



Glowing products for science

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

## 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 June 3, 2026 at 06:49:34 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/cfq520-maleimide/>