

**Methods for preparing labeled hybridization probes**

The following chart summarizes the common methods of DNA and RNA probe generation and their relative advantages. These methods are most commonly used with our <sup>32</sup>P and <sup>35</sup>S nucleotides, but can also be utilized with <sup>3</sup>H nucleotides.

Method	Enzyme used	Specific activity (dpm/μg)	Features	Optimal nucleotides
Nick translation	DNA Pol I	10 <sup>8</sup> -10 <sup>9</sup>	-High specific activity -Trouble-free -Inexpensive *Hairpin loops of DNA can be formed	dCTP, dATP, dTTP
Random Priming	Klenow	10 <sup>8</sup> -10 <sup>9</sup>	-Highest specific activity *Need primer	dCTP, dATP, dTTP
Primer Extension	Klenow	10 <sup>7</sup> -10 <sup>8</sup>	-Partially single-stranded probe	dCTP, dATP, dTTP
3' End Labeling	Terminal transferase	10 <sup>7</sup>	-Highly specific -Oligo probes	3'-deoxy-ATP (cordycepin)
5' End Labeling	T4 Polynucleotide Kinase (T4 PNK)	10 <sup>7</sup>	-Highly specific -Oligo probes	gamma phosphate labeled-ATP
Fill-in	Klenow	10 <sup>7</sup>	-Highly specific -For 5' or 3' overhangs after restriction enzyme digest	dCTP, dATP, dTTP
T4 DNA labeling	T4 DNA Polymerase	10 <sup>7</sup>	End labeling for blunt ends. Can be made strand-specific by restriction enzyme digestion	dCTP, dATP, dTTP
RNA Probes	SP6 RNA polymerase	10 <sup>7</sup> -10 <sup>8</sup>	DNA/RNA has high binding coefficient	UTP, CTP, GTP
RNA Probes	T7 RNA polymerase	10 <sup>7</sup> -10 <sup>8</sup>	DNA/RNA has high binding coefficient	UTP, CTP, GTP
cDNA synthesis	Reverse transcriptase	10 <sup>7</sup>	High specificity for probing recombinants	dCTP, dATP, dTTP