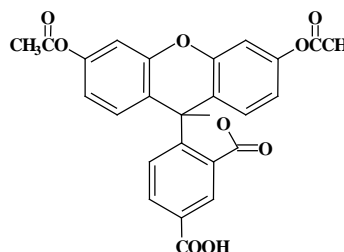


## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** 5-CFDA (5-Carboxyfluorescein diacetate, single isomer)**CATALOG #:** 51018**MOLECULAR INFORMATION:** C<sub>25</sub>H<sub>16</sub>O<sub>9</sub>  
Mwt: 460  
[79955-27-4]**PROPERTIES:**

<b>Color &amp; Form</b>	Off-white solid
<b>Purity</b>	≥ 99% by HPLC
<b>Solubility</b>	Soluble in DMSO
$\lambda_{ex} \backslash \lambda_{em}$	492nm \ 517nm (pH 9, after hydrolysis)
<b>Extinction Coefficient (<math>\epsilon</math>)</b>	78,000 (pH 9, after hydrolysis)

**STORAGE AND HANDLING:** Stored desiccated at 4°C upon receipt. Protect from light, especially when in solution.**APPLICATION:** 5-CFDA is membrane-permeant and thus can be loaded into cells via incubation. Once inside the cells, CFDA is hydrolyzed by intracellular esterases to carboxyfluorescein (**51013**).

Ref.: Boitano, S., et al. *J. Cell Sci.* **98**, 343(1991); Goodall, H., et al. *Nature* **295**, 524(1982); Hansson, Y., et al. *J. Immunol. Meth.* **100**, 261(1987); Bruning, J.W., et al. *J. Immunol. Meth.* **33**, 33(1980).

**TOXICITY:** Unknown

<b>FIRST AID:</b>	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.
-------------------	---

<b>Disclaimer:</b> 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.
---