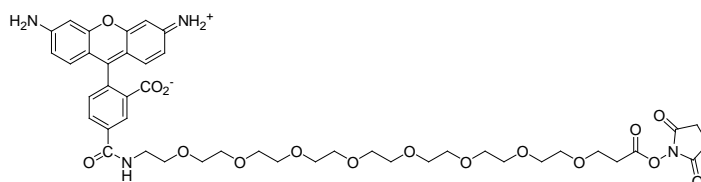


## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** **5-Carboxyrhodamine 110-PEO<sub>8</sub>-propionate, succinimidyl ester, single isomer** (5-CR110-PEO<sub>8</sub>, SE)

**CATALOG #** 90084

**MOLECULAR INFORMATION:** C<sub>44</sub>H<sub>54</sub>N<sub>4</sub>O<sub>16</sub>  
Mwt: 894.92

**PROPERTIES:**

**Color & Form** Orange red solid  
**Purity** ≥ 90% by TLC  
**Solubility** Soluble in DMF or DMSO  
**Absorption/Emission** 502nm/524nm (MeOH)  
**Extinction Coefficient** 39,600

**STORAGE AND HANDLING:** Store desiccated at -20 °C. Protect from light, especially when in solution.

**APPLICATION:** 5-CR110-PEO<sub>8</sub> SE is similar to 5-CR110-PEO<sub>4</sub> SE except that it has an even longer PEG spacer.

**RELATED PRODUCTS** *5-Carboxyrhodamine 110-PEO<sub>4</sub>-propionate, succinimidyl ester, single isomer (#90081)*

**TOXICITY:** Not established. Not listed by NTP, IARC or OSHA.

<b>FIRST AID:</b>	Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.
-------------------	---

<b>Disclaimer:</b> <i>Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.</i>
--