



Glowing products for science

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

## DMSO, Anhydrous

Anhydrous DMSO (dimethyl sulfoxide; methyl sulfoxide) is recommended for preparing stock solutions of AM ester dyes, reactive dyes and other related moisture-sensitive products.



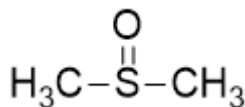
### Product attributes

CAS number	67-68-5
Storage Conditions	Store at room temperature, Desiccate

## Product Description

Anhydrous DMSO (dimethyl sulfoxide; methyl sulfoxide) is recommended for preparing stock solutions of AM ester dyes, reactive dyes and other related moisture-sensitive products.

- Colorless liquid at room temperature
- Store desiccated at room temperature
- C<sub>2</sub>H<sub>6</sub>OS
- MW: 78.13
- [67-68-5]



This datasheet was generated on January 21, 2026 at 12:26:45 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/dmsO-anhydrous-dimethyl-sulfoxide-or-methyl-sulfoxide/>