# **MTSES**

MTSES (2-Sulfonatoethyl methanethiosulfonate sodium salt) is a negatively charged MTS reagent.



#### **Product attributes**

Call us: 800-304-5357

## **Product Description**

MTSES (2-Sulfonatoethyl methanethiosulfonate sodium salt) is a negatively charged MTS reagent.

- White solid soluble in DMSO
- Store desiccated at -20°C
- C<sub>3</sub>H<sub>7</sub>NaO<sub>5</sub>S<sub>3</sub>
- MW: 242.21

### References

- 1. Biochemistry 33, 6840 (1994).
- 2. Science 258, 307 (1992).

This datasheet was generated on November 25, 2025 at 07:44:27 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/mtses-2-sulfonatoethyl-methanethiosulfonate-sodium-salt/