

## CERTIFICATE OF ANALYSIS

**Catalog Number** C037  
**Product Name** CellView Blue CMAC, (7-Amino-4-chloromethylcoumarin)  
**Molecular Formula** C<sub>10</sub>H<sub>8</sub>NO<sub>2</sub>Cl  
**Molecular Weight** 209.63  
**Appearance** yellow solid  
**Lot Number**

	LOT DATA	SPECIFICATION
<b>Absorption</b> <sup>1</sup> Maximum Extinction	354 nm 14,000 cm <sup>-1</sup> M <sup>-1</sup>	354 ± 3 nm 14,000 ± 1,000 cm <sup>-1</sup> M <sup>-1</sup>
<b>Fluorescence</b> <sup>1</sup> Emission Maximum	465 nm	464 ± 4 nm
<b>HPLC</b> Purity	96 % at 254 nm	≥ 90 % at 254 nm
<b>NMR</b> Result	meets specification	consistent with structure

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