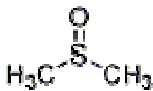


Instruction Manual

Catalog Number	PK-CA707-90082
Description	Anhydrous DMSO is recommended for preparing stock solutions of AM ester dyes, reactive dyes and other related moisture-sensitive products.
Quantity	10 ml
Excitation / Emission Maxima	NA
Molecular Structure	
Molecular Weight / Molecular Formula	78.13 Da; C ₂ H ₆ OS
Purity	>99.9%; Titration: 0.005% H ₂ O (maximum); Residual on Evaporation: 0.0005% (maximum)
Appearance / Formulation / Solubility	Colorless liquid.
Storage & Stability	Store desiccated at room temperature.
Applications	See Description
References	NA
Caution	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.

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.