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

## **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<sub>28</sub>H<sub>41</sub>N<sub>6</sub>O<sub>18</sub>P<sub>3</sub>SLi<sub>4</sub>
- MW: 902.5

This datasheet was generated on November 3, 2025 at 12:11:33 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/biotin-11-utp-10-mm-in-ph-7-5-tris-hcl-buffer/">https://biotium.com/product/biotin-11-utp-10-mm-in-ph-7-5-tris-hcl-buffer/</a>