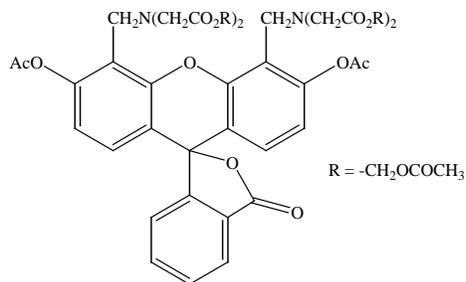


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: Calcein AM, 4mM in anhydrous DMSO**CATALOG #** 80011-1**MOLECULAR INFORMATION:** C₄₆H₄₆N₂O₂₃
MWt: 995**PROPERTIES:**

Color & Form	Off-white liquid
Purity	≥90% by HPLC
Solubility	N/A
Absorption/Emission	$\lambda_{\text{abs}} < 300\text{nm}$ (before hydrolysis); $\lambda_{\text{abs}} \backslash \lambda_{\text{em}} = 494/517\text{ nm}$ (pH 9, after hydrolysis).
Extinction Coefficient	Not determined.

STORAGE AND HANDLING:

Stored desiccated at -20 °C upon receiving and protect from light.

APPLICATION:

Calcein AM is a widely used green fluorescent cell marker. Calcein AM is membrane-permeant and thus can be introduced into cells via incubation. Once inside the cells, calcein AM is hydrolyzed by endogenous esterase into the highly negatively charged green fluorescent calcein, which is retained in the cytoplasm. Calcein AM is an excellent tool for the studies of cell membrane integrity and for long term cell tracing.

Ref.: 1) *J. Immunol. Meth.* **177**, 101(1994); 2) *J. Immunol. Meth.* **172**, 227(1994); 3) *J. Immunol. Meth.* **172**, 115(1994); 4) *Human Immunol.* **37**, 264(1993).

TOXICITY: Unknown

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.