

Product Information

NFQ1

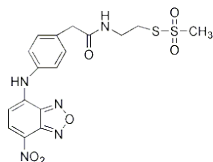
Catalog Number: 80501

Unit Size: 5 mg

Molecular Information

$C_{17}H_{17}N_5O_6S_2$

Molecular weight: 451



Properties

Color and form: Orange solid

Solubility: 20 mM in DMSO

Purity: >95% by TLC

Absorption: 474 nm

Extinction coefficient: 26,297

Storage and Handling

Store desiccated at -20°C and protect from light. Product is stable for at least 6 months from date of receipt when stored as recommended.

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.

References

Erbse, A.H. et al. (2011). *Biochemistry* 50 (4):451-7.

Please visit our website at www.biotium.com for information on our life science research products, including environmentally friendly EvaGreen® qPCR master mixes, fluorescent CF™ dye antibody conjugates and reactive dyes, apoptosis reagents, fluorescent probes, and kits for cell biology research.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.