266 (2013).
2. Naidoo, V.B., and Rautenbach, M. Bidirectional solid phase synthesis of a model oligoglycine bolaamphiphile and purification by rapid self-assembly. *J. Pept. Sci.* **18(5)**, 317-325 (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, 04/29/2024

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