

Biotin-20-dUTP, 1 mM Solution

This product is similar to biotin-11-dUTP (40029) and biotin-16-dUTP (40022) except that it has a longer and water-soluble spacer group, which should facilitate interaction between the biotin group and avidin or streptavidin.

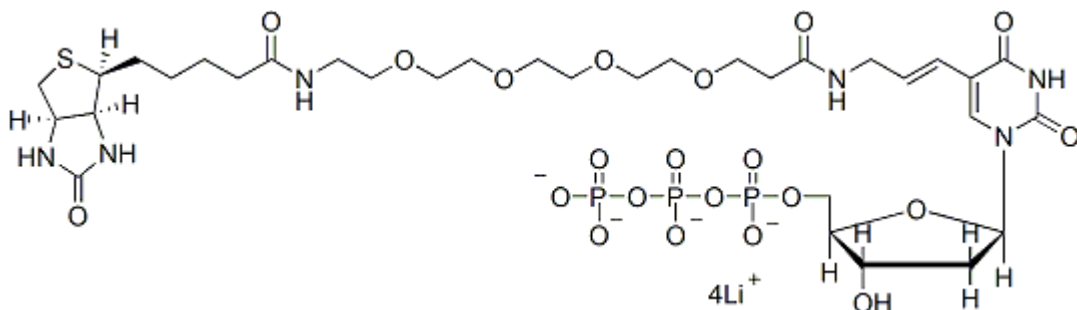


Product attributes

Product Description

This product is similar to biotin-11-dUTP ([40029](#)) and biotin-16-dUTP ([40022](#)) except that it has a longer and water-soluble spacer group, which should facilitate interaction between the biotin group and avidin or streptavidin. Biotin-20-dUTP is also available as a lyophilized solid (catalog no. [40030-1](#)).

- 1 mM solution in pH 7.5 Tris-HCl buffer
- Store at -20 °C
- C₃₃H₅₁N₆O₂₁P₃S Li₄
- MW: 1020.54



This datasheet was generated on June 11, 2026 at 08:25:52 PM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/biotin-20-dutp-1-mm-in-ph-7-5-tris-hcl-buffer/>