

NFQ1

NFQ1, a non-fluorescent quencher, is used for a new type of One Sample Fluorescence Resonance Energy Transfer (OS-FRET) method employing a novel, non-fluorescent methanethiosulfonate-linked acceptor that can be reversibly coupled to a target sulfhydryl residue via a disulfide bond.

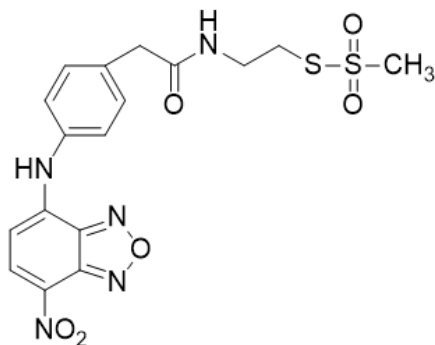


Product attributes

Product Description

NFQ1, a non-fluorescent quencher, is used for a new type of One Sample Fluorescence Resonance Energy Transfer (OS-FRET) method employing a novel, non-fluorescent methanethiosulfonate-linked acceptor that can be reversibly coupled to a target sulfhydryl residue via a disulfide bond. After the quenched donor emission is quantitated the acceptor is removed by reduction, enabling measurement of unquenched donor emission in the same sample.

- λ_{Abs} : 474 nm
- Orange solid soluble in DMSO at 20 mM
- Store desiccated at -20 °C and protect from light
- $\text{C}_{17}\text{H}_{17}\text{N}_5\text{O}_6\text{S}_2$
- MW: 451



References

Biochemistry 2010 doi: 10.1021/bi101188b

This datasheet was generated on December 22, 2025 at 12:22:48 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/nfq1/>