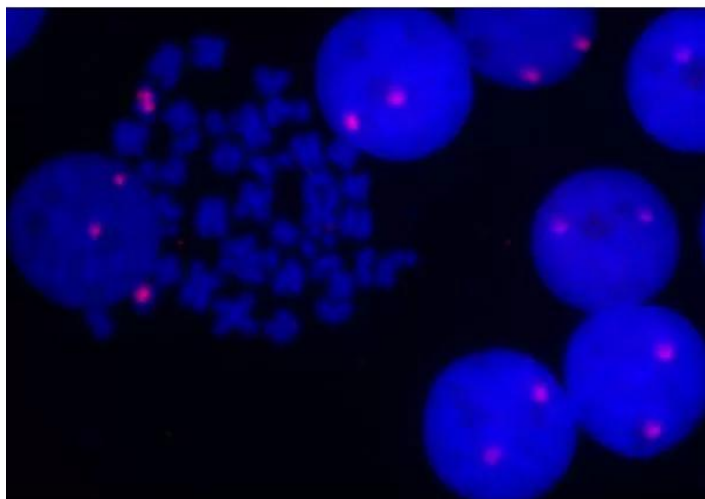




5' ENDTAG DNA/RNA LABELING KIT

SKU: MB-9001



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

SPECIFICATIONS

Unit Size

1 Kit

For research use only. Not intended for therapeutic or diagnostic use in animals or humans.



Species	DNA, Oligonucleotides, RNA
Tag/Group Incorporated	Thiol
Reactive Group	5'-OH group of nucleic acid
Safety Title	WARNING: NA -www.P65Warnings.ca.gov
Source Species	DNA, Oligonucleotides, RNA

KIT COMPONENTS

Materials suitable to perform 10 end labeling reactions of up to 0.6 nmols of 5' ends (e.g. about 5 µg of a 25 base oligo) per reaction. This includes:

- T4 polynucleotide kinase
- 10x reaction buffer
- ATPγS
- Precipitant
- Alkaline Phosphatase

***Note: Thiol-reactive labeling reagent is not included in the kit and must be purchased separately.** Labels suitable for this kit include:

- [Biotin \(Long Arm\) Maleimide](#)
- DNP Maleimide
- Fluorescein (Long Arm) Maleimide
- Fucose Maleimide
- Texas Red® Maleimide

Note regarding shipping: This item is shipped on wet ice and therefore requires additional shipping charges.

TECHNICAL INFORMATION

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

5' or 3' EndTag labeled nucleic acids can be used for applications such as DNA hybridization, PCR, *in situ* hybridization, the binding of capture probes to affinity matrices, or electrophoretic mobility shift assays (EMSA). Short oligonucleotides are labeled more efficiently with these

For research use only. Not intended for therapeutic or diagnostic use in animals or humans.



systems than with other methods. In addition, end labeling of oligonucleotides is an economical alternative to having labels inserted during synthesis.

Both the 5' EndTag and 3' EndTag Nucleic Acid Labeling Systems enable the covalent attachment of a variety of fluorescent dyes, haptens, or affinity tags to the respective ends of a nucleic acid using thiol-specific chemistry. Labels containing thiol-reactive groups (maleimides, iodoacetamides, etc.) can easily be incorporated.

Labeling time is about 1 hour with very little hands-on time.

5' EndTag Labeling Kit

The 5' EndTag System labels 5' ends of DNA, RNA, or unmodified oligonucleotides. This kit is ideal for labeling PCR primers because a label is attached only to the 5' end, leaving the 3' end available for polymerization.

Labeling with the 5' EndTag Kit is achieved in two steps:

1. T4 polynucleotide kinase transfers a thiophosphate from ATP γ S to the 5' -OH group of the nucleic acid. (5' phosphorylated ends are converted to 5' -OH groups with the included alkaline phosphatase).
2. Thiolated sample is coupled to a thiol-reactive label. The 5' EndTag Kit is designed to perform 10 labeling reactions of up to 0.6 nmols of 5' ends (e.g. about 5 μ g of a 25 base oligo) per reaction. Thiol-reactive label is not included in the kit and should be selected separately.

For research use only. Not intended for therapeutic or diagnostic use in animals or humans.

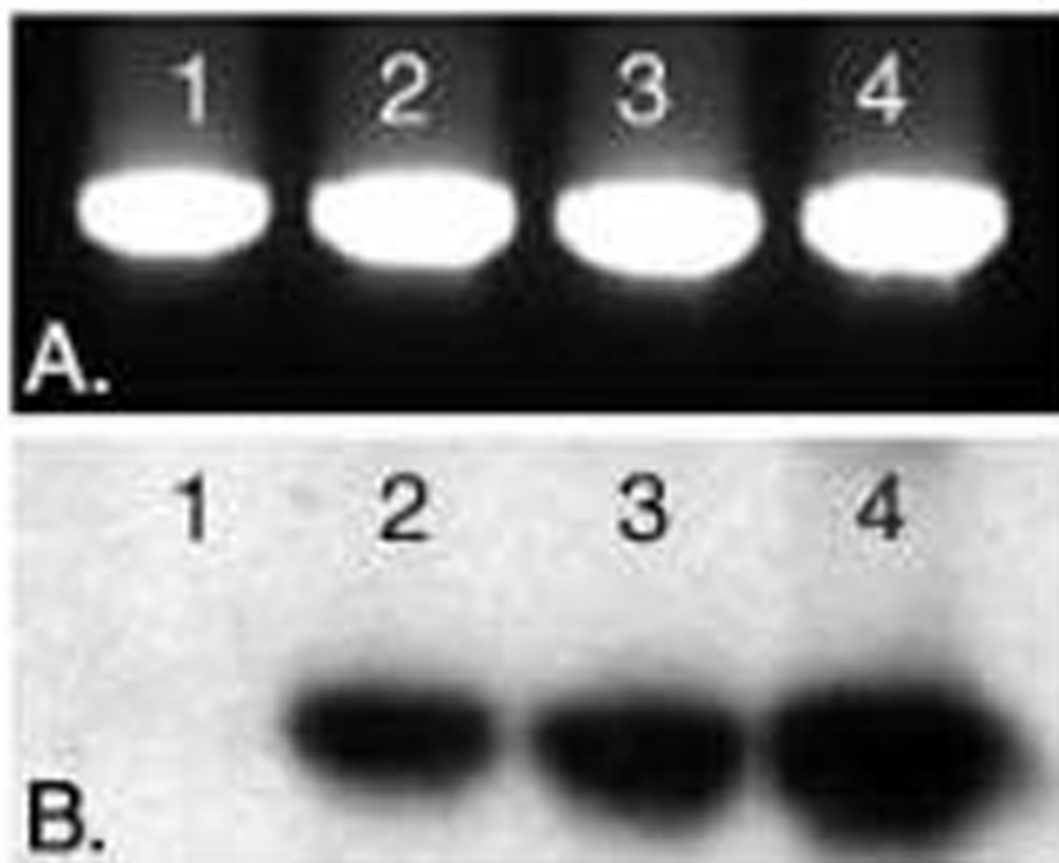


Figure 1: EtBr-stained gel of a 1002 bp PCR product amplified using:

Ln 1) Unlabeled primers

Ln 2) EndTag biotinylated forward primer and unlabeled reverse primer

Ln 3) Unlabeled forward primer and EndTag biotinylated reverse primer

Ln 4) EndTag biotinylated forward and reverse primers.

For research use only. Not intended for therapeutic or diagnostic use in animals or humans.

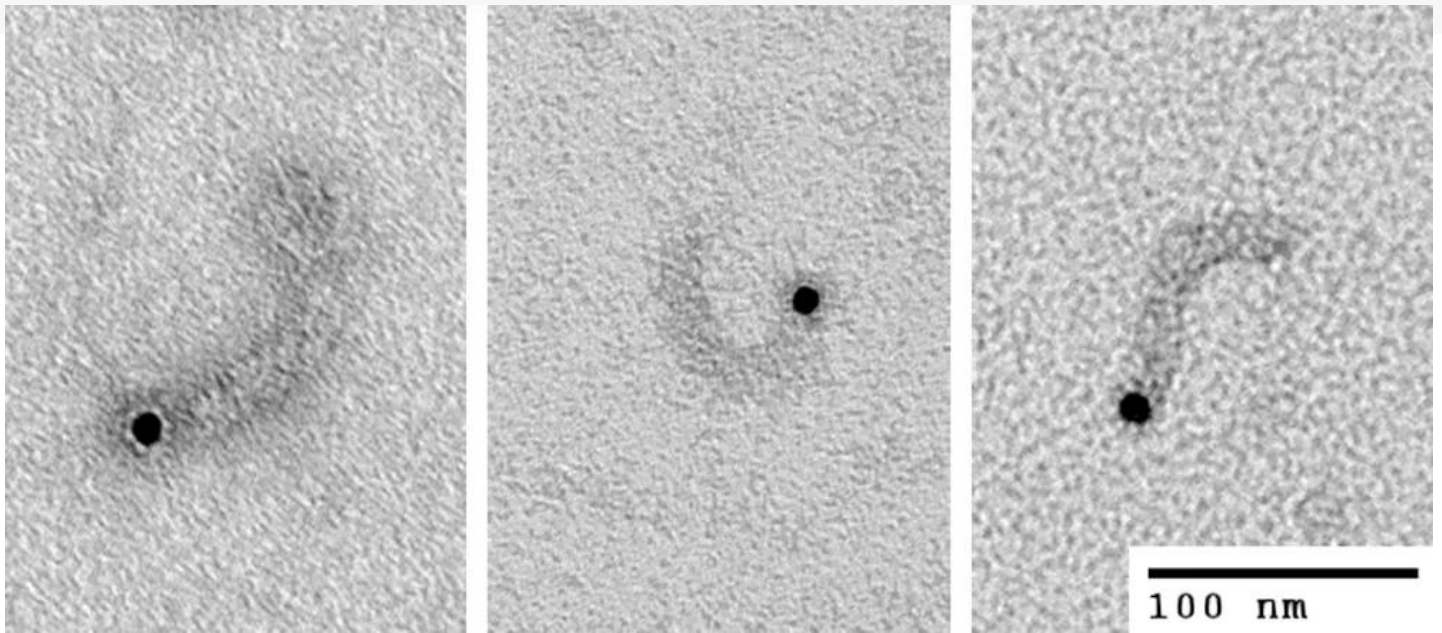
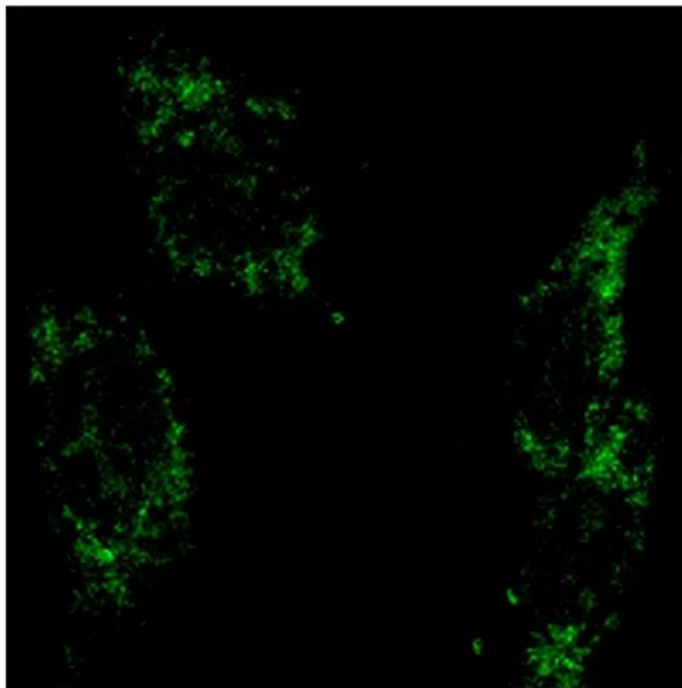
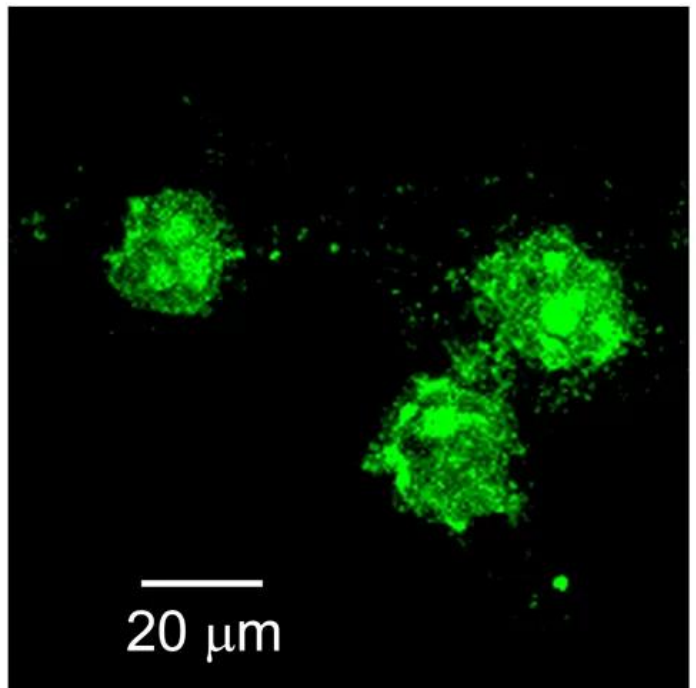


Figure 2: Transmission electron micrographs of molecules of influenza A viral ribonucleoprotein particles (vRNPs) labeled at the 5' end of the vRNA with biotin using 5' EndTag Kit, and further labeled with streptavidin gold. vRNPs from influenza A were purified according to Kemler et al., on a glycerol gradient with modifications as described in Wu et al. Courtesy of Drs. Winco WH Wu and Nelly Panté, University of British Columbia, Vancouver BC, Canada.

For research use only. Not intended for therapeutic or diagnostic use in animals or humans.



- energy
- cytosol



+ energy
+ cytosol

Figure 3: Nuclear import assay in digitonin-permeabilized HeLa cells of biotinylated vRNPs. vRNPs were labeled first at the 5' end of the vRNA with biotin using the 5' EndTag Kit and then with **Vector Fluorescein Streptavidin**. This allowed for direct fluorescence visualization of the vRNPs on a confocal fluorescence microscope. Nuclear import assays were carried out as described in Wu et al. The negative control consists of vRNPs added to the cells in the absence of energy and exogenous cytosol. In the presence of energy and cytosol, the fluorescein-labeled vRNPs successfully enter the nucleus, with a high degree of nucleolar staining. Courtesy of Drs. Winco WH Wu and Nelly Panté, University of British Columbia, Vancouver BC, Canada.

For research use only. Not intended for therapeutic or diagnostic use in animals or humans.

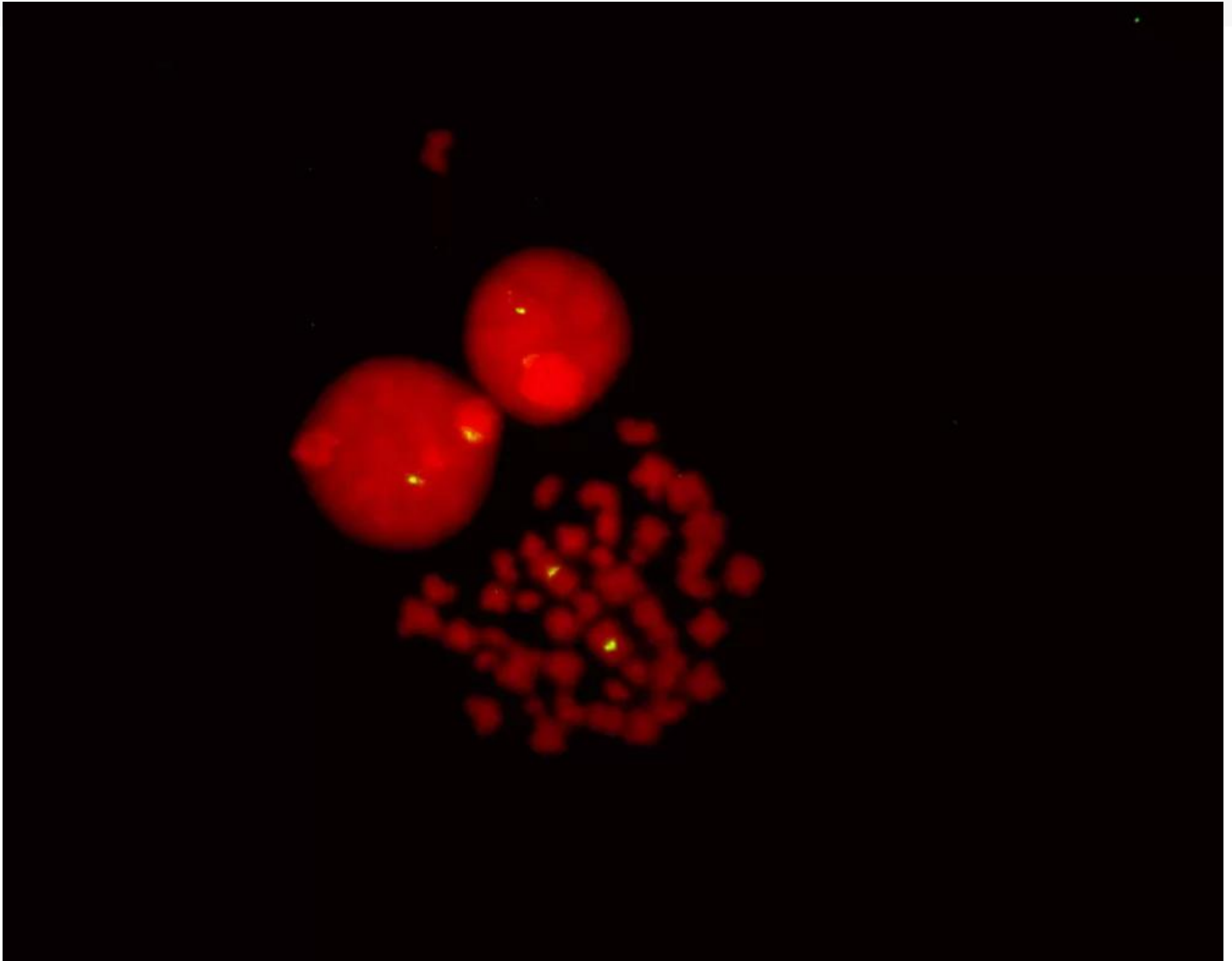


Figure 4: Fluorescence in situ hybridization of human chromosomes using 5' EndTag Fluorescein-labeled pUC1.77 detected directly and mounted in **Vectashield Mounting Media with PI**.

For research use only. Not intended for therapeutic or diagnostic use in animals or humans.

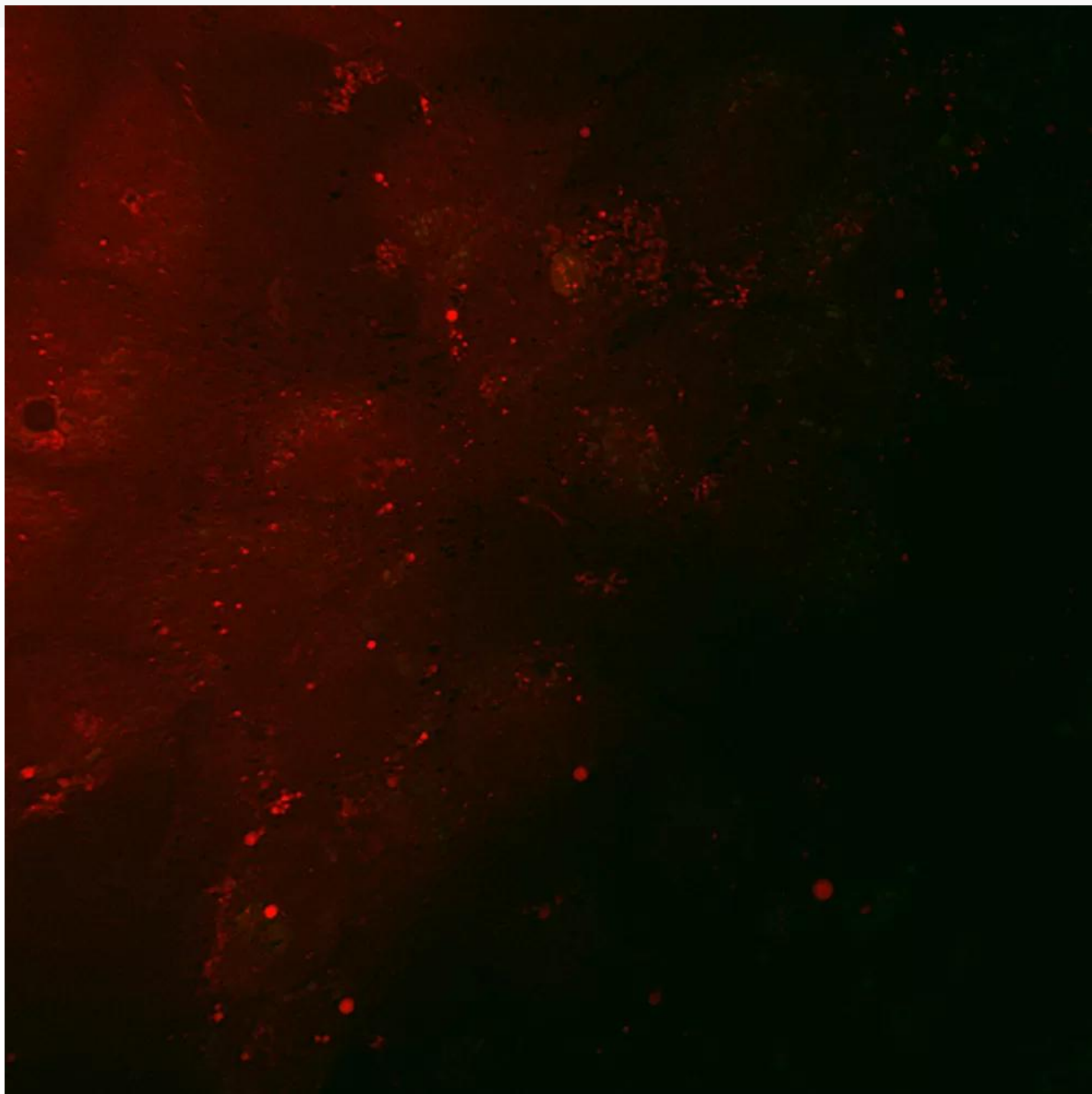


Figure 4: Coherent anti-Stokes Raman scattering (CARS) and two-photon fluorescence (TPF) imaging of fluorescein-labeled HCV RNA inside Huh-7 cells. HCV RNA in-vitro transcripts were labeled with a fluorescein-maleimide label using the 5' EndTag Nucleic Acid Labeling System. Huh-7 were transfected with 3 μ g of 5'-fluorescein-labeled HCV RNA and imaged at 4 h post-transfection. CARS and TPF of Mock transfected Huh-7 cells. Figure courtesy of Jennifer Haley,

For research use only. Not intended for therapeutic or diagnostic use in animals or humans.



Sylvie Belanger, Adrian Pegoraro, Albert Stolow, and John P. Pezacki, The Steacie Institute for Molecular Sciences, The National Research Council of Canada. Unpublished data.

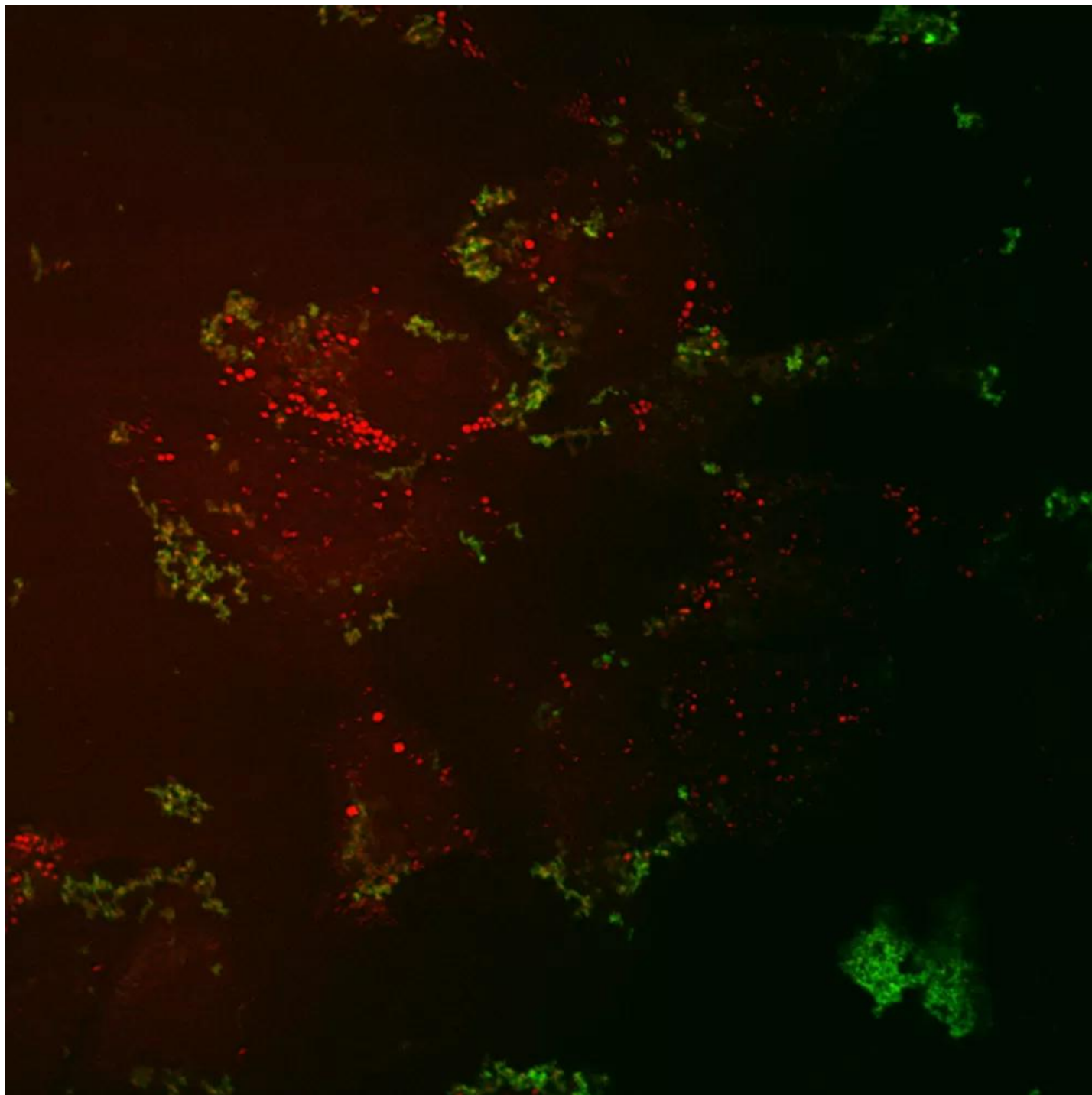


Figure 5: Coherent anti-Stokes Raman scattering (CARS) and two-photon fluorescence (TPF) imaging of fluorescein-labeled HCV RNA inside Huh-7 cells. HCV RNA in-vitro transcripts were

For research use only. Not intended for therapeutic or diagnostic use in animals or humans.



labeled with a fluorescein-maleimide label using the 5' EndTag Nucleic Acid Labeling System. Huh-7 were transfected with 3µg of 5'-fluorescein-labeled HCV RNA and imaged at 4 h post-transfection. Here, CARS and TPF of Huh-7 cells transfected with fluorescein-labeled HCV RNA. Figure courtesy of Jennifer Haley, Sylvie Belanger, Adrian Pegoraro, Albert Stolow, and John P. Pezacki, The Steacie Institute for Molecular Sciences, The National Research Council of Canada. Unpublished data.

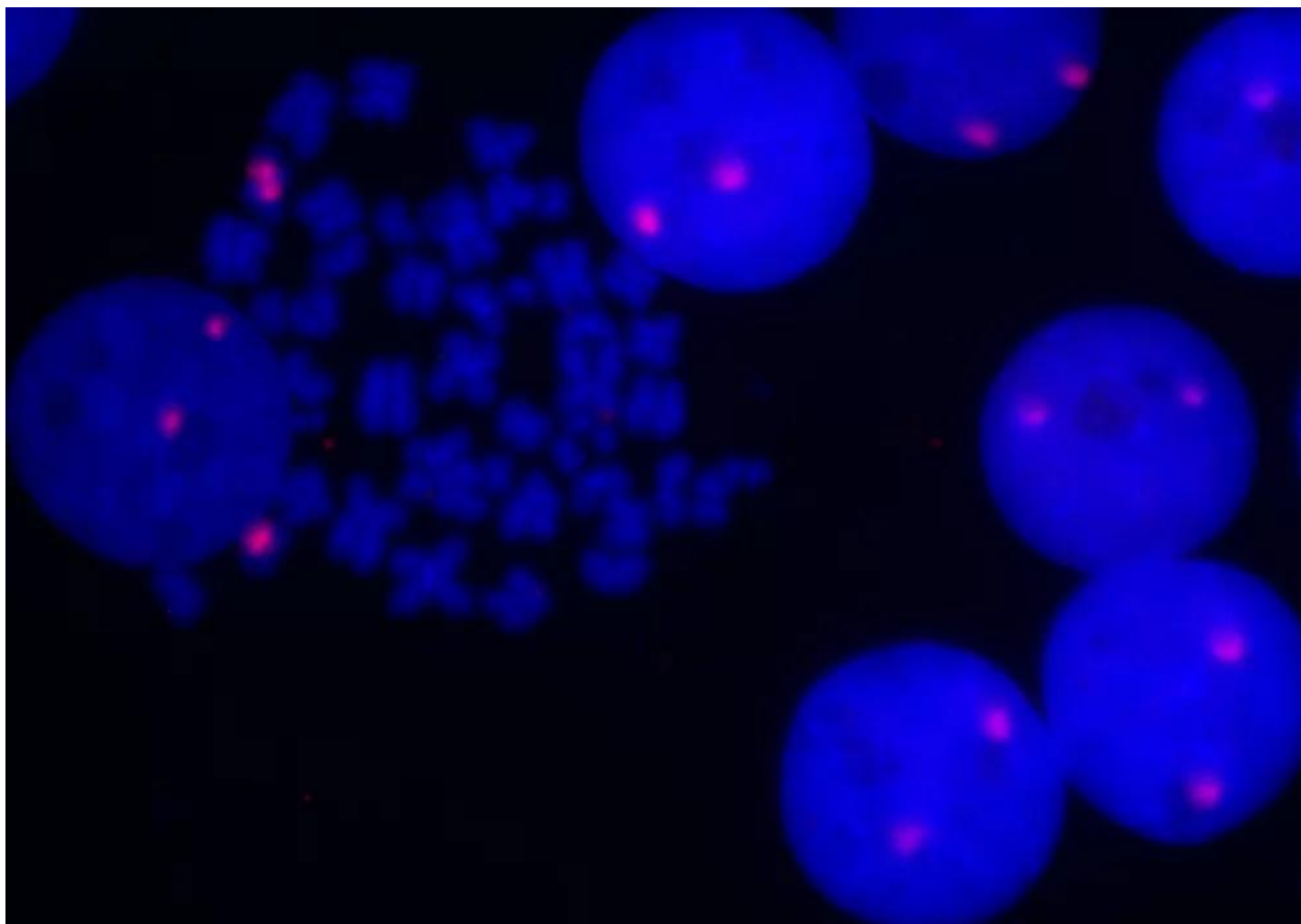


Figure 7: Fluorescence in situ hybridization of human chromosomes using 5' EndTag Texas Red-labeled pHuR 98 detected directly and mounted in **Vectashield Mounting Media with DAPI**.

CITATIONS

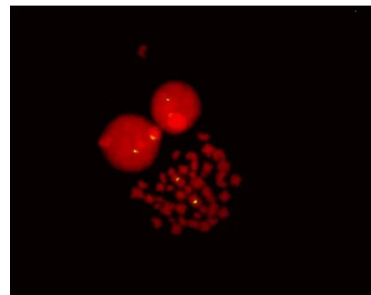
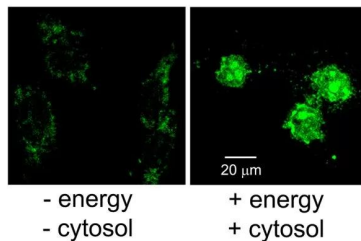
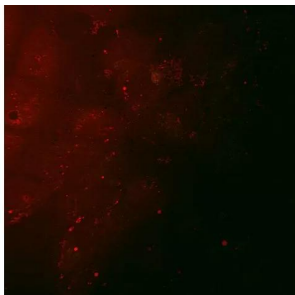
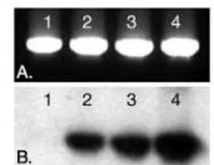
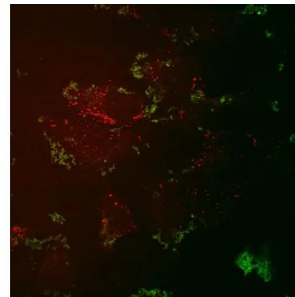
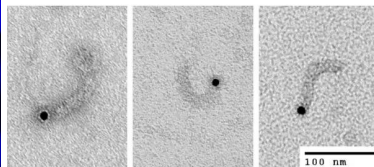
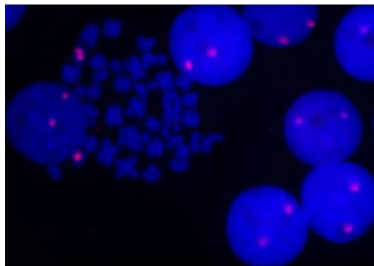
For research use only. Not intended for therapeutic or diagnostic use in animals or humans.



DOCUMENTS

- [User Guide](#)
- [Safety Data Sheet](#)
- [Download CoA](#)
- [Datasheet](#)

GALLERY IMAGES



For research use only. Not intended for therapeutic or diagnostic use in animals or humans.