

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 aminooxy 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 aminooxy compounds.

This datasheet was generated on November 14, 2025 at 03:12:39 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/aniline-10x-acetate-buffer/