MTS-ATFB

2-((4-Azido-2,3,5,6-tetrafluorobenzoyl)amino)ethyl methanethiosulfonate is a cross-linker containing a thiol-reactive MTS group and a photoaffinity label 4-azido-2,3,5,6-tetrafluorobenzene (ATFB). ATFB is known to be a superior photoaffinity label that undergoes highly efficient C-H bond insertion on UV photolysis.



Product attributes

Product Description

2-((4-Azido-2,3,5,6-tetrafluorobenzoyl)amino)ethyl methanethiosulfonate is a cross-linker containing a thiol-reactive MTS group and a photoaffinity label 4-azido-2,3,5,6-tetrafluorobenzene (ATFB). ATFB is known to be a superior photoaffinity label that undergoes highly efficient C-H bond insertion on UV photolysis.

- White solid soluble in DMF or DMSO
- Store at -20°C and protect from light
- C₁₀H₈F₄N₄O₃S₂
- MW: 372.32

References

1. J Org Chem 55, 3640 (1990).

This datasheet was generated on November 25, 2025 at 07:42:35 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/2-4-azido-2356-tetrafluorobenzoylaminoethyl-methanethiosulfonate/