

MTSES

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

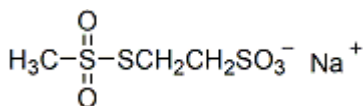


Product attributes

Product Description

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

- White solid soluble in DMSO
- Store desiccated at -20 °C
- C₃H₇NaO₅S₃
- MW: 242.21



References

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

This datasheet was generated on June 14, 2026 at 11:53:55 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/mtses-2-sulfonatoethyl-methanethiosulfonate-sodium-salt/>