

## MTSPT

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

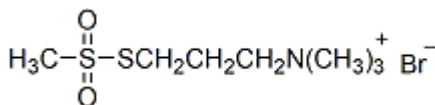


### Product attributes

## 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
- C<sub>7</sub>H<sub>18</sub>BrNO<sub>2</sub>S<sub>2</sub>
- MW: 292



This datasheet was generated on June 22, 2026 at 01:18:21 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/mtspt-3-trimethylammoniumpropyl-methanethiosulfonate-bromide-salt/>