

5-Aminoallyl-dUTP (AA-dUTP), Lyophilized Powder

Product attributes

CAS number 116840-18-7

Call us: 800-304-5357

5-Aminoallyl-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.

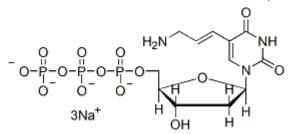
Product Description

5-Aminoallyl-dUTP (AA-dUTP) in lyophilized form suitable for long-term storage.

5-Aminoallyl-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.

- White solid soluble in water
- Store at -20°C
- C₁₂H₁₇N₃Na₃O₁₄P₃
- MW: 589.17
- [116840-18-7]

AA-dUTP is also available as a 10 mM solution in TE buffer (catalog no. 40020).



References

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

This datasheet was generated on November 25, 2025 at 11:30:23 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/5-aminoallyl-dutp-lyophilized-powder-aa-dutp/