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

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

side.

Product attributes

Excitation/Emission

375/508 nm

Call us: 800-304-5357 Email: btinfo@biotium.com

Product Description

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

- Off-white liquid soluble in DMF or water
- Store desiccated at -20°C
- C₈H₁₈O₆S₄
- MW: 338.49

This datasheet was generated on November 3, 2025 at 04:47:26 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/36-dioxaoctane-18-diyl-dimethanethiosulfonate/