

## CERTIFICATE OF ANALYSIS

**Catalog Number** C023  
**Product Name** calcein blue, AM, 1 mg/mL in DMSO  
**CAS Number** 168482-84-6  
**Molecular Formula** C<sub>21</sub>H<sub>23</sub>NO<sub>11</sub>  
**Molecular Weight** 465.41  
**Appearance** colorless solution  
**Lot Number**

	LOT DATA	SPECIFICATION
<b>Absorption</b> <sup>1</sup>		
Maximum	323 nm	322 ± 3 nm
Extinction	14,000 cm <sup>-1</sup> M <sup>-1</sup>	13,000 ± 1,000 cm <sup>-1</sup> M <sup>-1</sup>
<b>Fluorescence</b> <sup>1</sup>		
Emission Maximum	439 nm	437 ± 4 nm
<b>HPLC</b>		
Purity	95 % at 254 nm	≥ 80 % at 254 nm
<b>NMR</b>		
Result	meets specification	consistent with structure

1. Solvent: Methanol.