

3,6,9-Trioxaundecane-1,11-diyl dimethanethiosulfonate

3,6,9-Trioxaundecane-1,11-diyl dimethanethiosulfonate is a water soluble MTS cross-linker.

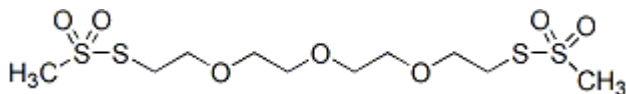


Product attributes

Product Description

3,6,9-Trioxaundecane-1,11-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
- $C_{10}H_{22}O_7S_4$
- MW: 382.52



This datasheet was generated on January 28, 2026 at 09:42:39 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/369-trioxaundecane-111-diyl-dimethanethiosulfonate/>