

Z-FR-AMC

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



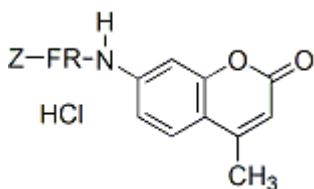
Product attributes

CAS number	65417-22-0
Excitation/Emission	342/441 nm (end product)

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
- $C_{33}H_{37}ClN_6O_6$
- 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 June 9, 2026 at 01:27:43 AM. 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-l-arginine-amide-hydrochloride/>