## 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

Email: btinfo@biotium.com

## **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
- C39H53N4O20P3Li4
- MW: 1018.54
- Store at -20°C

This datasheet was generated on November 16, 2025 at 07:24:13 PM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/digoxigenin-dutp-alkali-stable/">https://biotium.com/product/digoxigenin-dutp-alkali-stable/</a>