

# Bodi Fluor™ 488 amine

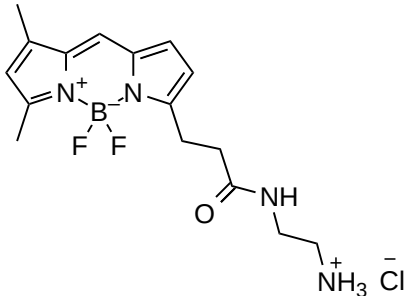
Catalog Number: 702

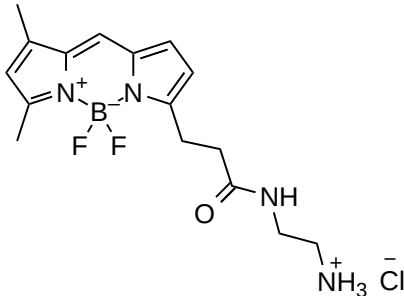
Unit Size: 5 mg

## Product Details

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

## Chemical Properties

|                    |                                                                                    |
|--------------------|------------------------------------------------------------------------------------|
| Appearance         | Solid orange-red                                                                   |
| Molecular Weight   | 370.64                                                                             |
| Soluble In         | DMSO                                                                               |
| Chemical Structure |  |



## Spectral Properties

|                       |     |
|-----------------------|-----|
| Excitation Wavelength | N/A |
| Emission Wavelength   | N/A |

## Applications

Bodi Fluor 488 amine is an excellent building block for developing fluorescent probes. Bodi Fluor 488 is the same fluorophore as BODIPY FL (4,4-difluoro-4-bora-3a,4a-diaza-s-indacene FL). It is a popular lipophilic fluorescent dye molecule used in biology, chemistry, and materials science. BODI Fluor 488 has high fluorescence quantum yields, photostability, and brightness. When excited with light in the blue to green range (around 488 to 515 nanometers), it emits around 510 to 520 nanometers. Researchers and scientists use BODIPY FL for various purposes, such as labeling biomolecules (proteins, nucleic acids, etc.) for fluorescence microscopy, flow cytometry, and other imaging techniques. It can also be employed as a probe for sensing and detecting various analytes or as a fluorescent marker in chemical and biological assays.