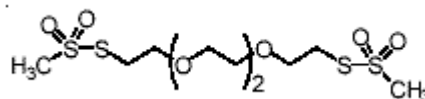


## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** 3,6,9-Trioxaundecane-1,11-diyl dimethanethiosulfonate**CATALOG #** 91043**MOLECULAR INFORMATION:** C<sub>10</sub>H<sub>22</sub>O<sub>7</sub>S<sub>4</sub>  
MWt: 382.52**PROPERTIES:**

<b>Color &amp; Form</b>	Of-white solid
<b>Purity</b>	≥ 95% by TLC
<b>Solubility</b>	DMF or water
<b>Absorption/Emission</b>	N/A
<b>Extinction Coefficient</b>	N/A

**STORAGE AND HANDLING:** Store the solid at -20 °C. Stock solution made in anhydrous DMF is good for at least two months if stored at -20 °C. Stock solution made in water or buffer should be used immediately.**APPLICATION:** 3,6,9-Trioxaundecane-1,11-diyl dimethanethiosulfonate is a water soluble MTS cross-linker.**TOXICITY:** Not established. Not listed by NTP, IARC or OSHA.

<b>FIRST AID:</b>	Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.
-------------------	---

<b>Disclaimer:</b> <i>Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.</i>
--