

CERTIFICATE OF ANALYSIS

Catalog Number C031
Product Name 5(6)-CFDA (5-(and-6)-carboxyfluorescein diacetate)
CAS Number 124387-19-5
Molecular Formula C₂₅H₁₆O₉
Molecular Weight 460.4
Appearance off-white solid
Lot Number

	LOT DATA	SPECIFICATION
Absorption ¹ Maximum Extinction	291 nm 5,900 cm ⁻¹ M ⁻¹	290 ± 4 nm 5,500 ± 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.