Z-FR-AMC

Z-FR-AMC is a substrate for serine proteases, including cathepsins, kallikrein and plasmin.



Product attributes

CAS number 65417-22-0

Excitation/Emission 342/441 nm (end product)

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

Product Description

Z-FR-AMC is a substrate for serine proteases, including cathepsins, kallikrein and plasmin.

- $\lambda_{Ex}/\lambda_{Em}$ of substrate = 330/390 nm (weak fluorescence)
- $\lambda_{Ex}/\lambda_{Em}$ of end product (AMC) = 342/441 nm (strong fluorescence)
- Off-white solid soluble in DMSO
- Store at 4°C and protect from light
- C₃₃H₃₇CIN₆O₆
- MW: 649.14

References

- 1. J Biol Chem 270, 558 (1995).
- 2. Arch Biochem Biophys 314, 171 (1994).
- 3. FEBS Lett 341, 197 (1994).
- 4. J Biochem (Tokyo) 113, 441 (1993).
- 5. Biochem J 201, 367 (1982).

This datasheet was generated on November 1, 2025 at 07:34:11 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/z-fr-amc-7-amino-4-methylcoumarin-n-cbz-l-phenylalanyl-larginine-amide-hydrochloride/