

CERTIFICATE OF ANALYSIS

Catalog Number C032
Product Name 5(6)-CFDA, SE; (5-(and-6)-carboxyfluorescein diacetate, succinimidyl ester)
CAS Number 150347-59-4
Molecular Formula C₂₉H₁₉NO₁₁
Molecular Weight 557.47
Appearance white solid
Lot Number

	LOT DATA	SPECIFICATION
Absorption ¹ Maximum Extinction	291 nm 6,400 cm ⁻¹ M ⁻¹	290 ± 4 nm 5,900 ± 900 cm ⁻¹ M ⁻¹
HPLC Purity	96 % at 254 nm	≥ 90 % at 254 nm
NMR Result	meets specification	consistent with structure

1. Solvent: Methanol.