

Biotin-11-CTP, 10 mM Solution

Biotin-11-CTP (Biotin-11-cytidine-5'-triphosphate, tetralithium salt) can be enzymatically incorporated into RNA.

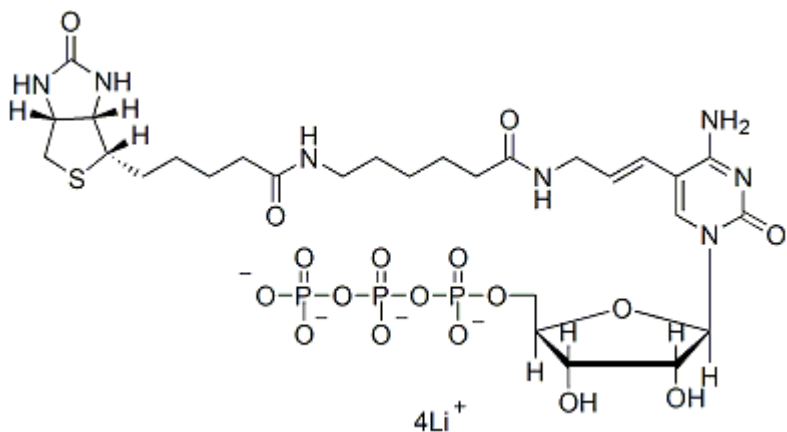


Product attributes

Product Description

Biotin-11-CTP (Biotin-11-cytidine-5'-triphosphate, tetralithium salt) can be enzymatically incorporated into RNA

- 10 mM solution in pH 7.5 Tris-HCl buffer
- Store at -20°C
- $C_{28}H_{42}N_7O_{17}P_3SLi_4$
- MW: 901.42



This datasheet was generated on June 10, 2026 at 10:29:27 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/biotin-11-ctp-10-mm-in-ph-7-5-tris-hcl-buffer-biotin-11-cytidine-5-triphosphate-tetralithium-salt/>