

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.

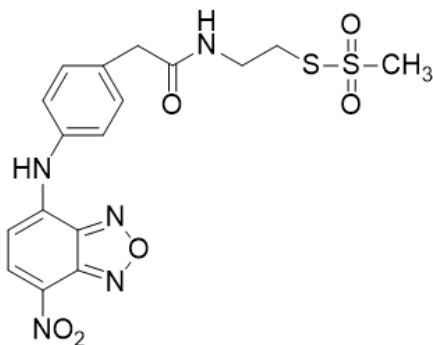


Product attributes

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
- $C_{17}H_{17}N_5O_6S_2$
- MW: 451



References

Biochemistry 2010 doi: 10.1021/bi101188b

This datasheet was generated on June 24, 2026 at 08:20:11 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/nfq1/>