

CERTIFICATE OF ANALYSIS

Catalog Number C027
Product Name 2',7'-bis-(2-carboxyethyl)-5-(and-6)-carboxyfluorescein (BCECF acid)
CAS Number 85138-49-4
Molecular Formula C₂₇H₂₀O₁₁
Molecular Weight 520.45
Appearance orange solid
Lot Number

	LOT DATA	SPECIFICATION
Absorption ¹		
Maximum	503 nm	503 ± 4 nm
Extinction	87,000 cm ⁻¹ M ⁻¹	83,000 ± 7,000 cm ⁻¹ M ⁻¹
Fluorescence ¹		
Emission Maximum	525 nm	528 ± 4 nm
HPLC		
Purity	96 % at 254 nm	≥ 90 % at 254 nm
NMR		
Result	meets specification	consistent with structure

1. Solvent: Buffer, 50 mM potassium phosphate (pH 9).