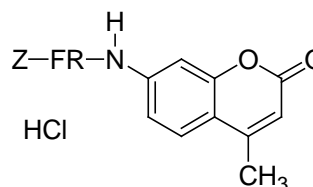


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: 7-Amino-4-methylcoumarin, *N*-CBZ-L-phenylalanyl-L-arginine amide, hydrochloride (Z-FR-AMC, or Z-FR-MCA)

CATALOG #: 10201

MOLECULAR INFORMATION: C₃₃H₃₇N₆O₆
MW: 649.14

**PROPERTIES:**

Color & Form Off-white solid
Purity ≥ 95% by HPLC
Solubility Soluble in DMSO
Absorption/Emission 325 nm/385 nm (weak fluorescence)
(of substrate)
Absorption/Emission 340 nm/440 nm (strong fluorescence)
(of end product AMC)

STORAGE AND HANDLING: Store desiccated at ≤ 4 °C. Protect from light.

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

Ref.: 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).

TOXICITY: Unknown.

FIRST AID: Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.

Disclaimer: Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.