



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

CFQ520 succinimidyl ester

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.



Product attributes

Product Description

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.

This datasheet was generated on January 16, 2026 at 09:31:01 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/cfq520-succinimidyl-ester/>