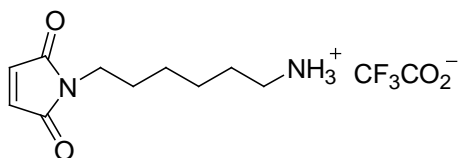


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: N-(6-aminohexyl)maleimide, trifluoroacetic acid salt**CATALOG #:** 91051**MOLECULAR INFORMATION:** C₁₂H₁₇F₃N₂O₄
Mwt: 310.27**PROPERTIES:**

Color & Form	White solid.
Purity	≥ 95% by TLC
Solubility	Soluble in DMF, DMSO or water
$\lambda_{ex} \backslash \lambda_{em}$	N/A
Extinction Coefficient	N/A

STORAGE AND HANDLING: Store desiccated at -20 °C and protect from light especially when in solution.**APPLICATION:** N-(6-aminohexyl)maleimide is a bifunctional cross-linker.**TOXICITY:** Unknown.

FIRST AID:	Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.
-------------------	---

Disclaimer: <i>Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.</i>
--