5-Aminoallyl-dUTP (AA-dUTP), 10 mM solution

alidum

CAS number 116840-18-7

Product attributes

Call us: 800-304-5357 Email: btinfo@biotium.com

5-Aminoallyl-dUTP (AA-dUTP) can be enzymatically incorporated into DNA. The resulting amine-containing DNA can be subsequently labeled with a fluorescent dye, biotin or other haptens via conventional peptide coupling method.

Product Description

5-Aminoallyl-dUTP (AA-dUTP) can be enzymatically incorporated into DNA. The resulting amine-containing DNA can be subsequently labeled with a fluorescent dye, biotin or other haptens via conventional peptide coupling method (1,2). This two-step method for labeling nucleic acids is considerably more economical than the one-step method using a prelabeled dUTP. AA-dUTP is also available in lyophilized form (catalog no. 40020-1).

- 10 mM solution in TE buffer
- Store at -20°C
- C₁₂H₁₇N₃Na₃O₁₄P₃
- MW: 589.17
- [116840-18-7]

References

- 1. Biol Chem Hoppe Seyler 371, 953 (1990).
- 2. Biotechniques 28, 518 (2000).

This datasheet was generated on November 4, 2025 at 03:24:10 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/5-aminoallyl-dutp10-mm-in-te-buffer-aa-dutp/