

Chitosan-Bodipy TMR

Catalog number: H-C00M

MW: 278 kDa

Solubility: 1 mg/mL in H₂O

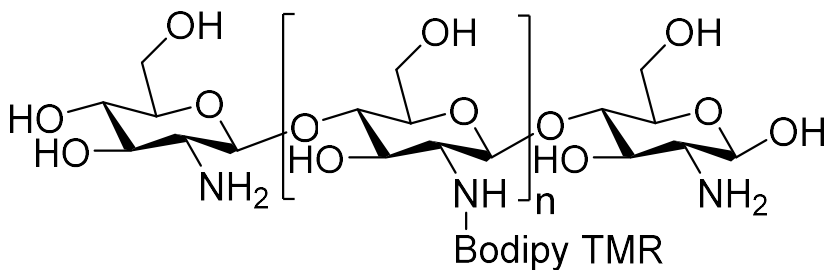
Storage and Handling: store in -20 °C, light sensitive

Background: Chitosan, the deacetylated derivative of chitin, has attracted considerable attention in recent years, because of its remarkable properties, which make it particularly suited in such applications as, e.g., biomedical aids, tissue engineering, drug delivery, packaging materials and microcapsules for diverse end-uses. Fluorescent labeled chitosan is used for tracing chitosan used in a biological system.

References:

Fernandes, S.C.M., et al, "a study of the distribution of chitosan onto and within a paper sheet using a fluorescent chitosan derivative" *Carbohydrate Polymers*, 2009, 78, 760-766

Kumar, M.N.V.R., et al, "Chitosan chemistry and pharmaceutical perspectives" *Chem. Rev.*, 2004, 104, 6017-6084



Hazardous Properties and Cautions: The toxicological and pharmacological properties of this compound are not fully known. For further information see the MSDS on request. This product is manufactured and shipped only in small quantities, intended for research and development in a laboratory utilizing prudent procedures for handling chemicals of unknown toxicity, under the supervision of persons technically qualified to evaluate potential risks and authorized to enforce appropriate health and safety measures. As with all research chemicals, precautions should be taken to avoid unnecessary exposures or risks.

Warranty and Disclaimer: Echelon warrants the product conforms to the specifications stated herein. In the event of nonconformity, Echelon will replace products or refund purchase price, at its sole option, and Echelon shall not be responsible for any other loss or damage, whether known or foreseeable to Echelon. No other warranties apply, express or implied, including but not limited to warranty of fitness for any purpose or implied warranty of merchantability. Purchaser is solely responsible for all consequences of its use of the product and Echelon assumes no responsibility therefore, including success of purchaser's research and development, or health or safety of any uses of the product.

Technical Data Sheet, Rev 1, 9/25/2024 – **For research use only.** Not intended for diagnostic or therapeutic use.