

Biotin-20-dUTP, Lyophilized Powder

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.



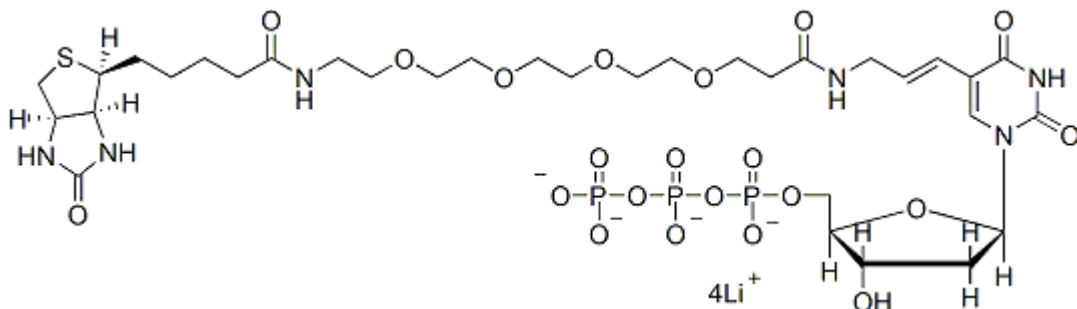
Product attributes

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.

- Lyophilized solid suitable for long term storage; contains lyophilized TE buffer
- Store at -20°C
- $C_{33}H_{51}N_6O_{21}P_3SLi_4$
- MW: 1020.54

Biotin-20-dUTP is also available as a 1 mM solution in Tris buffer ([40030](#)).



This datasheet was generated on January 12, 2026 at 01:47:19 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/biotin-20-dutp-lyophilized-powder/>