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

nicture 686

Product attributes

Call us: 800-304-5357 Email: btinfo@biotium.com

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

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