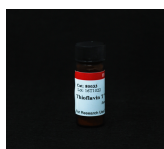


Thioflavin T, High Purity Grade

A cell-permeable dye that exhibits enhanced fluorescence upon binding to amyloid fibrils, and has been used in histology and for protein characterization.



Product attributes

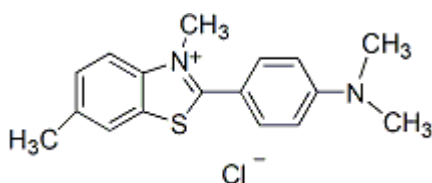
CAS number	2390-54-7
Apoptosis/viability marker	Amyloid stain
Colors	Blue
Excitation/Emission	450/482 nm

Product Description

A cell-permeable benzothiazole dye that exhibits enhanced fluorescence (Ex/Em 450/482 nm) upon binding to amyloid fibrils. Useful in monitoring stacked β -sheet aggregates, Thioflavin T has also been used in histology and for protein characterization.

Biotium offers high purity grade of Thioflavin T with HPLC purity greater than 97%.

- Yellow solid soluble in DMF and DMSO
- Stored desiccated at 4 °C
- C₁₇H₁₉ClN₂S
- MWt: 318.86
- [2390-54-7]



References

LeVine, H. *Protein Sci.* 1993 Mar; 2(3): 404–410. doi: [10.1002/pro.5560020312](https://doi.org/10.1002/pro.5560020312)

This datasheet was generated on June 6, 2026 at 07:26:46 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/thioflavin-t-high-purity-grade/>