

CERTIFICATE OF ANALYSIS

Catalog Number C035
Product Name CDCFDA (5-(and-6)-carboxy-2',7'-dichlorofluorescein diacetate)
CAS Number 127770-45-0
Molecular Formula C₂₅H₁₄Cl₂O₉
Molecular Weight 529.28
Appearance white solid
Lot Number

	LOT DATA	SPECIFICATION
Absorption ¹ Maximum Extinction	292 nm 6,400 cm ⁻¹ M ⁻¹	292 ± 3 nm 6,000 ± 500 cm ⁻¹ M ⁻¹
HPLC Purity	96 % at 254 nm	≥ 90 % at 254 nm
NMR Result	meets specification	consistent with structure

1. Solvent: Methanol.