

Thioflavin T *CAS 2390-54-7*

Catalog number: 23060

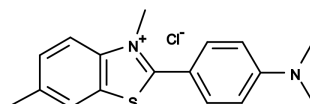
Unit size: 1 g

Product Details

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

Chemical Properties

Appearance	Yellow solid
Molecular Weight	318.87
Soluble In	DMSO
Chemical Structure	


Spectral Properties

Excitation Wavelength	349 nm
Emission Wavelength	454 nm

Applications

Thioflavin T (ThT), also called Basic Yellow 1 or CI 49005, is a benzothiazole salt. The dye is used to visualize plaques composed of beta-amyloid found in the brains of Alzheimer's disease patients. When it binds to beta sheets, such as those in amyloid oligomers, the dye undergoes a characteristic 120 nm red shift of its excitation spectrum that may be selectively excited at 450 nm, resulting in a fluorescence signal at 482 nm. ThT binds rapidly and specifically to the anti-parallel beta-sheet fibrils formed from synthetic beta-amyloid (1-40), but does not bind to monomer or oligomeric intermediates. The binding of ThT does not interfere with the aggregation of beta-amyloid peptide into amyloid fibrils.