

CERTIFICATE OF ANALYSIS

Catalog Number	C039
Product Name	CellView Green CMFDA, [5-Chloromethylfluorescein Diacetate]
CAS Number	136832-63-8
Molecular Formula	C ₂₅ H ₁₇ O ₇ Cl
Molecular Weight	464.86
Appearance	white solid
Lot Number	

	LOT DATA	SPECIFICATION
Absorption ¹ Maximum Extinction	290 nm 5,800 cm ⁻¹ M ⁻¹	290 ± 3 nm 5,500 ± 500 cm ⁻¹ M ⁻¹
HPLC Purity	96 % at 254 nm	≥ 95 % at 254 nm
NMR Result	meets specification	consistent with structure

1. Solvent: Acetonitrile.