

# HighFidelity Fluorescein PCR Labeling Kit, S pack

HighFidelity Fluorescein PCR Labeling Kit is designed to produce randomly Fluorescein-modified DNA probes by PCR. Such probes are ideally suited for Fluorescence in situ hybridization (FISH) and Northern Blot experiments. PCR-based labeling is superior to random-primed labeling with Klenow fragment if template amounts are limited or amplification of a specific DNA fragments is required. Amplification of probes up to 4kbp is feasible. Fluorescein-12-dUTP is efficiently incorporated into DNA as substitute for its natural counterpart dTTP using an optimized reaction buffer and a High Fidelity Polymerase blend consisting of Taq polymerase and a proofreading enzyme. 50 % Fluorescein-12-dUTP substitution typically results in an optimal balance between reaction and labeling efficiency. Individual optimization of Fluorescein-12-dUTP/dTTP ratio however, can easily be achieved with the single nucleotide format. The kit contains sufficient reagents for 10 labeling reactions (S-Pack) or 50 labeling reactions (L-Pack) of 20 µl each (50% Fluorescein-12-dUTP substitution, 100 µM dATP/dGTP/dCTP, 50 µM dTTP, 50 µM Fluorescein-12-dUTP).

## Ordering Information

[Order Online »](#)

JBS-APP-101-FAMX-S	10Reactions
--------------------	-------------

## 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 Fluorescein-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)