

## Biotin-20-UTP, 10 mM Solution

This product is similar to biotin-11-UTP (40033) and biotin-16-UTP (40023) except that it has a longer and water soluble spacer group, which should facilitate interaction between the biotin group and avidin or streptavidin.

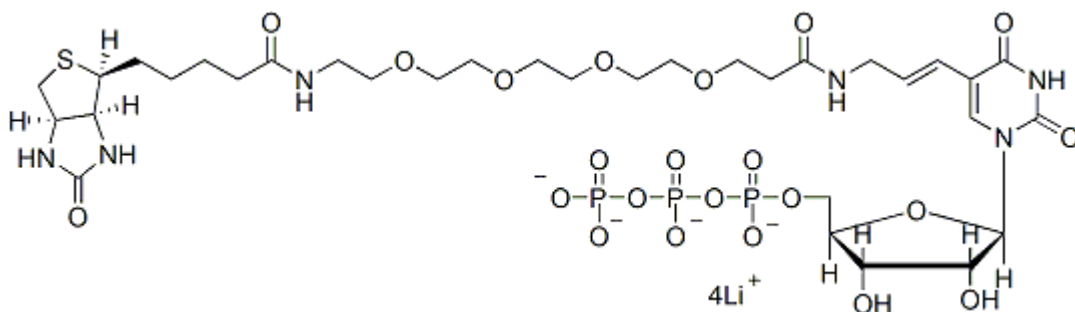


### Product attributes

### Product Description

This product is similar to biotin-11-UTP ([40033](#)) and biotin-16-UTP ([40023](#)) except that it has a longer and water soluble spacer group, which should facilitate interaction between the biotin group and avidin or streptavidin.

- 10 mM solution in pH 7.5 Tris-HCl buffer
- Store at -20 °C
- C<sub>33</sub>H<sub>51</sub>N<sub>6</sub>O<sub>22</sub>P<sub>3</sub>SLi<sub>4</sub>
- MW: 1036.26



This datasheet was generated on June 10, 2026 at 08:59:15 PM. Visit product page to check for updated information before use.  
 Product link: <https://biotium.com/product/biotin-20-utp-10-mm-in-ph-7-5-tris-hcl-buffer/>