

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



Product attributes

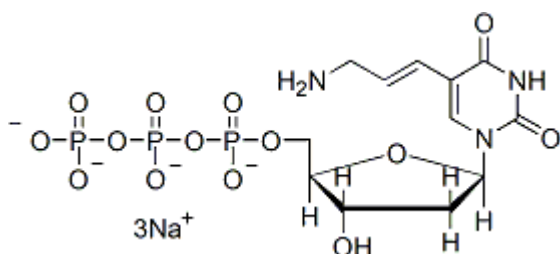
CAS number	116840-18-7
------------	-------------

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 June 3, 2026 at 08:07:27 PM. 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/>