

N-(6-Aminohexyl)maleimide, trifluoroacetic acid salt

N-(6-Aminohexyl)maleimide is a cross-linking reagent that can react with an amine-reactive group and a thiol group.

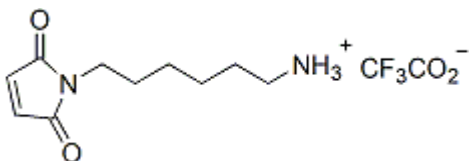


Product attributes

Product Description

N-(6-Aminohexyl)maleimide is a cross-linking reagent that can react with an amine-reactive group and a thiol group.

- White solid soluble in DMSO, DMF, or H₂O
- Store at -20 °C
- C₁₂H₁₇F₃N₂O₄
- MW: 310



This datasheet was generated on June 22, 2026 at 11:22:21 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/n-6-aminohexylmaleimide-trifluoroacetic-acid-salt/>