



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

N-(5-Aminopentyl)maleimide, trifluoroacetic acid salt

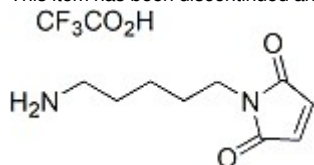
This item has been discontinued and replaced by N-(6-aminoheptyl)maleimide ([91051](#)).



Product attributes

Product Description

This item has been discontinued and replaced by N-(6-aminoheptyl)maleimide ([91051](#)).



This datasheet was generated on June 4, 2026 at 12:33:11 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/n-5-aminopentylmaleimide-trifluoroacetic-acid-salt/>