

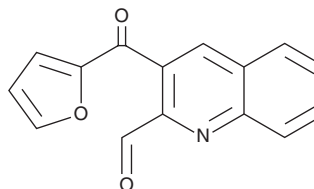
PRODUCT INFORMATION



3-(2-Furoyl)quinoline-2-carboxaldehyde

Item No. 44747

CAS Registry No.: 126769-01-5
Formal Name: 3-(2-furanylcarbonyl)-2-quinolinecarboxaldehyde
Synonym: FQCA
MF: C₁₅H₉NO₃
FW: 251.2
Purity: ≥95%
Ex./Em. Max: 490/590 nm
Supplied as: A 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

3-(2-Furoyl)quinoline-2-carboxaldehyde is supplied as a solid. A stock solution may be made by dissolving the 3-(2-furoyl)quinoline-2-carboxaldehyde in water. 3-(2-Furoyl)quinoline-2-carboxaldehyde is slightly soluble (0.1-1 mg/ml) in water. We do not recommend storing the aqueous solution for more than one day.

Description

3-(2-Furoyl)quinoline-2-carboxaldehyde is a reactive probe for the derivatization of primary amines.¹ 3-(2-Furoyl)quinoline-2-carboxaldehyde reacts with primary amines to form fluorescent isoindoles, which display excitation/emission maxima of 490/590 nm, respectively. 3-(2-Furoyl)quinoline-2-carboxaldehyde has been used for protein labeling.²

References

1. Beale, S.C., Hsieh, Y.-Z., Wiesler, D., *et al.* Application of 3-(2-furoyl)quinoline-2-carbaldehyde as a fluorogenic reagent for the analysis of primary amines by liquid chromatography with laser-induced fluorescence detection. *J. Chromatogr.* **499**, 579-587 (1990).
2. Stoyanov, A.V., Ahmadzadeh, H., and Krylov, S.N. Heterogeneity of protein labeling with a fluorogenic reagent, 3-(2-furoyl)quinoline-2-carboxaldehyde. *J. Chromatogr. B Analyt. Technol. Biomed. Life Sci.* **780(2)**, 283-287 (2002).

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, 02/13/2026

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