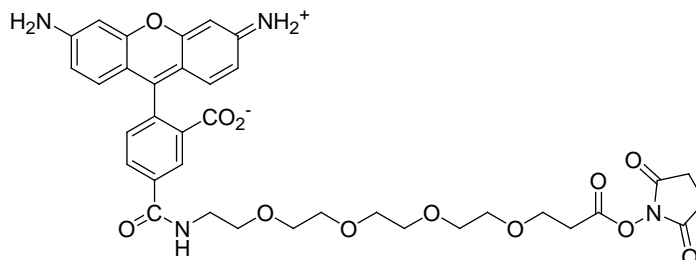


## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** 5-CR110-PEO<sub>4</sub>-propionate SE**CATALOG #** 90081**MOLECULAR INFORMATION:** C<sub>36</sub>H<sub>38</sub>N<sub>4</sub>O<sub>12</sub>  
MWt: 718.71**PROPERTIES:**

**Color & Form** Orange red solid  
**Solubility** Soluble in DMSO or DMF  
**Absorption/Emission** 502/524 nm  
**Extinction Coefficient**

**STORAGE AND HANDLING:** Store desiccated at -20°C and protect from light

**APPLICATION:** 5-CR110-PEO<sub>4</sub>-propionate SE is a derivative of the green fluorescent 5-carboxyrhodamine 110 with a polyethylene oxide spacer between the dye and the amine-reactive group. The spacer group increases the water solubility of the dye as well as the distance between the dye and the molecule to be labeled.

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

**FIRST AID:** 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.

**Disclaimer:** 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.