

PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: CFQ520 succinimidyl ester**CATALOG #:** 80500**MOLECULAR WEIGHT:** 675**PROPERTIES:**

| | |
|-------------------------------|-------------------------|
| Color & Form | Orange red solid |
| Purity | ≥ 95% by TLC |
| Solubility | Soluble in DMF and DMSO |
| Absorption/Emission | 521 nm / --- (MeOH) |
| Extinction Coefficient | 65,000 |

STORAGE AND HANDLING: Store at -20°C. Protect material from light, especially in solution.**APPLICATION:** The amine-reactive quencher, CFQ520 succinimidyl ester, has a broad and intense absorption (approximately 520 nm) but no fluorescence, making it useful as an acceptor in fluorescence resonance energy transfer (FRET) applications.**TOXICITY:** Unknown**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.*