Call us: 800-304-5357 Email: btinfo@biotium.com

CFQ520 maleimide

The thiol-reactive quencher, CFQ520 maleimide, 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 thiol-reactive quencher, CFQ520 maleimide, 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 November 4, 2025 at 08:57:25 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/cfq520-maleimide/