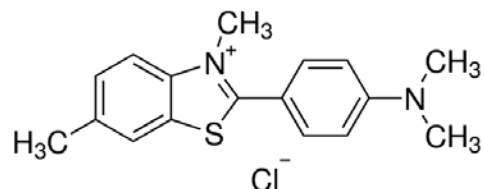


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: Thioflavin T "high purity grade"**CATALOG NUMBER:** 80033**MOLECULAR INFORMATION:** C₁₇H₁₉ClN₂S
Mwt: 318.86
[2390-54-7]**PROPERTIES:****Color & Form** Yellow solid
Purity ≥ 97% by HPLC
Solubility Soluble in DMF and DMSO**STORAGE AND HANDLING:** Stored desiccated at 4 °C upon receipt.**APPLICATION:** A cell-permeable benzothiazole dye that exhibits enhanced fluorescence upon binding to amyloid fibrils. Useful in monitoring stacked β sheet aggregates. Thioflavin T has also been used in histology and for protein characterization.**TOXICITY:** Unknown**FIRST AID:** Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.

Disclaimer: Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.