

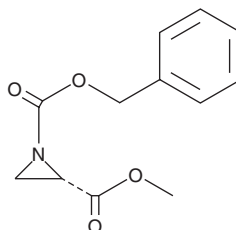
# PRODUCT INFORMATION



## Methyl (S)-(-)-N-Z-Aziridine-2-carboxylate

Item No. 20731

**CAS Registry No.:** 104597-98-0  
**Formal Name:** (2S)-1,2-aziridinedicarboxylic acid, 2-methyl 1-(phenylmethyl) ester  
**MF:** C<sub>12</sub>H<sub>13</sub>NO<sub>4</sub>  
**FW:** 235.2  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 258 nm  
**Supplied as:** A neat oil  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

Methyl (S)-(-)-N-Z-aziridine-2-carboxylate is supplied as a neat oil. A stock solution may be made by dissolving the methyl (S)-(-)-N-Z-aziridine-2-carboxylate in the solvent of choice, which should be purged with an inert gas. Methyl (S)-(-)-N-Z-aziridine-2-carboxylate is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of methyl (S)-(-)-N-Z-aziridine-2-carboxylate in ethanol and DMSO is approximately 20 mg/ml and approximately 30 mg/ml in DMF.

Methyl (S)-(-)-N-Z-aziridine-2-carboxylate is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, methyl (S)-(-)-N-Z-aziridine-2-carboxylate should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Methyl (S)-(-)-N-Z-aziridine-2-carboxylate has a solubility of approximately 0.125 mg/ml in a 1:7 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Methyl (S)-(-)-N-Z-aziridine-2-carboxylate is a chiral building block that is commonly used in the organic stereoselective synthesis of various compounds, including anticancer agents, antibiotics, and enzyme inhibitors.<sup>1,2</sup>

### References

- Li, Y., Hayman, E., Plesescu, M., *et al.* Synthesis of potent BCRP inhibitor - Kol143. *Tetrahedron Lett.* **49(9)**, 1480-1483 (2008).
- Bartocchini, F., Venturi, S., Retini, M., *et al.* Total synthesis of (-)-clavicipitic acid via γ,γ-dimethylallyltryptophan (DMAT) and chemoselective C-H hydroxylation. *J. Org. Chem.* **84(12)**, 8027-8034 (2019).

#### 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, 11/12/2022

#### 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