

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

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

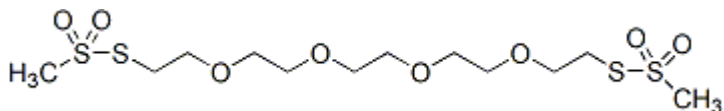


Product attributes

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
- $C_{12}H_{26}O_8S_4$
- MW: 426.57



This datasheet was generated on January 28, 2026 at 08:15:05 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/36912-tetraoxatetradecane-114-diyl-dimethanethiosulfonate/>