

HighFidelity Biotin16 PCR Labeling Kit

HighFidelity Biotin16 PCR Labeling Kit is designed to produce randomly Biotin16-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. Biotin-16-dUTP is efficiently incorporated into DNA as substitute for its natural counterpart dTTP using an optimized reaction buffer and a High Fidelity Polymerase. 50 % Biotin-16-dUTP substitution typically results in an optimal balance between reaction and labeling efficiency. Individual optimization of Biotin-16-dUTP/dTTP ratio however, can easily be achieved with the single nucleotide format. The resulting Biotin16-modified DNA probe can subsequently be detected by HRP- or AP-modified Streptavidin. The kit contains sufficient reagents for 175 labeling reactions of 20 µl each (50 % Biotin16-dUTP substitution, 100 µM dATP/dGTP/dCTP, 50 µM dTTP, 50 µM Biotin16-dUTP).

Ordering Information

[Order Online »](#)

JBS-APP-101-BIO16	175Reactions
-------------------	--------------

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 Biotin16-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

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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