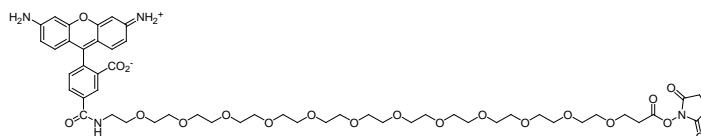


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: 5-Carboxyrhodamine 110-PEO₁₂-propionate, succinimidyl ester, single isomer
(5-CR110-PEO₁₂, SE)

CATALOG # 90086

MOLECULAR INFORMATION: C₅₂H₇₀N₄O₂₀
MWt: 1071.13

**PROPERTIES:**

Color & Form Orange red solid
Purity ≥ 70% reactivity by TLC (remaining amount is free acid)
Solubility Soluble in DMSO or DMF
Absorption/Emission 502/524 nm (MeOH)
Extinction Coefficient >65,000

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

APPLICATION: 5-CR110-PEO₁₂, SE has a long, water-soluble PEG spacer between the green fluorescent rhodamine dye and the amine-reactive SE group. See Figure 7.5 for spectra and Figure 7.6 for photostability comparison with other dyes.

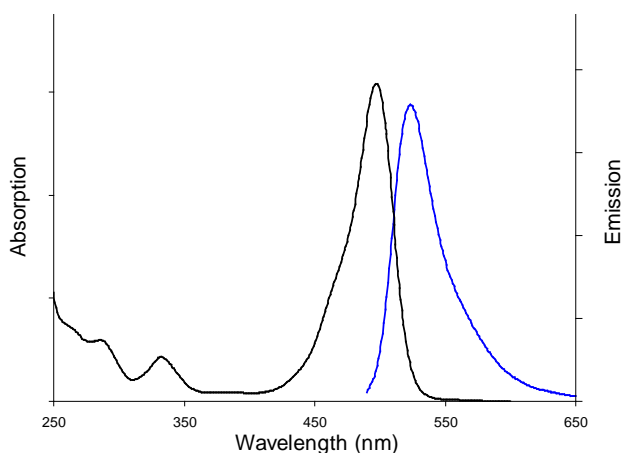


Figure 7.5 Absorption and emission spectra of CR110 and its derivatives in pH7 buffer.

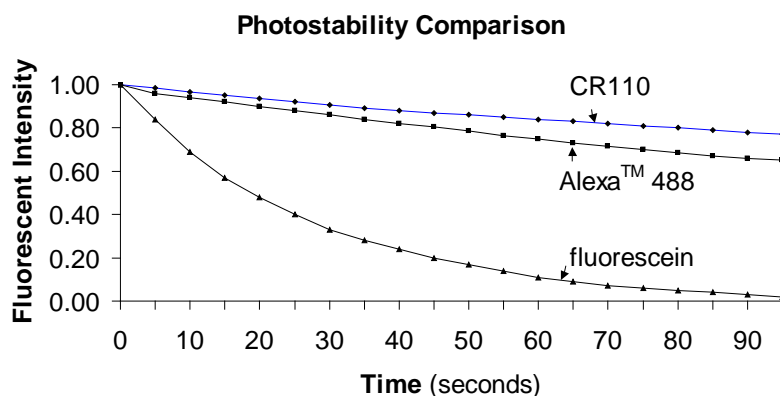


Figure 7.6 Photostability comparison of carboxyrhodamine 110-, fluorescein- and Alexa™ 488. Carboxyrhodamine 110 is more photostable than fluorescein and Alexa™ 488.

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.*