

---

## Product Information

### CDCFDA, SE

Cat. No.	Product Name	Unit Size
C036	CDCFDA, SE	25 mg

**Storage upon receipt:**

- -20°C
- Protect from light

### Product Description

**5(6)-CDCFDA, SE** is a fixable-cell-permeant, fluorescein-based tracer for very long-term cell labeling.

### Specifications:

**Chemical name:** 5-(and-6)-Carboxy-2',7'-dichlorofluorescein Diacetate, succinimidyl ester

**Molecular Formula:** C<sub>25</sub>H<sub>17</sub>Cl<sub>2</sub>NO<sub>11</sub>

**Molecular Weight:** 626.36

**CAS Number:** 147265-60-9