

## Gln-AMC

Catalog Number: 13495

Unit Size: 5 mg

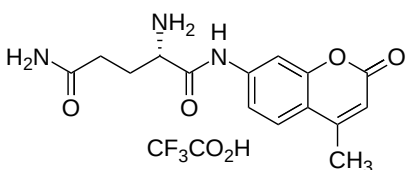
### Product Details

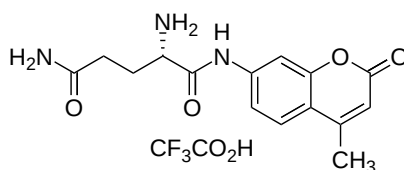
---

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

### Chemical Properties

---

|                    |                                                                                    |
|--------------------|------------------------------------------------------------------------------------|
| Appearance         | N/A                                                                                |
| Molecular Weight   | 417.34                                                                             |
| Soluble In         | DMSO                                                                               |
| Chemical Structure |  |



### Spectral Properties

---

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

### Applications

---

Gln-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.