

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

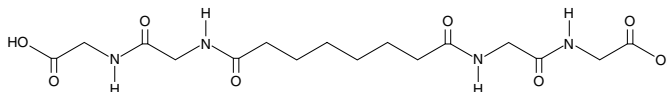
MF: C₁₆H₂₆N₄O₈
FW: 402.4

Purity: ≥95%

Stability: ≥2 years at -20°C

Supplied as: A crystalline solid

Special Conditions: DMSO should be warmed to solubilize



Laboratory Procedures

For long term storage, we suggest that CAY10679 be stored as supplied at -20°C. It should be stable for at least two years.

CAY10679 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10679 in the solvent of choice. CAY10679 is soluble in DMSO at a concentration of approximately 1 mg/ml.

CAY10679 is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

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).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/9001475

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 12/20/2013

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com