Burd of all 2

Biotin-11-CTP, 10 mM Solution

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



Product attributes

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

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
- C₂₈H₄₂N₇O₁₇P₃SLi₄
- MW: 901.42

This datasheet was generated on November 2, 2025 at 09:10:07 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/