

Fluorescamine

Fluorescamine is nonfluorescent until it reacts with primary amines. The reaction is completed in miniseconds and can be used to detect amine-containing compounds including peptides and proteins.



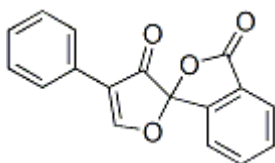
Product attributes

Excitation/Emission	381/470 nm (reaction product)
---------------------	-------------------------------

Product Description

Fluorescamine is nonfluorescent until it reacts with primary amines. The reaction is completed in miniseconds and can be used to detect amine-containing compounds including peptides and proteins. Unreacted fluorescamine is rapidly converted to a non-fluorescent product by reaction with water, making amine detection highly specific.

- $\lambda_{Ex}/\lambda_{Em}(\text{CH}_3\text{CN}) = 381/470 \text{ nm}$ (for reaction product with amines)
- Off-white solid soluble in DMF or DMSO
- Store at -20°C and protect from light
- $\text{C}_{17}\text{H}_{10}\text{O}_4$
- MW: 278.26



References

1. Clin Chim Acta 157, 73 (1986).
2. J Lipid Res 27, 792 (1986).
3. J Chromatog 502, 427 (1990).

This datasheet was generated on January 11, 2026 at 07:35:14 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/fluorescamine/>