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

Call us: 800-304-5357 Email: btinfo@biotium.com

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
- C₁₇H₁₉CIN₂S
- MWt: 318.86
- [2390-54-7]

References

LeVine, H. Protein Sci. 1993 Mar; 2(3): 404-410. doi: 10.1002/pro.5560020312

This datasheet was generated on November 6, 2025 at 09:18:14 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/thioflavin-t-high-purity-grade/