



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

## MTSEA

MTSEA (2-Aminoethyl methanethiosulfonate hydrobromide) is a positively charged MTS reagent.

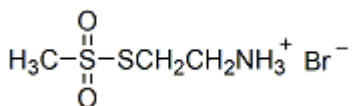


### Product attributes

## Product Description

MTSEA (2-Aminoethyl methanethiosulfonate hydrobromide) is a positively charged MTS reagent.

- Off-white solid soluble in water, DMSO or DMF
- Store desiccated at -20 °C
- $C_3H_{10}BrNO_2S_2$
- MW: 236.15



## References

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

This datasheet was generated on January 9, 2026 at 07:16:37 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/mtsea-2-aminoethyl-methanethiosulfonate-hydrobromide/>