

CERTIFICATE OF ANALYSIS

Catalog Number C036
Product Name CDCFDA, SE (5-(and-6)-carboxy-2',7'-dichlorofluorescein diacetate, succinimidyl ester)
CAS Number 147265-60-9
Molecular Formula C₂₉H₁₇Cl₂NO₁₁
Molecular Weight 626.36
Appearance white solid
Lot Number

	LOT DATA	SPECIFICATION
Absorption ¹ Maximum Extinction	301 nm 6,600 cm ⁻¹ M ⁻¹	300 ± 4 nm 6,500 ± 600 cm ⁻¹ M ⁻¹
HPLC Purity	96 % at 254 nm	≥ 90 % at 254 nm
NMR Result	meets specification	consistent with structure

1. Solvent: Methanol.