CPM

CPM is a widely used blue fluorescent thiol-reactive dye that is essentially nonfluorescent until it reacts with thiols, making it possible to quantify thiols without a separation step.



Excitation/Emission 384/469 nm (reaction product)

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

CAS number 76877-33-3

Product attributes

Product Description

CPM (full name: 7-diethylamino-3-(4'-maleimidylphenyl)-4-methylcoumarin) is probably the most widely used blue fluorescent thiol-reactive dye. This maleimide derivative of coumarin is essentially nonfluorescent until it reacts with thiols, making it possible to quantify thiols without a separation step. CPM is a good energy acceptor from tryptophan and a good energy donor to fluorescein. Please also see our MTS products.

- Yellow solid soluble in anhydrous DMSO
- MW: 402.45
- [76877-33-3]

References

- 1. Biochemistry 22, 5369 (1983).
- 2. Anal Lett 18, 393 (1985).
- 3. J Biol Chem 265, 14796 (1990).

This datasheet was generated on November 4, 2025 at 10:31:36 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/cpm/