

AMCA-X, SE

AMCA-X SE is an amine-reactive, UV-excitable, blue fluorescent dye.



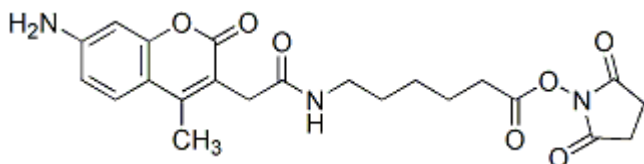
Product attributes

Excitation/Emission	353/442 nm
---------------------	------------

Product Description

AMCA-X SE is an amine-reactive, UV-excitable, blue fluorescent dye.

- $\lambda_{Ex}/\lambda_{Em}(\text{MeOH}) = 353/442 \text{ nm}$
- White solid soluble in DMSO or DMF
- Store at -20°C and protect from light
- $\text{C}_{22}\text{H}_{25}\text{N}_3\text{O}_7$
- MW: 443.45



References

1. Angew Chem Int Ed Engl 40, 402 (2001).
2. J Biol Chem 276, 25236 (2001).
3. J Immunol 164, 5236 (2000).

This datasheet was generated on June 15, 2026 at 01:07:05 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/amca-x-se-6-7-amino-4-methylcoumarin-3-acetyl-amino-hexanoic-acid-succinimidyl-ester/>