

## CERTIFICATE OF ANALYSIS

**Catalog Number** C044  
**Product Name** Rhodamine 123  
**CAS Number** 62669-70-9  
**Molecular Formula** C<sub>21</sub>H<sub>17</sub>ClN<sub>2</sub>O<sub>3</sub>  
**Molecular Weight** 380.83  
**Appearance** red solid  
**Lot Number**

|   | LOT DATA  | SPECIFICATION  |
|---|---|--|
| <b>Absorption</b> <sup>1</sup><br>Maximum<br>Extinction | 507 nm<br>97,000 cm <sup>-1</sup> M <sup>-1</sup> | 507 ± 3 nm<br>101,000 ± 7,000 cm <sup>-1</sup> M <sup>-1</sup> |
| <b>Fluorescence</b> <sup>1</sup><br>Emission Maximum    | 530 nm  | 530 ± 4 nm   |
| <b>HPLC</b><br>Purity                                   | 95 % at 254 nm                                    | ≥ 90 % at 254 nm   |
| <b>NMR</b><br>Result                                    | meets specification                               | consistent with structure                                      |

1. Solvent: Methanol.