



Glowing products for science

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

Aniline, 10X in Acetate Buffer

Aniline, 10X in acetate buffer is designed to use as catalytic reagent for the optimal coupling of protein carbonyls with aminoxy compounds.



Product attributes

Product Description

Aniline, 10X in acetate buffer is designed to be used as a catalytic reagent for the optimal coupling of protein carbonyls with aminoxy compounds.

This datasheet was generated on January 12, 2026 at 08:13:51 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/aniline-10x-acetate-buffer/>