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

Call us: 800-304-5357

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 November 25, 2025 at 11:35:35 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/