

## 5(6)-CR110, SE [5-(and-6)-Carboxy-rhodamine 110, succinimidyl ester]

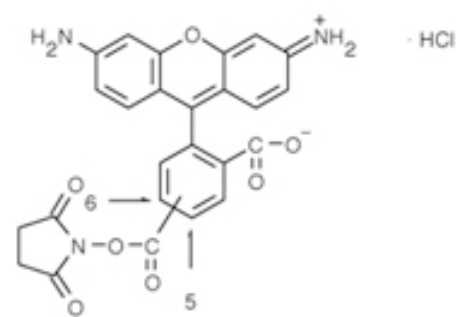
| Catalog Number | Packaging Size |
|----------------|----------------|
| C128           | 5 mg           |

**Storage upon receipt:** -20°C, protect from light

### Introduction

**5(6)-CR110, SE** is a amine-reactive fluorescent probe for preparation of bioconjugates and it is more photo-stable and less pH sensitive (between 4-9) than fluorescein. The labeled conjugated probes can be used for Fluorescence Correlation Spectroscopy (FCS).

### Specifications

|                            |                                                                 |                                                                                     |
|----------------------------|-----------------------------------------------------------------|-------------------------------------------------------------------------------------|
| <b>Label:</b>              | Rhodamine 110                                                   |  |
| <b>Ex/Em:</b>              | 505/525 nm                                                      |                                                                                     |
| <b>Detection Method:</b>   | Fluorescent                                                     |                                                                                     |
| <b>Solubility:</b>         | DMSO, DMF                                                       |                                                                                     |
| <b>Molecular Formula:</b>  | C <sub>25</sub> H <sub>18</sub> ClN <sub>3</sub> O <sub>7</sub> |                                                                                     |
| <b>Molecular Weight:</b>   | 507.89                                                          |                                                                                     |
| <b>CAS Number:</b>         | 254732-34-8                                                     |                                                                                     |
| <b>Storage Conditions:</b> | -20°C, protect from light                                       |                                                                                     |
| <b>Shipping Condition:</b> | Room Temperature                                                |                                                                                     |

### Applications

Fluorescent labeling

### References:

1. Single-molecule detection technologies in miniaturized high-throughput screening: fluorescence intensity distribution analysis.  
Haupts U, Rudiger M, Ashman S, Turconi S, Bingham R, Wharton C, Hutchinson J, Carey C, Moore KJ, Pope AJ  
J Biomol Screen (2003) 0:19-19
2. A single-shot approach for measuring two-photon action cross-section of fluorescent markers  
Shi K, Heikal AA, Liu Z  
Opt Express (2006) 14:8722-8749
3. [Frequency and voltage dependence of the dielectrophoretic trapping of short lengths of DNA and dCTP in a nanopipette.](#)  
Ying L, White SS, Bruckbauer A, Meadows L, Korchev YE, Klenerman D  
Biophys J (2004) 86:1018-1027