

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

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 attributes

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

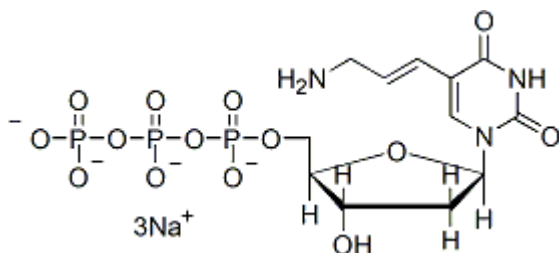
Product Description

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
- $C_{12}H_{17}N_3Na_3O_{14}P_3$
- 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 June 10, 2026 at 08:57:26 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/5-aminoallyl-dutp-lyophilized-powder-aa-dutp/>