

# iFluor® 750-beta-Amyloid (1-28) [beta-Amyloid (1-28), iFluor® 750-Labeled]

Catalog Number: 81033

Unit Size: 100 ug

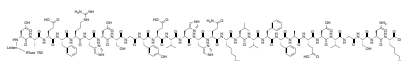
## Product Details

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

## Chemical Properties

|                  |         |
|------------------|---------|
| Appearance       | Solid   |
| Molecular Weight | 4941.46 |
| Soluble In       | DMSO    |

Chemical Structure



## Spectral Properties

|                       |        |
|-----------------------|--------|
| Excitation Wavelength | 757 nm |
| Emission Wavelength   | 779 nm |

## Applications

iFluor® 750-beta-Amyloid (1-28) is an iFluor® 750-labeled peptide used extensively in neuroscience research, particularly in studies related to Alzheimer's disease. iFluor® 750 tag is conjugated to the N-termini of beta-Amyloid (1-42) peptide with sequence: iFluor® 750-Asp-Ala-Glu-Phe-Arg-His-Asp-Ser-Gly-Tyr-Glu-Val-His-His-Gln-Lys-Leu-Val-Phe-Phe-Ala-Glu-Asp-Val-Gly-Ser-Asn-Lys-OH (DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVG). It has a near-IR fluorescence. The spectral properties of iFluor® 750 label allow researchers to visualize and track the fluorescent peptide in biological systems using the most common settings either in a fluorescence microscope or flow cytometer. AAT Bioquest's iFluor® 750-beta-Amyloid (1-28) material is thoroughly analyzed to ensure its quality by HPLC, electrophoresis and cell functional analysis. While the full-length Aβ peptides like Aβ(1-40) and Aβ(1-42) are more commonly associated with plaque formation, Aβ(1-28) is often used for specific mechanistic and biochemical studies. It is used as a monomeric starting material in seeding and fibril formation experiments.