

iFluor® 647-beta-Amyloid (1-42) [beta-Amyloid (1-42), iFluor® 647-Labeled]

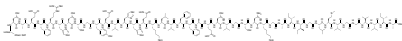
Catalog Number: 81140

Unit Size: 100 ug

Product Details

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

Chemical Properties

Appearance	Solid blue
Molecular Weight	6373.34
Soluble In	DMSO
Chemical Structure	

Spectral Properties

Excitation Wavelength	656 nm
Emission Wavelength	670 nm

Applications

iFluor® 647-beta-Amyloid (1-42) is an iFluor® 647-labeled peptide used extensively in neuroscience research, particularly in studies related to Alzheimer's disease. iFluor® 647 tag is conjugated to the N-termini of beta-Amyloid (1-42) peptide with sequence: iFluor® 647-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-Gly-Ala-Ile-Ile-Gly-Leu-Met-Val-Gly-Gly-Val-Val-Ile-Ala-OH (DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVVIA). It has a far-red fluorescence. The spectral properties of iFluor® 647 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® 647-beta-Amyloid (1-42) material is thoroughly analyzed to ensure its quality by HPLC, electrophoresis and cell functional analysis. Beta-Amyloid (1-42) is a peptide fragment derived from the amyloid precursor protein (APP). It is the primary component of amyloid plaques found in the brains of individuals with Alzheimer's disease. Beta-Amyloid (1-42) is used in neurotoxicity, neuroinflammation and drug screening studies. Aβ(1-42) aggregates into fibrils and oligomers that are neurotoxic.