

46117 Landing Parkway, Fremont, CA 94538 U.S.A. Tel: 1-510-265-1027; Fax: 1-510-265-1352

www.biotium.com

PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: Dimethyl sulfoxide, anhydrous (DMSO)

CATALOG # 90082

MOLECULAR $(CH_3)_2SO$ **INFORMATION:** MWt: 78.13

[67-68-5]

H₃C^SCH₃

PROPERTIES:

Color & Form Colorless liquid

Solubility N/A **Purity** $\geq 99.9\%$

Titration 0.005% H₂O (maximum) **Residual on** 0.0005% (maximum)

Evaporation

STORAGE AND Store desiccated at room temperature

HANDLING:

APPLICATION: Dimethyl sulfoxide, anhydrous is an organic solvent used for dissolving moisture-

sensitive compounds.

TOXICITY: Not established. Not listed by NTP, IARC or OSHA.

FIRST AID: Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with

plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing,

move individual to fresh air and seek medical advice immediately.

Disclaimer: Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.