

**Ac-LKKR-AMC**

Catalog number: 13446

Unit size: 1 mg

**Product Details**

---

Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

**Chemical Properties**

---

Appearance	Solid
Molecular Weight	1084.99
Soluble In	DMSO

**Spectral Properties**

---

Excitation Wavelength	341 nm
Emission Wavelength	441 nm

**Applications**

---

Ac-Leu-Lys-Lys-Arg-AMC is a fluorogenic aminopeptidase and endopeptidase substrate. It might be used as a fluorogenic probe for detecting thrombin and metacaspases. Metacaspases are found in plants, fungi, and protozoa. They are cysteine-proteases and are distantly related to metazoan caspases. The protease activity is quantified by release of free fluorescent 7-amino-4-methylcoumarin (AMC) which excites at 360-380 nm and emits at 440-460 nm.