

5-Aminoallyl-UTP, Sodium Salt, 10 mM Solution



Product attributes

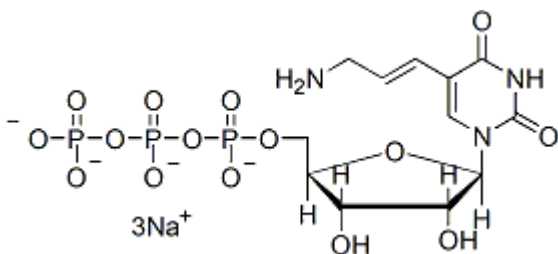
CAS number	112131-73-4
------------	-------------

5-Aminoallyl-UTP can be enzymatically incorporated into RNA for subsequent labeling with an amine-reactive fluorescent dye, biotin, or other hapten.

Product Description

5-Aminoallyl-UTP (5-(3-Aminoallyl)uridine-5'-triphosphate, trisodium salt; AA-UTP) can be enzymatically incorporated into RNA. The resulting amine-containing RNA can be subsequently labeled with a fluorescent dye, biotin, or other hapten via conventional peptide coupling methods. Biotium also offers AA-UTP in lyophilized form (catalog no. [40021-1](#)).

- 10 mM solution in TE buffer
- Store at -20°C
- $C_{12}H_{17}N_3Na_3O_{15}P_3$
- MW: 605.17
- [112131-73-4]



References

1. PNAS 90, 4206 (1993).
2. J Clin Microbiol 29, 583 (1991).
3. Histochemistry 93, 191 (1989).

This datasheet was generated on June 9, 2026 at 08:16:19 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/5-aminoallyl-utp-sodium-salt-10-mm-in-te-buffer-5-3-aminoallyluridine-5-triphosphate-trisodium-salt-aa-utp/>