

# HighFidelity Digoxigenin PCR Labeling Kit, S pack

HighFidelity Digoxigenin PCR Labeling Kit is designed to produce randomly Digoxigenin-modified DNA probes by PCR. Such probes are ideally suited for in situ hybridization and Northern Blot experiments. PCR-based labeling is superior to random-primed labeling by Klenow fragment if template amounts are limited or amplification of a specific DNA fragments is required. The labeling principle is similar to the underlying labeling principles of PCR DIG Labeling Mix (Roche) and PCR DIG Probe Synthesis Kit (Roche) with the exception that an alkali-stable version of DIG-11-dUTP is used (DIG label removal not feasible). DIG-11-dUTP is efficiently incorporated into DNA as substitute for its natural counterpart dTTP using an optimized reaction buffer and a High Fidelity Polymerase. 35 % DIG-11-dUTP substitution typically results in an optimal balance between reaction and labeling efficiency. Individual optimization of DIG-11-dUTP/dTTP ratio however, can easily be achieved with the single nucleotide format. The resulting Digoxigenin-modified DNA probe can subsequently be detected by HRP- or AP-modified Digoxigenin antibody. The kit contains sufficient reagents for 35 labeling reactions (S-Pack) or 175 labeling reactions (L-Pack) of 20 µl each (35 % DIG-11-dUTP substitution, 100 µM dATP/dGTP/dCTP, 65 µM dTTP, 35 µM DIG-11-dUTP).

## Ordering Information

[Order Online »](#)

JBS-APP-101-DIGX-S	35Reactions
--------------------	-------------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage      -20°C

Shipping                      Blue Ice

**Regulatory Status**      RUO - Research Use Only

## Product Details

Application Notes              Preparation of Digoxigenin-labeled DNA probes by PCR

Technical Info / Product Notes      For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)