

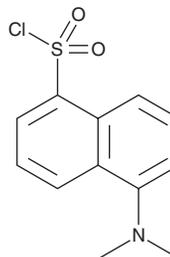
PRODUCT INFORMATION



Dansyl Chloride

Item No. 27458

CAS Registry No.: 605-65-2
Formal Name: 5-(dimethylamino)-1-naphthalenesulfonyl chloride
Synonyms: DNSCl, NSC 83616
MF: C₁₂H₁₂ClNO₂S
FW: 269.7
Purity: ≥95%
UV/Vis.: λ_{max}: 210, 265, 369 nm
Ex./Em. Max: 340/535 nm, respectively, in acetone
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

Dansyl chloride is supplied as a solid. A stock solution may be made by dissolving the dansyl chloride in the solvent of choice, which should be purged with an inert gas. Dansyl chloride is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of dansyl chloride in these solvents is approximately 1.6, 10, and 3 mg/ml, respectively.

Dansyl chloride is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, dansyl chloride should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Dansyl chloride has a solubility of approximately 0.02 mg/ml in a 1:40 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Dansyl chloride is a reactive probe for the derivatization of primary amines, including those on amino acids, peptides, and polyamines, for detection by HPLC.¹⁻³ It has also been used for the derivatization of compounds containing phenol groups, such as steroids, for detection by MS/MS.⁴ Dansyl chloride is a fluorescent probe for proteins and enzymes.⁵ It displays excitation/emission maxima of 340/535 nm, respectively, in acetone.

References

1. Walker, J.M. The dansyl method for identifying N-terminal amino acids. *Methods in molecular biology: Basic protein and peptide protocols*. Walker, J.M., editor, Humana Press, Inc. (1994).
2. Takeuchi, T. HPLC of amino acids as dansyl and dabsyl derivatives. *J. Chromatogr. Lib.* **70**, 229-241 (2005).
3. Hunter, K.J. A dansyl chloride-HPLC method for the determination of polyamines. *Methods in molecular biology: Polyamine protocols*. Morgan, D.M.L., editor, Humana Press, Inc. (1998).
4. Santa, T. Derivatization reagents in liquid chromatography/electrospray ionization tandem mass spectrometry. *Biomed. Chromatogr.* **25**(1-2), 1-10 (2011).
5. Chen, R.F. and Scott, C.H. Atlas of fluorescence spectra and lifetimes of dyes attached to protein. *Anal. Lett.* **18**(4), 393-421 (1985).

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, 10/19/2022

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