Call us: 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 November 25, 2025 at 07:39:32 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/mtspt-3-trimethylammoniumpropyl-methanethiosulfonate-bromide-salt/