

Biotin-16-UTP

Biotin-16-UTP can be enzymatically incorporated into RNA. The number 16 indicates the number of atoms between biotin and UTP.

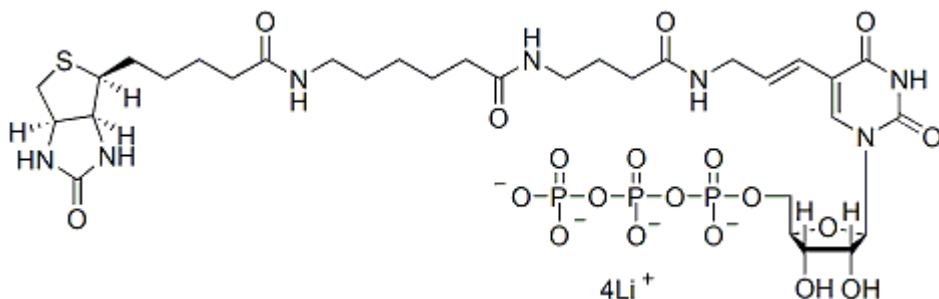


Product attributes

Product Description

Biotin-16-UTP (Biotin-16-uridine-5'-triphosphate, tetralithium salt) can be enzymatically incorporated into RNA. The number 16 indicates the number of carbons between biotin and UTP.

- Supplied in pH 7.5 Tris-HCl buffer
- $C_{32}H_{48}N_7O_{19}P_3SLi_4$
- MW: 987.51
- Store at $-20^{\circ}C$



This datasheet was generated on January 1, 2026 at 11:29:46 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/biotin-16-utp-10-mm-in-ph-7-5-tris-hcl-buffer-biotin-16-uridine-5-triphosphate-tetralithium-salt/>