

3,6-Dioxaoctane-1,8-diyl dimethanesulfonate

3,6-dioxaoctane-1,8-diyl dimethanesulfonate is a water soluble MTS cross-linker.



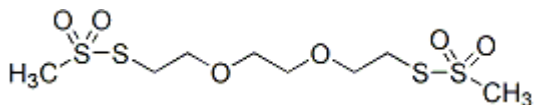
Product attributes

Excitation/Emission	375/508 nm
---------------------	------------

Product Description

3,6-dioxaoctane-1,8-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 a buffer should be used immediately.

- Off-white liquid soluble in DMF or water
- Store desiccated at -20 °C
- $C_8H_{18}O_6S_4$
- MW: 338.49



This datasheet was generated on June 3, 2026 at 11:33:30 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/36-dioxaoctane-18-diyl-dimethanesulfonate/>