

Biotin-11-UTP

Biotin-11-UTP can be enzymatically incorporated into RNA. The number 11 is the number of atoms in the linker between biotin and UTP.

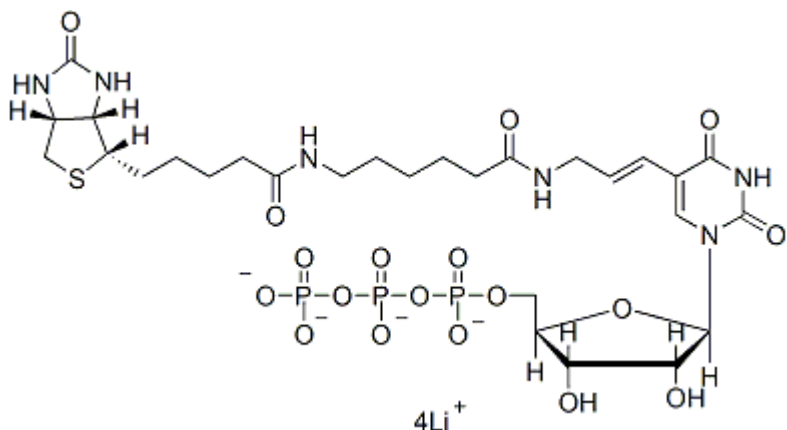


Product attributes

Product Description

Biotin-11-UTP (biotin-11-uridine-5'-triphosphate, tetralithium salt) can be enzymatically incorporated into RNA. The number 11 is the number of atoms in the linker between biotin and UTP.

- Supplied in pH 7.5 Tris-HCl buffer
- Store at -20 °C
- $C_{28}H_{41}N_6O_{18}P_3SLi_4$
- MW: 902.5



This datasheet was generated on August 10, 2026 at 09:53:31 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/biotin-11-utp-10-mm-in-ph-7-5-tris-hcl-buffer/>