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

Call us: 800-304-5357

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
- C33H51N6O22P3SLi4
- MW: 1036.26

This datasheet was generated on December 2, 2025 at 05:39:38 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/