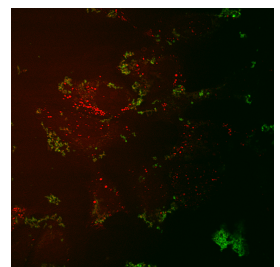
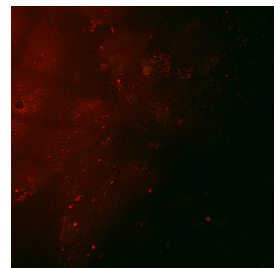
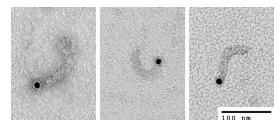
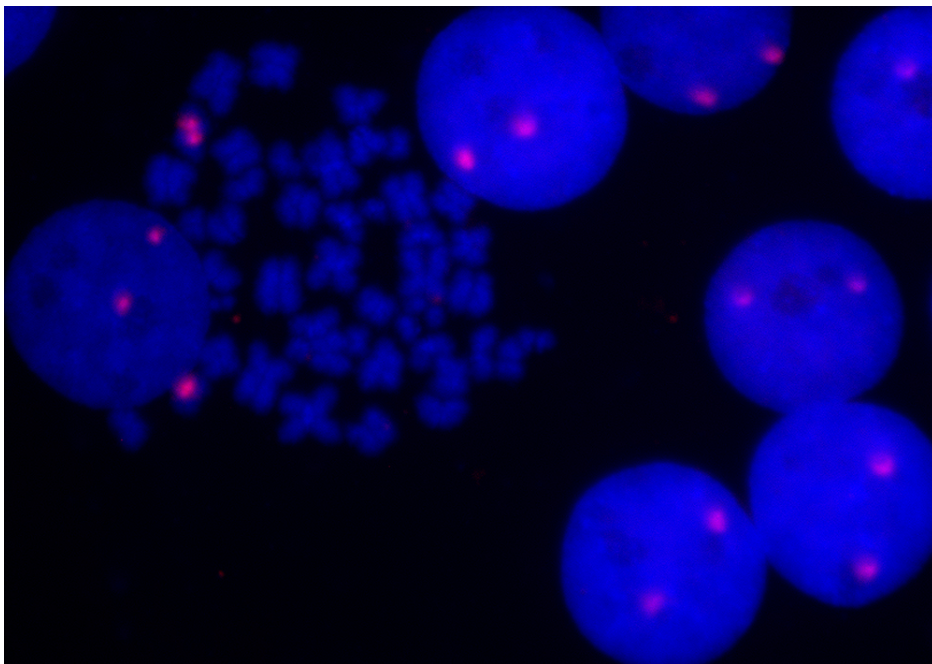


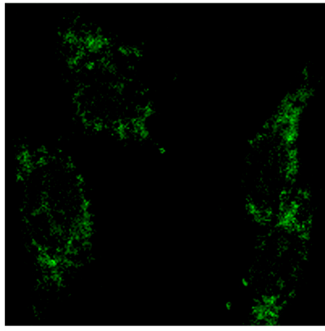


5' EndTag™ Nucleic Acid Labeling Kit

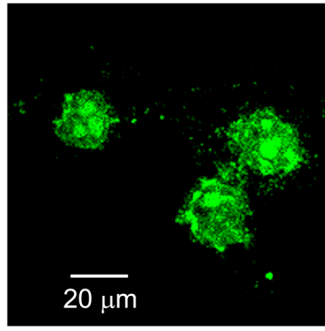
MB-9001

Product Images

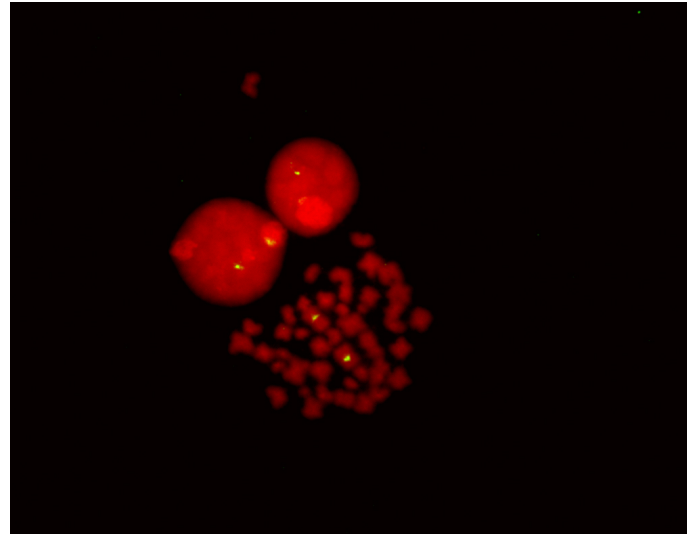




- energy
- cytosol



+ energy
+ cytosol



Short Description

End labeling is a favored method for applications in which an internal label might interfere with hybridization or sequence specific protein binding.

Features of the 5' EndTag DNA/RNA Labeling Kit:

- Enables simple and uniform labeling of the 5' end of a DNA, RNA or unmodified oligonucleotide molecule.
- Preserves the unmodified 3' end of the oligonucleotide (ideal for the labeling of PCR primers as a result).
- Compatible with a choice of diverse labels - fluorophores, biotin, haptens, or affinity ligands*

Additional Information

Unit Size	1 kit
Target for Labeling	DNA, Oligonucleotides, RNA
Required Reactive Group	5 '-OH group of nucleic acid
Tag/Group Incorporated	Thiol
Safety Title	NA

