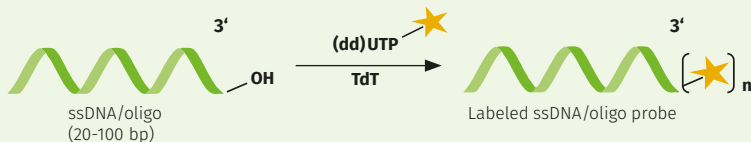
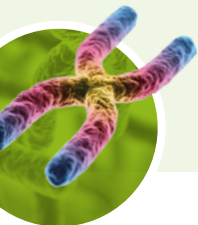


# 3' End Oligonucleotide Labeling Kits

for preparation of Biotin & Digoxigenin-labeled ssDNA/oligo probes



**TdT:** Terminal deoxynucleotidyl Transferase

$n = 1$  for Biotin-11-UTP

$n = 1 - 3$  (average) for Digoxigenin-11-ddUTP

★: Biotin or Digoxigenin

Kit	Cat.-No.	Label	# of incorporated label	Labeling principle similar to ...	Price (€)
BIO 3'-End Oligonucleotide Labeling Kit with Biotin-11-UTP	APP-001	Biotin	1-3	Biotin 3'-End DNA Labeling Kit (Pierce/ Thermo Scientific)	275,00
DIG 3'-End Oligonucleotide Labeling Kit with Digoxigenin-11-ddUTP	APP-002	Digoxigenin	1	DIG Oligonucleotide 3' End Labeling Kit, 2nd generation (Roche)	315,00



Biotin & Digoxigenin  
3' End Oligonucleotide  
Labeling Kits

# 3' End Oligonucleotide Labeling Kits for preparation of Biotin & Digoxigenin-labeled ssDNA/oligo probes

## Comparable labeling efficiency of BIO 3' End Oligonucleotide Labeling Kit (JBS) and Biotin 3' End DNA Labeling Kit (competitor P/TS)

### Labeling reaction:

5 pmol oligonucleotide (60 bp), 50 pmol UTP (negative control) or 50 pmol Biotin-11-UTP (Biotin labeling reaction)\*, 1x TdT Reaction Buffer\*, 20 U TdT\* ad 50 µl PCR-grade H<sub>2</sub>O. Incubation for 30 min at 37°C.

(\*derived from JBS or P/TS Kit)

**Dot Blot analysis of labeling efficiency:** Oligonucleotide sample dilution series (50–3.125 fmol) and subsequent Biotin detection with Streptavidin-Alkaline Phosphatase.

**P/TS:** Biotin 3' End DNA Labeling Kit (competitor P/TS)

**JBS:** BIO 3'-End Oligonucleotide Labeling Kit (JBS)

-: negative control (labeling reaction with UTP)

+: Biotin-labeled oligonucleotide (labeling reaction with Biotin-11-UTP)

