Call us: 800-304-5357 Email: btinfo@biotium.com

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
- C₃₂H₄₈N₇O₁₉P₃SLi₄
- MW: 987.51
- Store at -20°C

This datasheet was generated on November 2, 2025 at 07:41:01 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/