



Glowing products for science

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

MTSPT

MTSPT (3-trimethylammoniumpropyl methanethiosulfonate, bromide salt) is a positively charged MTS reagent.

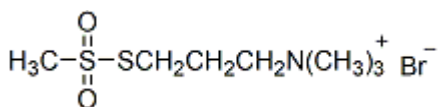


Product attributes

Product Description

MTSPT (3-trimethylammoniumpropyl methanethiosulfonate, bromide salt) is a positively charged MTS reagent.

- White solid soluble in water, DMSO or DMF
- Store desiccated at -20°C
- C₇H₁₈BrNO₂S₂
- MW: 292



This datasheet was generated on January 28, 2026 at 08:15:10 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/mtspt-3-trimethylammoniumpropyl-methanethiosulfonate-bromide-salt/>