



Glowing products for science

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

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

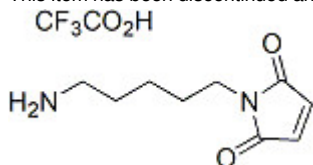
This item has been discontinued and replaced by N-(6-aminoheptyl)maleimide ([91051](#)).



Product attributes

Product Description

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



This datasheet was generated on December 16, 2025 at 08:17:55 PM. Visit product page to check for updated information before use.

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