

# PRODUCT INFORMATION



## DAUDA

Item No. 10005188

**CAS Registry No.:** 73025-02-2  
**Formal Name:** 11-[[[5-(dimethylamino)-1-naphthalenyl]sulfonyl]amino]-undecanoic acid

**Synonym:** 11-(Dansylamino)undecanoic Acid

**MF:** C<sub>23</sub>H<sub>34</sub>N<sub>2</sub>O<sub>4</sub>S

**FW:** 434.6

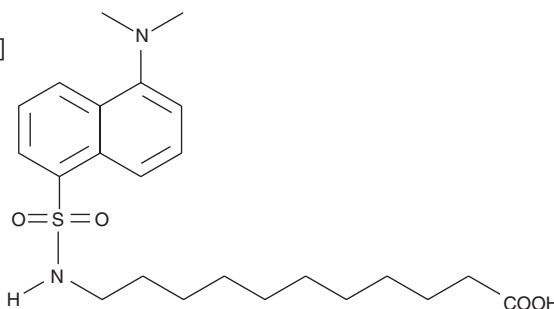
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 217, 251, 336 nm

**Supplied as:** A crystalline 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

DAUDA is supplied as a crystalline solid. A stock solution may be made by dissolving the DAUDA in the solvent of choice, which should be purged with an inert gas. DAUDA is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of DAUDA in ethanol is approximately 30 mg/ml and approximately 25 mg/ml in DMSO and DMF.

DAUDA is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, DAUDA should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. DAUDA has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

DAUDA is an environment-sensitive fluorescent fatty acid analogue that alters its fluorescent emission spectra and intensities on entry into binding proteins. DAUDA has been used to determine the relative affinity of natural fatty acids for polymorphs of the *Schistosoma mansoni* Sm14 fatty acid-binding protein by monitoring excitation and emission at 345 and 543 nm, respectively.<sup>1</sup>

### Reference

1. Ramos, C.R.R., Figueredo, R.C.R., Pertinhez, T.A., *et al.* Gene structure and M20T polymorphism of the *Schistosoma mansoni* Sm14 fatty acid-binding protein. Molecular, functional, and immunoprotection analysis. *J. Biol. Chem.* **278**(15), 12745-12751 (2003).

#### 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, 11/09/2023

#### 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