

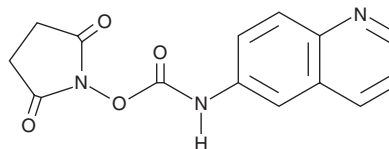
PRODUCT INFORMATION



AQC

Item No. 30877

CAS Registry No.: 148757-94-2
Formal Name: N-6-quinolinyl-carbamic acid, 2,5-dioxo-1-pyrrolidinyl ester
Synonym: 6-Aminoquinolyl-N-hydroxysuccinimidyl Carbamate
MF: $C_{14}H_{11}N_3O_4$
FW: 285.3
Purity: $\geq 95\%$
UV/Vis.: λ_{max} : 211, 242 nm
Ex./Em. Max: 248/398 nm, respectively, for derivatized polyamines
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Special Conditions: Hygroscopic



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

Laboratory Procedures

AQC is supplied as a solid. A stock solution may be made by dissolving the AQC in the solvent of choice, which should be purged with an inert gas. AQC is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of AQC in these solvents is approximately 5 and 1 mg/ml, respectively.

AQC is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AQC should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. AQC has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

AQC is a reactive probe for the pre-column derivatization of primary and secondary amines, including those found in amino acids, peptides, proteins, and polyamines.^{1,2} It forms stable and highly fluorescent derivatives and has been used for chromatographic separation and analysis of derivatized amino acids and polyamines.¹⁻³ AQC-derivatized polyamines display excitation/emission maxima of 248/398 nm, respectively.²

References

1. Cohen, S.A. Quantitation of amino acids as 6-aminoquinolyl-N-hydroxysuccinimidyl carbamate derivatives. *J. Chromatogr. Lib.* **70**, 242-267 (2005).
2. Weiss, T.S. HPLC of biogenic amines as 6-aminoquinolyl-N-hydroxysuccinimidyl derivatives. *J. Chromatogr. Lib.* **70**, 502-523 (2005).
3. Pawlowska, M., Chen, S., and Armstrong, D.W. Enantiomeric separation of fluorescent, 6-aminoquinolyl-N-hydroxysuccinimidyl carbamate, tagged amino acids. *J. Chromatogr. A* **641**(2), 257-265 (1993).

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, 07/25/2023

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