

## ULYSIS<sup>®</sup> Nucleic Acid Labeling Kits

Table 1. Contents and storage

Material	Amount	Storage	Stability
ULS <sup>®</sup> labeling reagent (Component A)	1 vial for 20 labelings or 5 vials for 4 labelings each (for Cat. no. U21650)	<ul style="list-style-type: none"> <li>• ≤-20°C</li> <li>• Desiccate</li> <li>• Protect from light</li> </ul>	When stored as directed, kit components are stable for at least 6 months.
Dimethylformamide (DMF, 50% solution in water) or Dimethylsulfoxide (DMSO) (Component B)	200 µL		
Labeling buffer (Component C)	600 µL		
Deoxyribonuclease I (DNase I) (Component D)	100 µg		
DNase I storage buffer (Component E)	200 µL		
10X DNase I reaction buffer (Component F)	500 µL		
DNA from calf thymus (Component G)	100 µL of a 0.1 mg/mL solution in TE buffer		
Nuclease-free H <sub>2</sub> O (Component H)	5 mL		

**Number of assays:** Sufficient materials are supplied for 20 labelings of 1 µg DNA each.

## Introduction

ULYSIS<sup>®</sup> Nucleic Acid Labeling kits provide a non-enzymatic method for chemically labeling purine bases in nucleic acids with fluorescent dyes. The method, the Universal Linkage System (ULS<sup>®</sup>), is based on the use of a platinum dye complex that forms a stable adduct with the N7 position of guanine and, to a lesser extent, adenine bases in DNA, RNA, PNA, and oligonucleotides (Figure 1, page 2). The labeling reaction takes only 15 minutes and separation of the labeled nucleic acids from the unreacted ULS<sup>®</sup> complex can be accomplished through the use of a simple spin-column procedure (Figure 2, page 2). The ULS<sup>®</sup> method has been used to prepare labeled probes for dot, Southern and Northern blot analysis, RNA and DNA *in situ* hybridization, multicolor fluorescence *in situ* hybridization (mFISH), and comparative genome hybridization (CGH). The ULYSIS<sup>®</sup> kits allow researchers to label DNA with our exceptionally bright and photostable fluorescent dyes, including the proprietary Alexa Fluor<sup>®</sup> dyes (Table 2, page 2).

This user guide describes how to prepare fluorescent DNA hybridization probes optimized for chromosome *in situ* hybridization and dot blot hybridization. Suggestions for labeling RNA are also provided.

For Research Use Only. Not for use in diagnostic procedures.

Table 2. Spectral characteristics of the fluorescent dyes available in the ULYSIS® Nucleic Acid Labeling Kits

Cat. no.	Fluorescent Dye	$\lambda_{\text{max}}$ (nm) *	Em (nm) †	$\epsilon_{\text{dye}}$ (cm <sup>-1</sup> M <sup>-1</sup> ) ‡	CF <sup>260</sup> §	Spectrally Similar Dyes
U21650	Alexa Fluor® 488	492	520	62,000	0.30	Fluorescein (FITC)
U21652	Alexa Fluor® 546	555	570	104,000	0.21	Cy®3, tetramethylrhodamine (TRITC)
U21654	Alexa Fluor® 594	588	615	80,400	0.43	Texas Red®
U21660	Alexa Fluor® 647	650	670	239,000	0.00	Cy®5

\* Absorbance maximum for the fluorophore; † Emission maximum for the fluorophore; ‡ Extinction coefficient for the dye;

§ Correction factor =  $A_{260}$  for the free dye /  $A_{\text{max}}$  for the free dye.

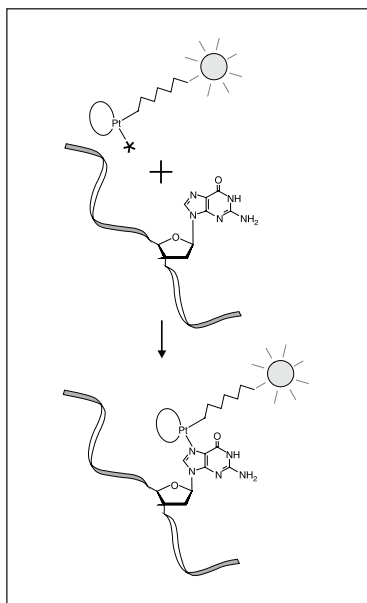


Figure 1. The ULS® reagent in the ULYSIS® Nucleic Acid Labeling Kits reacts with the N7 of guanine residues to provide a stable coordination complex between the nucleic acid and the fluorophore label.

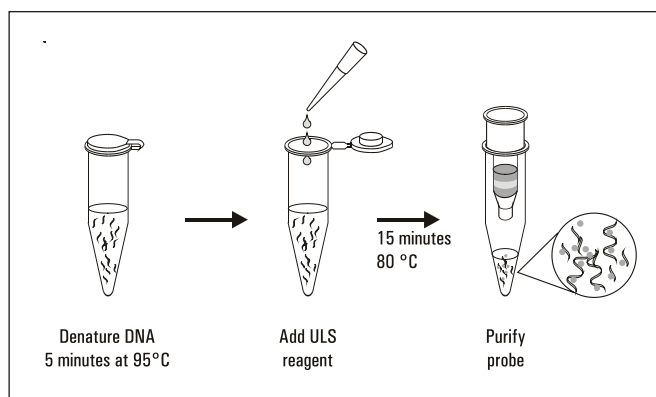


Figure 2. ULYSIS® nucleic acid labeling method.

## Before you begin

### Storage and handling

Upon receipt, store the kit at  $\leq -20^{\circ}\text{C}$  in a desiccator, protected from light. When stored properly, the kit should be stable for at least six months.

### Caution

No data are available addressing the mutagenicity or toxicity of the ULS® labeling reagents. Because these reagents bind to nucleic acids, they should be handled with appropriate care. ULS® labeling reagents should be disposed of safely and in accordance with applicable regulations. ULS® labeling reagents can be removed from aqueous solutions by filtration through activated charcoal. The charcoal and adsorbed dye must then be disposed of in a safe and appropriate manner.

**Materials required but not provided**

- Ethanol, absolute
- Sodium acetate, 3 M (pH 5.2)
- Spin column (see step 2.5)

**Spectral characteristics**

For the best results, match the light source, excitation filters, and emission filters to the spectral characteristics of the dye. Refer to Table 2 (page 2) for this information.

## Sample preparation

---

The ULS<sup>®</sup> labeling reagent will label double-stranded or single-stranded DNA of any length. However, DNA longer than about 1,000 bp may aggregate when labeled and precipitate out of solution, most likely due to hydrophobic interactions between closely spaced dye molecules. If such a probe is used for *in situ* hybridization, large aggregates may form over the sample, obscuring the signal. To produce optimally labeled probes, DNA longer than 1,000 bp should be fragmented by DNase I digestion prior to labeling as described below. Alternatively, it is possible to fragment the DNA by sonication or by digestion with a restriction enzyme that has a four-basepair recognition sequence. The fragmentation step is not needed for labeling DNA shorter than ~1,000 bp, or for labeling RNA samples (proceed to step 2.1).

**DNase I stock solution**

Centrifuge the vial of DNase I (Component D) briefly in a microcentrifuge to deposit the solids in the bottom of the tube. Make a 1 mg/mL DNase I stock solution by adding 100  $\mu$ L of chilled DNase I storage buffer (Component E) to the vial of DNase I (Component D). MIX GENTLY by inversion to dissolve the DNase I. Do not vortex, as the DNase I is unusually sensitive to physical denaturation. Store the DNase I stock solution at  $\leq -20^{\circ}\text{C}$  for up to six months.

**DNase I digestion protocol**

The following DNase I digestion protocol for 1  $\mu$ g of DNA results in fragments ~100 bp to ~1,000 bp in length. This protocol has been successfully applied to a wide variety of DNA species and preparations with excellent results when subsequently labeled and used in FISH experiments. However, because DNase I preparations show some variation, we recommend that its activity be titrated before using it with your samples.

- 1.1** Prepare 200 $\mu$ L of 1X DNase I reaction buffer by adding 20  $\mu$ L of 10X DNase I reaction buffer (Component F) to 180  $\mu$ L of nuclease-free H<sub>2</sub>O (Component H). Chill the solution on ice.
- 1.2** Immediately before use, make a DNase I working solution by diluting the DNase I stock solution into chilled 1X DNase I reaction buffer (from step 1.1). To find the optimal concentration of DNase I to use, test 1,650-, 2,500-, 5,000-, and 10,000-fold dilutions of DNase I using either the DNA from calf thymus (Component G) provided with the kit or a small portion of sample DNA. It is most convenient to perform the dilutions in two steps. First, dilute 1  $\mu$ L of DNase I stock solution into 49  $\mu$ L of 1X DNase I reaction buffer. Mix thoroughly by gently flicking the tube. Next, dilute 0.5, 1, 2, or 3  $\mu$ L of the first dilution into 1X DNase I reaction buffer to a final volume of 100  $\mu$ L. Mix thoroughly by gently flicking the tube. Leave DNase I working solutions on ice until ready to use. The working solutions are not stable and should be used only on the day prepared.

- 1.3 Add the following to a microfuge tube on ice in the order indicated, starting with the volume of nuclease-free H<sub>2</sub>O (Component H) necessary to achieve a final volume of 25  $\mu$ L. The following example is for 1  $\mu$ g of DNA, from a 0.1 mg/mL solution:
- 9.5  $\mu$ L nuclease-free H<sub>2</sub>O (Component H)
  - 2.5  $\mu$ L 10X DNase I reaction buffer (Component F)
  - 10  $\mu$ L 0.1 mg/mL sample DNA in TE buffer (10 mM Tris, 1 mM EDTA, pH 8.0)
  - 3  $\mu$ L DNase I working solution (from step 1.2)
- 1.4 Incubate at 37°C for 10 minutes. Stop the reaction by plunging the reaction tube into an ice bath.
- 1.5 To check the extent of DNase I digestion, take 2.5  $\mu$ L from the reaction mixture, heat it at 65°C for 10 minutes and then analyze it by agarose gel electrophoresis. Confirm that the DNA has been digested to fragments between ~100 bp and ~1,000 bp in length.

## Labeling reaction

---

The optimal amount of ULS<sup>®</sup> labeling reagent to use depends upon the application. Table 3 shows the optimal amount of ULS<sup>®</sup> labeling reagent required to label 1  $\mu$ g of DNA for hybridization to human metaphase chromosome spreads (FISH) or dot blots. To label more than 1  $\mu$ g of DNA, we recommend using proportionally more ULS<sup>®</sup> reagent and scaling up the reaction. Although we have not fully optimized the reaction for RNA labeling, preliminary experiments show good results using the same amounts of ULS<sup>®</sup> reagent as for DNA labeling.

### ULS<sup>®</sup> reagent stock solution

To prepare the ULS<sup>®</sup> labeling reagent stock solutions, except for Alexa Fluor<sup>®</sup> 488 ULS<sup>®</sup> reagent, add 100  $\mu$ L of 50% DMF (Component B) or 100  $\mu$ L of DMSO (Component B), depending upon the kit, to the vial containing the ULS<sup>®</sup> reagent (Component A). Vortex mix until all of the ULS<sup>®</sup> labeling reagent has dissolved and no particulate matter remains. In order to completely dissolve the ULS<sup>®</sup> reagents, vigorous vortexing, followed by pipetting up and down, may be required. DO NOT HEAT the solution. ULS<sup>®</sup> reagent stock solutions may be stored at 4°C for up to six months.

For the Alexa Fluor<sup>®</sup> 488 ULS<sup>®</sup> reagent (Cat. no. U21650), add 5  $\mu$ L of DMSO (Component B) to one vial of the ULS<sup>®</sup> reagent (Component A). Vortex mix until all of the ULS<sup>®</sup> labeling reagent has dissolved and no particulate matter remains. The stock solution will be nearly colorless. The Alexa Fluor<sup>®</sup> 488 ULS<sup>®</sup> reagent stock solution is stable for at least one month at 4°C.

**Note:** The ULS<sup>®</sup> Alexa Fluor<sup>®</sup> 488 labeling reagent has an absorption maximum of 257 nm and almost no absorption at 492 nm. Lack of fluorescence under 492 nm light does not mean there is no reagent present.

### Labeling protocol

- 2.1 Precipitate 1  $\mu$ g of DNA by adding 1/10 volume of 3 M sodium acetate (pH 5.2) and two volumes of absolute ethanol, freeze at -70°C for 30 minutes and then centrifuge for 15 minutes at 12,000 rpm. Wash the pellet with 70% ethanol and allow it to air dry. Resuspend the pellet in 20  $\mu$ L of the labeling buffer (Component C). DO NOT use ammonium acetate for DNA precipitation, as the residual ammonium ions will interfere with the ULS<sup>®</sup> labeling reaction. DO NOT use carrier DNA to assist in the ethanol precipitation, as this will also interfere with the ULS<sup>®</sup> labeling reaction.

- 2.2 Denature the DNA (from step 2.1) at 95°C for 5 minutes and then snap cool on ice. Centrifuge the tube briefly to redeposit the sample to the bottom of the tube.

**Note:** Denaturation is not required for labeling; however, it improves labeling efficiency by 20–40%.

- 2.3 Referring to Table 3, below, add the appropriate volume of ULS<sup>®</sup> labeling reagent stock solution to the tube containing the denatured sample DNA. If necessary, add labeling buffer (Component C) to bring the final volume to 25 µL.

**Table 3.** Amount of ULS<sup>®</sup> reagent solution required to label 1 µg of DNA for different applications.

ULS <sup>®</sup> Labeling Reagent	Application		Acceptable Labeling Ratios *
	FISH	Dot Blots	
Alexa Fluor <sup>®</sup> 488	1.0 µL	1.0 µL	43–67
Alexa Fluor <sup>®</sup> 546	5 µL	4 µL	44–89
Alexa Fluor <sup>®</sup> 594	5 µL	2.5 µL	40–98
Alexa Fluor <sup>®</sup> 647	5 µL	5 µL	30–40

\* Base:dye ratio for the labeled nucleic acid. See *Calculating the Labeling Efficiency and Concentration of Nucleic Acid* on page 5 for instructions on how to determine this ratio.

- 2.4 Incubate the reaction at 80°C for 15 minutes (for RNA, incubate at 90°C for 10 minutes). Stop the reaction by plunging the reaction tube into an ice bath. Centrifuge the tube briefly to redeposit the sample to the bottom of the tube.
- 2.5 The DNA must now be purified from the excess ULS<sup>®</sup> labeling reagent. We recommend purifying the labeled DNA by using a gel filtration–based spin column, according to the manufacturer’s protocol. For example, BioRad<sup>®</sup> Micro Bio-Spin<sup>®</sup> P-30 or Princeton Separations Centri-Sep<sup>™</sup> columns give good results. Avoid using silica-based separation techniques.

## Calculating the labeling efficiency and concentration of nucleic acid

The relative efficiency of a labeling reaction can be evaluated by calculating the approximate ratio of bases to dye molecules. This ratio can be determined by measuring the absorbance of the nucleic acid at 260 nm and the absorbance of the dye at its absorbance maximum ( $\lambda_{\max}$ ) and by using the Beer-Lambert law:

$$A = \epsilon \times \text{path length (cm)} \times \text{concentration (M)},$$

where  $\epsilon$  is the extinction coefficient in  $\text{cm}^{-1}\text{M}^{-1}$ . The absorbance measurements can also be used to determine the concentration of nucleic acid in the sample. Values needed for these calculations are found in Table 2 (page 2) and Table 4 (page 6). Acceptable labeling ratios for ULS-labeled nucleic acids are listed in Table 3 (page 5).

## Measuring the base:dye ratio

3.1 Measure the absorbance of the nucleic acid–dye conjugate at 260 nm ( $A_{260}$ ) and at the  $\lambda_{\text{max}}$  for the dye ( $A_{\text{dye}}$ ). Measure the background absorbance at 260 nm and  $\lambda_{\text{max}}$  using buffer alone, and subtract these numbers from the raw absorbance values for the sample. The  $\lambda_{\text{max}}$  values for the fluorophores used in the ULYSIS® kits are given in Table 2 (page 2).

- To perform these measurements, the nucleic acid–dye conjugate should be at a concentration of at least 5  $\mu\text{g}/\text{mL}$ . Depending on the dye used and the degree of labeling, a higher concentration may be required.
- For most applications, it will be necessary to measure the absorbance of the entire sample using either a conventional spectrophotometer with a 100 or 200  $\mu\text{L}$  cuvette or an absorbance microplate reader with a microplate.
- Use a cuvette or microplate that does not block UV light and that is clean and nuclease-free. Note that most plastic disposable cuvettes and microplates have significant absorption in the UV.

3.2 Correct for the contribution of the dye to the  $A_{260}$  reading. Most fluorescent dyes absorb light at 260 nm as well as at their  $\lambda_{\text{max}}$ . To obtain an accurate absorbance measurement for the nucleic acid, it is therefore necessary to account for the dye absorbance using a correction factor ( $\text{CF}_{260}$ ). Use the  $\text{CF}_{260}$  values given in Table 2 (page 2) in the following equation:

$$A_{\text{base}} = A_{260} - (A_{\text{dye}} \times \text{CF}_{260})$$

3.3 Calculate the ratio of bases to dye molecules using the following equation:

$$\text{base dye} = (A_{\text{base}} \times \epsilon_{\text{dye}}) / (A_{\text{dye}} \times \epsilon_{\text{base}})$$

where  $\epsilon_{\text{dye}}$  is the extinction coefficient for the fluorescent dye (found in Table 2, page 2) and  $\epsilon_{\text{base}}$  is the average extinction coefficient for a base in double stranded DNA (dsDNA), long single-stranded DNA (ssDNA), or RNA (found in Table 4, below). Note that since the calculation is a ratio, the path length has canceled out of the equation.

**Table 4.** Average values for bases in different nucleic acids.

Nucleic Acid	$\epsilon_{\text{base}} \text{ (cm)}^{-1} \text{ (M)}^{-1} *$	$\text{MW}_{\text{base}} \dagger$
dsDNA	6,600	330
ssDNA	8,919	330
RNA	8,250	340

\* Average extinction coefficient for a base; † Average molecular weight for a base (g/mol).

## Measuring the concentration of nucleic acid

The absorbance values,  $A_{260}$  and  $A_{\text{dye}}$ , and the Beer-Lambert law may also be used to measure the concentration of nucleic acid in the sample ([N.A.]). In order to obtain an accurate measurement for a dye-labeled nucleic acid, a dye-corrected absorbance value ( $A_{\text{base}}$ ) must be used, as explained in step 3.2. In addition, for concentration measurements, the path length (in cm) is required. If the path length of the cuvette or of the solution in a microplate well is unknown, consult the manufacturer. Follow steps 3.1 and 3.2 above and then use the following equation:

$$[\text{N.A.}] \text{ (mg/mL)} = (A_{\text{base}} \times \text{MW}_{\text{base}}) / (\epsilon_{\text{base}} \times \text{path length})$$

## Product list

Current prices may be obtained from our website or from our Customer Service Department.

Cat. no.	Product Name	Unit Size
U21650	ULYSIS® Alexa Fluor® 488 Nucleic Acid Labeling Kit *20 labelings* .....	1 kit
U21652	ULYSIS® Alexa Fluor® 546 Nucleic Acid Labeling Kit *20 labelings* .....	1 kit
U21654	ULYSIS® Alexa Fluor® 594 Nucleic Acid Labeling Kit *20 labelings* .....	1 kit
U21660	ULYSIS® Alexa Fluor® 647 Nucleic Acid Labeling Kit *20 labelings* .....	1 kit

## Purchaser notification

These high-quality reagents and materials must be used by, or directly under the supervision of, a technically qualified individual experienced in handling potentially hazardous chemicals. Read the Safety Data Sheet provided for each product; other regulatory considerations may apply.

### Obtaining Support

For the latest services and support information for all locations, go to [www.lifetechnologies.com](http://www.lifetechnologies.com).

At the website, you can:

- Access worldwide telephone and fax numbers to contact Technical Support and Sales facilities
- Search through frequently asked questions (FAQs)
- Submit a question directly to Technical Support ([techsupport@lifetech.com](mailto:techsupport@lifetech.com))
- Search for user documents, SDSs, vector maps and sequences, application notes, formulations, handbooks, certificates of analysis, citations, and other product support documents
- Obtain information about customer training
- Download software updates and patches

### SDS

Safety Data Sheets (SDSs) are available at [www.lifetechnologies.com/sds](http://www.lifetechnologies.com/sds).

### Certificate of Analysis

The Certificate of Analysis provides detailed quality control and product qualification information for each product. Certificates of Analysis are available on our website. Go to [www.lifetechnologies.com/support](http://www.lifetechnologies.com/support) and search for the Certificate of Analysis by product lot number, which is printed on the product packaging (tube, pouch, or box).

### Limited Product Warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at [www.lifetechnologies.com/termsandconditions](http://www.lifetechnologies.com/termsandconditions). If you have any questions, please contact Life Technologies at [www.lifetechnologies.com/support](http://www.lifetechnologies.com/support).

### Disclaimer

LIFE TECHNOLOGIES CORPORATION AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

### Important Licensing Information

These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

All trademarks are the property of Thermo Fisher Scientific and its subsidiaries, unless otherwise specified. ULS is a trademark of Kreotech Biotechnology B.V. Cy is a trademark of GE Healthcare UK. Lissamine is a trademark of PerkinElmer, Inc. SpectrumAqua is a trademark of Abbott Molecular, Inc.

©2015 Thermo Fisher Scientific Inc. All rights reserved.