

## CERTIFICATE OF ANALYSIS

**Catalog Number** C029  
**Product Name** 2',7'-bis-(2-carboxyethyl)-5-(and-6)-carboxyfluorescein, acetoxymethyl ester (BCECF, AM), 1 mg/mL in DMSO  
**CAS Number** 117464-70-7  
**Appearance** yellow solution  
**Lot Number**

	LOT DATA	SPECIFICATION
<b>Absorption</b> <sup>1</sup> Maximum Extinction	227 nm 73,000 cm <sup>-1</sup> M <sup>-1</sup>	226 ± 4 nm 68,000 ± 6,000 cm <sup>-1</sup> M <sup>-1</sup>
<b>NMR</b> Result	meets specification	consistent with structure

1. Solvent: Acetonitrile.