3,6,9,12-Tetraoxatetradecane-1,14-diyl dimethanethiosulfonate

Bictum

Product attributes

Call us: 800-304-5357

3,6,9,12-Tetraoxatetradecane-1,14-diyl dimethanethiosulfonate is a water soluble MTS cross-linker.

Product Description

3,6,9,12-Tetraoxatetradecane-1,14-diyl dimethanethiosulfonate is a water soluble MTS cross-linker. Stock solutions made in anhydrous DMF are stable for at least two months if stored at -20°C. However, stock solutions made in water or buffer should be used immediately.

- Off-white solid soluble in DMF or water
- Store desiccated at -20°C
- C₁₂H₂₆O₈S₄
- MW: 426.57

This datasheet was generated on November 25, 2025 at 07:41:24 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/36912-tetraoxatetradecane-114-diyl-dimethanethiosulfonate/