

## Val-AMC

Catalog Number: 13469

Unit Size: 5 mg

### Product Details

---

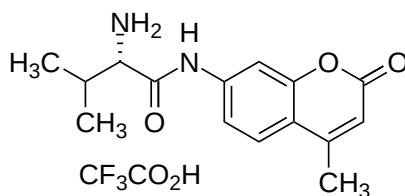
|                    |                                            |
|--------------------|--------------------------------------------|
| Storage Conditions | Freeze (< -15 °C), Minimize light exposure |
| Expiration Date    | 12 months upon receiving                   |

### Chemical Properties

---

|                  |        |
|------------------|--------|
| Appearance       | N/A    |
| Molecular Weight | 388.34 |
| Soluble In       | DMSO   |

Chemical Structure



### Spectral Properties

---

|                       |        |
|-----------------------|--------|
| Excitation Wavelength | 341 nm |
| Emission Wavelength   | 441 nm |

### Applications

---

Val-AMC is a sensitive fluorogenic protease substrate. The non-fluorescent substrate generates a bright blue fluorescent AMC product that has an emission spectra that can be easily detected with the DAPI filter set. It can be used for monitoring protease activity, making it a valuable research tool in enzymology, inhibitor screening, and mechanistic protease studies.