

iFluor® 488-beta-Amyloid (1-28) [beta-Amyloid (1-28), iFluor® 488-Labeled]

Catalog Number: 81010

Unit Size: 100 ug

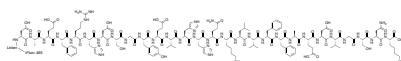
Product Details

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

Chemical Properties

Appearance	Solid
Molecular Weight	4570.89
Soluble In	DMSO

Chemical Structure



Spectral Properties

Excitation Wavelength	491 nm
Emission Wavelength	516 nm

Applications

iFluor® 488-beta-Amyloid (1-28) is an iFluor® 488-labeled peptide used extensively in neuroscience research, particularly in studies related to Alzheimer's disease. iFluor® 488 tag is conjugated to the N-termini of beta-Amyloid (1-42) peptide with sequence: iFluor® 488-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 green fluorescence. The spectral properties of iFluor® 488 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® 488-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.