

## CERTIFICATE OF ANALYSIS

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

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

1. Solvent: Acetonitrile.