

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

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

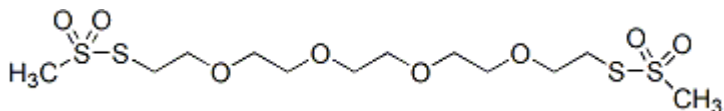


Product attributes

Product Description

3,6,9,12-Tetraoxatetradecane-1,14-diyl dimethanesulfonate 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
- $\text{C}_{12}\text{H}_{26}\text{O}_8\text{S}_4$
- MW: 426.57



This datasheet was generated on June 16, 2026 at 11:34:47 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/36912-tetraoxatetradecane-114-diyl-dimethanesulfonate/>