

## Product Information

### CFDA, SE

Cat. No.	Product Name	Unit Size
C032	CFDA, SE	25 mg

**Storage upon receipt:**

- -20°C
- Protect from light

### Product Description

**5(6)-CFDA, SE** is a amine-reactive probe which passively diffuse into cells and it is nonfluorescent until the acetate groups are cleaved by intracellular esterases to yield the highly fluorescent, amine-reactive fluorophore. Upon reaction with amine-containing residues of intracellular proteins, these probes form dye protein adducts that are well retained in cells as they move and divide during embryonic development.

### Specifications:

**Chemical name:** 5-(and-6)-Carboxyfluorescein Diacetate

**Molecular Formula:** C<sub>29</sub>H<sub>19</sub>NO<sub>11</sub>

**Molecular Weight:** 557.47

**CAS Number:** 150347-59-4