

6-Carboxyfluorescein diacetate (single isomer)

Instruction Manual

| | |
|---------------------------------------|---|
| Catalog Number | PK-CA707-51021 |
| Description | 6-CFDA is membrane-permeant and thus can be loaded into cells via incubation. Once inside the cells, 6-CFDA is hydrolyzed by intracellular esterases to 6-carboxyfluorescein. |
| Quantity | 100 mg |
| Excitation / Emission Maxima | $\lambda_{ex}/\lambda_{em} = 492/517$ nm (pH 9.0, after hydrolysis); Extinction coefficient = 78,000 (pH 9.0, after hydrolysis) |
| Molecular Structure | |
| Molecular Weight / Molecular Formula | 460 Da; C ₂₅ H ₁₆ O ₉ |
| Purity | >99% (as determined by TLC) |
| Appearance / Formulation / Solubility | Off-white solid; soluble in DMSO. |
| Storage & Stability | Store at $\leq 4^{\circ}\text{C}$ and protect from light |
| Applications | Fluorescent pH indicator |
| References | <ol style="list-style-type: none"> 1) Boitano, S., et al. J. Cell Sci. 98, 343(1991) 2) Goodall, H., et al. Nature 295, 524(1982) 3) Hansson, Y., et al. J. Immunol. Meth. 100, 261(1987) 4) Bruning, J.W., et al. J. Immunol. Meth. 33, 33(1980) |
| 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.