



# Label Your Own FISH Probes

## Nick Translation DNA Labeling System & Fluorescent dye-dUTPs

### Ensure Peak Labeling Performance for FISH Probe Synthesis

Fluorescent dye-dUTPs are well recognized as superior to analogous but cumbersome indirect two-step labeling methods. When coupled with the Nick Translation DNA Labeling System, this direct approach provides a simple and efficient method to label DNA for FISH, suitable for a wide range of molecular biology and cytogenetics applications.

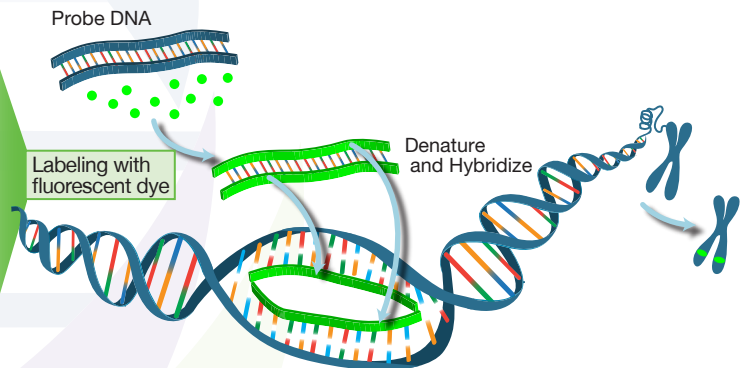
#### Nick Translation DNA Labeling System (ENZ-42910)

- Simple and efficient method for generating labeled DNA
- Allows the user to optimize incorporation and product size by adjusting the ratio of components
- Compatible with both biotin-labeled and fluorophore-labeled nucleotides

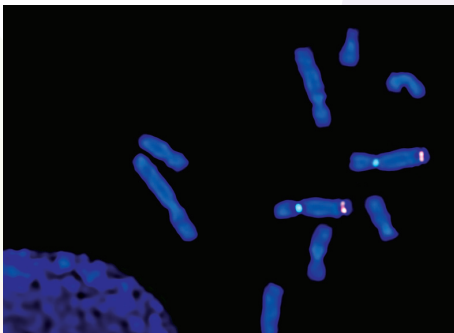
#### Fluorescent dye-dUTPs

- Eight distinct colors to choose from, spanning the visible light spectrum
- High signal intensity and good photostability
- Available in aqueous format or new lyophilized sizes

#### Labeling Probes for FISH Analysis



#### Total FISH Solution: Nick Translation Kit + Fluorescent dye-dUTPs



Composite fluorescent image using the Nick Translation Kit and BAC Target DNA labeled with Enzo Orange dUTP (TAMRA) and Centromere BAC Probe labeled with ENZO Green 496 dUTP (FITC). Metaphase chromosome spreads were counterstained with DAPI.

| PRODUCT NAME                         | PRODUCT # |
|--------------------------------------|-----------|
| Nick Translation DNA Labeling System | ENZ-42910 |
| Gold 550 dUTP                        | ENZ-42521 |
| Red 650 dUTP                         | ENZ-42522 |
| Green 496 dUTP                       | ENZ-42831 |
| Orange 552 dUTP                      | ENZ-42842 |
| Gold 525 dUTP                        | ENZ-42843 |
| Red 580 dUTP                         | ENZ-42844 |
| Green 500 dUTP                       | ENZ-42845 |
| Aqua 431 dUTP                        | ENZ-42853 |

Fuel Your Research:  
[www.enzolifesciences.com](http://www.enzolifesciences.com)

#### GLOBAL HEADQUARTERS

Enzo Life Sciences Inc.  
10 Executive Boulevard  
Farmingdale, NY 11735  
Phone: 800.942.0430  
Fax: 631.694.7501  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

#### EUROPE

Enzo Life Sciences (ELS) AG  
Industriestrasse 17  
CH-4415 Lausen, Switzerland  
Phone: +41 61 926 8989  
Fax: +41 61 926 8979  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)