## **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

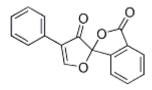
381/470 nm (reaction product)

Call us: 800-304-5357 Email: btinfo@biotium.com

## **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.

- $\lambda_{Ex}/\lambda_{Em}(CH3CN) = 381/470$  nm (for reaction product with amines)
- Off-white solid soluble in DMF or DMSO
- Store at -20°C and protect from light
- C<sub>17</sub>H<sub>10</sub>O<sub>4</sub>
- 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 November 11, 2025 at 11:56:36 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/fluorescamine/">https://biotium.com/product/fluorescamine/</a>