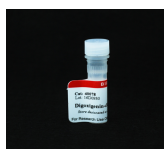


Digoxigenin-dUTP, Alkali Stable

Digoxigenin-dUTP can be enzymatically incorporated into DNA via nick translation, random priming, or 3'-end terminal labeling to synthesize labeled DNA probes for in-situ hybridization, microarray or blotting techniques. The digoxigenin labeled probe could be detected by using fluorescent labeled or enzyme labeled anti-digoxigenin antibody.

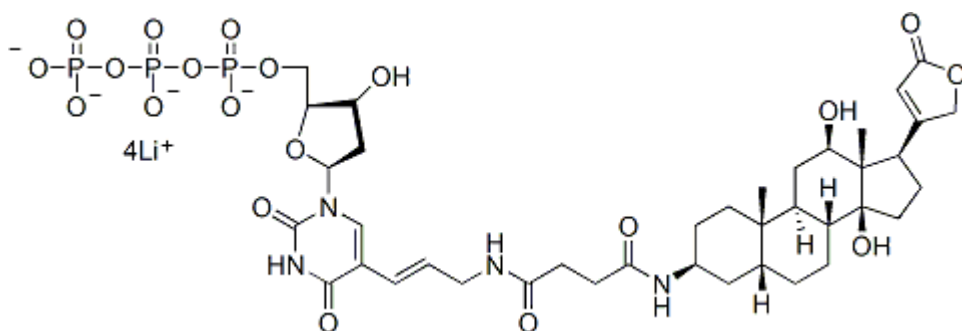


Product attributes

Product Description

Digoxigenin-dUTP can be enzymatically incorporated into DNA via nick translation, random priming, or 3'-end terminal labeling to synthesize labeled DNA probes for in-situ hybridization, microarray or blotting techniques. The digoxigenin labeled probe could be detected by using fluorescent labeled or enzyme labeled anti-digoxigenin antibody.

- Colorless solid soluble in water
- $C_{39}H_{53}N_4O_{20}P_3Li_4$
- MW: 1018.54
- Store at $-20^{\circ}C$



This datasheet was generated on January 16, 2026 at 02:36:38 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/digoxigenin-dutp-alkali-stable/>