

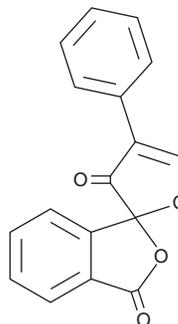
# PRODUCT INFORMATION



## Fluorescamine

Item No. 34286

**CAS Registry No.:** 38183-12-9  
**Formal Name:** 4-phenyl-spiro[furan-2(3H),1'(3'H)-isobenzofuran]-3,3'-dione  
**Synonyms:** Fluram, Ro 20-7234  
**MF:** C<sub>17</sub>H<sub>10</sub>O<sub>4</sub>  
**FW:** 278.3  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 222, 237 nm  
**Ex./Em. Max:** 370/488 nm, in acetone  
**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

Fluorescamine is supplied as a crystalline solid. A stock solution may be made by dissolving the fluorescamine in the solvent of choice, which should be purged with an inert gas. Fluorescamine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of fluorescamine in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Fluorescamine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, fluorescamine should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Fluorescamine has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Fluorescamine is an amine-reactive fluorescent probe.<sup>1</sup> It has commonly been used as a reagent in fluorometric protein assays.<sup>2,3</sup> It displays excitation/emission maxima of 370/488 nm, respectively, in acetone.<sup>1</sup>

### References

1. 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).
2. Udenfriend, S., Stein, S., Böhlen, P., et al. Fluorescamine: A reagent for assay of amino acids, peptides, proteins, and primary amines in the picomole range. *Science* **178(4063)**, 871-872 (1972).
3. Tsukagoshi, K., Tanaka, A., Nakajima, R., et al. On-line capillary zone electrophoretic separation-chemiluminescence detection of protein labeled with fluorescamine. *Anal. Sci.* **12**, 525-528 (1996).

#### 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, 10/17/2022

#### 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