

**DELFLIA[®] Tb-N1 ITC Chelate &
Terbium Standard****AD0009****For Research Use Only****INTRODUCTION**

DELFLIA[®] Tb-N1 ITC Chelate is optimized for the terbium labelling of proteins and peptides for use in dissociation-enhanced time-resolved fluorometric assays (1). The reagent is the Tb³⁺-chelate of N¹-(p-isothiocyanatobenzyl)diethylenetriamine-N¹, N², N³, N³-tetraacetic acid. The DTTA group (diethylenetriaminetetraacetic acid) forms a stable complex with Tb³⁺ and the isothiocyanato group reacts with free amino groups on the proteins and peptides, forming a stable, covalent thiourea bond.

PACKAGE CONTENTS

1 vial (1 mg, 1.5 µmol) of Tb-N1 ITC Chelate
1 vial (0.5 mL) of 100 nmol/L Terbium Standard

STORAGE

The manufacturing date of the chelate is stated on the vial label. Store the chelate with a desiccant at -20°C before reconstitution. Store the standard at +2 - +8°C.

REAGENT RECONSTITUTION

Dissolve the chelate in distilled water (e.g. in 100 µL giving 15 mmol/L solution of the chelate) for immediate use. Keep at 0°C (icebath).

If all chelate is not used during the same day, it can be dissolved either in 10 mmol/L sodium succinate (pH 5.0) or in 10 mmol/L sodium acetate (pH 4.8) (pH of the reconstituted chelate should be below 7 for storage purposes). Keep at 0°C (icebath).

RECONSTITUTED STABILITY

The chelate reconstituted in water should be used the same day.

For a longterm storage at -20°C, the chelate should be dissolved in succinate or acetate buffer and aliquoted.

WARNINGS AND PRECAUTIONS

This labelling reagent is intended for research use only.

The handling of concentrated Tb³⁺-solutions constitutes a contamination risk, which may cause elevated backgrounds in an assay based on time-resolved fluorometry. Keep the labelling reagents and required accessories separated from the place and accessories where the actual assay is performed.

Disposal of all waste should be in accordance with local regulations.

LABELLING OF PROTEINS

Tb-N1 ITC chelate has an aromatic isothiocyanato group as a reactive arm. Isothiocyanato group reacts with primary aliphatic amino groups at alkaline pH.

1. Parameters of labelling reaction

Parameters of labelling reaction include protein concentration, pH, temperature, reaction time and molar excess of chelate over protein.

2. Labelling

The recommended reaction conditions for labelling of proteins are pH 9 - 9.3, +4°C and overnight incubation. Under these conditions, the following calculations are valid for labelling of a protein with an isoelectric point (pI) between 4 and 7.

Protein concentration (mg/mL)	Percentage of chelate reacted
5	20 %
2.5	10 %
1	4 %

Table 1. The effect of protein concentration on the percentage of Tb-N1 ITC chelate reacting with the protein.

For example, if a protein (pI around 6, molecular weight 160 000) is reacted at a concentration of 5 mg/mL under the conditions described above, a 40-fold molar excess of chelate over protein would give a labelling degree of about 8 Tb-N1 ITC chelates per protein.

If the protein to be labelled is not stable in the labelling conditions (+4°C, pH 9 - 9.3, overnight incubation), it is possible to run a 4-hour reaction (+4°C, pH 9 - 9.3) by increasing the molar excess of chelate over protein. A suitable amount of chelate is three times higher for the 4-hour reaction than for the overnight reaction. For example, if a protein (5 mg/mL in the labelling reaction) requires 40-fold molar excess of chelate during overnight reaction for the introduction of 8 chelates per protein, 120-fold molar excess of Tb-N1 ITC chelate is needed to obtain the same label incorporation during a 4-hour reaction. Note that the separation of unreacted chelate from labelled protein is more difficult with four hour labelling because of the greater amount of chelate in the reaction.

If the protein to be labelled is not stable at pH 9 – 9.3 labelling can also be done at pH 8.4 – 8.5. Reactivity of Tb-N1 chelate at pH 8.4 – 8.5 is only about 25 % of the reactivity at pH 9 – 9.3. For example, if a protein (pI around 6, molecular weight

160 000) is reacted at a concentration of 5 mg/mL, pH 8.4 – 8.5, +4°C and overnight reaction, a 160-fold molar excess of chelate over protein would give a labelling degree of about 7 – 9 Tb-N1 ITC chelates per protein.

Suitable number of Tb-N1 ITC chelates coupled to a protein depends on the molecular weight (MW). When the MW of a protein is higher than 100 000, 4 - 15 chelates per protein is a good labelling yield. For proteins with a MW in the range of 30 000 - 100 000 the preferred number of coupled chelates is 2 - 10. Proteins with a MW less than 30 000 should be labelled with 1 - 3 chelates. The given values may be higher for basic proteins (pI between 8 and 10).

LABELLING OF PEPTIDES

Peptides (size up to about 40 amino acids) are labelled like proteins except that the molar excess of chelate over peptide is lower than in protein labelling. Recommended molar excess of chelate over peptide is 3 - 6 (peptide concentration 5 - 20 mg/mL), 5 - 10 (peptide concentration 2.5 - 5 mg/mL) or 8 - 30 (peptide concentration 1 - 2.5 mg/mL). Labelling is usually performed at +4°C but, if a peptide to be labelled is very stable, it can be labelled at room temperature (+20 - +25°C). Suitable number of chelates coupled to a peptide is 1 - 2 depending on the peptide.

LABELLING PROCEDURE

The protein or peptide to be labelled must not be stabilized with a protein (e.g. bovine serum albumin (BSA), casein or gelatin).

1. Pretreatment

If the buffer including the protein or peptide to be labelled contains primary amines (e.g. Tris, ammonium ions), sulfhydryl groups (e.g. mercaptoethanol) or sodium azide, a pretreatment step is necessary. The above mentioned compounds interfere with labelling. Suitable methods for removing interfering compounds include gel filtration (e.g. NAP¹ and PD-10 columns), dialysis and reverse phase HPLC (RP-HPLC).

2. Concentrating protein and peptide

If a protein is too dilute (less than 1 mg/mL) or it is preferable to use less chelate to facilitate purification after labelling, a concentration step is necessary. Suitable concentrators are e.g. Centricon and Centriprep concentrators².

If the concentration of a peptide is too low for an efficient labelling reaction (less than 1 mg/mL), a vacuum centrifuge can be used to concentrate a peptide solution.

3. Reconstitution of chelate

After calculating the amount of chelate needed in the reaction, Tb-N1 ITC chelate is dissolved either in water (for immediate use only) or in 10 mmol/L sodium succinate pH 5.0 or in 10 mmol/L sodium acetate pH 4.8 (in case some of the chelate will be

¹ NAP is a trademark of Amersham Pharmacia Biotech.

² Centricon and Centriprep are registered trademarks of Millipore Corp.

stored for future purposes). Suitable concentration for reconstituted Tb-N1 ITC is 5 - 30 mmol/L (1mg of Tb-N1 ITC is 1.5 μ mol). For example, dissolving 1 mg of Tb-N1 ITC chelate in 200 μ L gives a concentration of 7.5 mmol/L. After dissolving the chelate is kept on ice for immediate use.

4. Labelling

If the protein or peptide is already in a labelling buffer (50 - 100 mmol/L sodium carbonate, pH 9 - 9.3) after the pretreatment or reconstitution, the calculated amount of chelate is added into the protein (peptide) solution on ice.

If the protein (peptide) is not in a labelling buffer, 1 mol/L sodium carbonate (pH 9 - 9.3) is added to adjust buffer concentration to 50 - 100 mmol/L followed by the calculated amount of reconstituted chelate.

In both cases, pH is checked after adding the chelate by applying a 0.5 μ L sample on a pH-paper or pH-stick. It is advisable to check the performance of the pH-paper or pH-stick with 50 - 100 mmol/L sodium carbonate buffer of known pH. A suitable pH-paper is Spezial Indikatorpapier pH 8.2 - 10.0 (Merck Art. No. 9558). If necessary, pH of the reaction mixture is adjusted to 9 - 9.3 using either 0.5 - 1 mol/L HCl or 0.5 - 1 mol/L NaOH.

After adding all necessary components and checking pH, incubate at +4°C overnight (or for 4 hours).

5. Purification

Separation of the labelled protein from unreacted chelate is performed by gel filtration. Elution buffer should be Tris-HCl based, e.g. 50 mmol/L Tris-HCl (pH 7.8) containing 0.9 % NaCl and 0.05 % sodium azide (TSA buffer). Proteins with a molecular weight over 100 000 can be purified using Superdex 200 column or a combination of Sephadex G-50 (e.g. 1 x 10 cm) layered on Sepharose 6B (e.g. 1 x 30 - 40 cm)³. Proteins with a MW in the range of 30 000 - 100 000 are best purified using Superdex 75 or Sephadex G-50. Sephadex G-50 is suitable also for purification of proteins with a MW between 15 000 and 30 000.

In gel filtration the collected fraction size should be about 1 mL. The gel filtration eluate can be monitored by UV-absorbance at 280 nm. The first peak contains the labelled protein and the second peak unreacted chelate. The Tb concentration of the fractions can be measured by making a 1:1000 - 1:10000 dilution in DELFIA Enhancement Solution (prod. no. 1244-105) and in DELFIA Enhancer (prod. no. C500-100). DELFIA Enhancer must equal 1/5 of the total volume (e.g. 0.2 mL if the total volume is 1 mL). The dilutions should be mixed gently and let stand for about 2 minutes before measuring in a time-resolved fluorometer. The fractions from the first peak with the highest Tb counts are pooled and characterized. Collecting of fractions should be stopped at least two fractions before the signal of unreacted chelate starts to rise.

When labelling only a small amount of antibody (< 0.5 mg) the purification can be done with a PD-10 column by applying the reaction mixture in the equilibrated column and

³ Superdex, Sephadex and Sepharose are trademarks of Amersham Pharmacia Biotech.

collecting 0.5 mL fractions. The fractions from the first peak with the highest Tb counts should be pooled and characterized.

Peptides having at least about 25 amino acid residues can be purified from the unreacted chelate on Sephadex G-25 or alternatively using reverse phase HPLC. Small peptides (less than 25 amino acid residues) can be purified from the unreacted chelate and at least in some cases also from the unlabelled peptide by using reverse phase HPLC. The labelled peptide is eluted from the column in acetonitrile gradient in 0.02 - 0.1 mol/L triethylammonium acetate (pH 7.5). After collecting the labelled peptide acetonitrile is evaporated. It is advisable to add 50 µL/mL 1 mol/L Tris-HCl (pH 8.5) before evaporation of acetonitrile to make sure that pH stays neutral.

Proteins MW above 100 000	Proteins MW 30-100 000	Proteins MW 15-30 000	Proteins and peptides MW 2500-15 000	Peptides MW below 2500 (< 25 aa)
Superdex 200	Superdex 75	Sephadex G-50	Sephadex G-25	RP-HPLC
Sephadex G-50 /Sephacrose 6B	Sephadex G-50		RP-HPLC	

Table 2. Recommended columns for purification of proteins and peptides after labelling with Tb-N1 ITC chelate.

There should be dedicated columns for each lanthanide (europium, terbium, samarium, dysprosium) used for labelling. After purification, the gel filtration column should be decontaminated by washing with 10 mmol/L phthalate buffer (pH 4.1) containing 0.01 % DTPA. Before each run, it is advisable to saturate and further purify the gel filtration column on the previous day by applying concentrated BSA solution of high purity (e.g. 0.5 mL 7.5 % BSA; for suitable BSA see section STORAGE OF LABELLED COMPOUNDS). After adding the BSA the column should be equilibrated overnight. RP-HPLC columns can be washed using the phthalate buffer described above.

CHARACTERIZATION OF LABELLED PROTEINS AND PEPTIDES

To determine Tb³⁺ concentration in the labelled proteins (peptides), the labelled protein (peptide) is diluted in DELFIA Enhancement Solution (prod. no. 1244-105) and mixed gently. Wait for 5 minutes to allow Tb to dissociate and then add 1/5 of the total volume (e.g. 0.2 mL into 0.8 mL to give a total volume 1 mL; final dilution 1:10 000) DELFIA Enhancer (prod. no. C500-100). After mixing and incubating for 2 minutes measure in a time-resolved fluorometer against 100 nmol/L Tb standard diluted 1:100 in the same way as the sample (1 nmol/L Tb in Enhancement Solution plus DELFIA Enhancer in a clear 96-well plate, 250 µL per well, gives about 500 000 cps in 1420 VICTOR™ Multilabel Counter).

The protein (peptide) concentration can be measured with a suitable method (e.g. Lowry) or calculated from absorbance at 280 nm. The molar absorptivity of reacted Tb-N1 chelate is 8000 at 280 nm (1 µmol/L reacted chelate gives an absorbance of 0.008 at 280 nm).

FILTRATION

To remove particles and possible aggregates the labelled protein should be filtered through a 0.22 µm low protein binding membrane.

STORAGE OF LABELLED COMPOUNDS

To ensure stability, the labelled proteins and peptides should be stored at a high concentration and in the absence of chelators or competing metals in the buffer. A concentrated solution (0.1 mg/mL or higher) can be stored without any stabilizer. With lower concentrations, the stability can be increased by adding purified BSA (a component in the DELFIA Eu-labelling kit, prod. no. 1244-302, and Sm-labelling kit, prod. no. 1244-303; available also as a specialty product from PerkinElmer Life Sciences, prod. no. CR84-100 Stabilizer) to a final concentration of 0.1 %. Temperature during storage is determined by the stability of the protein. Suitable temperatures are e.g. +4°C, -20°C and -70°C.

USE OF LABELLED PROTEINS

The amount of proteins, incubation time, temperature and the buffers used must be optimized for each particular analyte. As a general rule about 5 - 100 ng of the labelled proteins per tube or well can be used. The DELFIA Assay Buffer (product no. 1244-106) is optimal for most assays. It contains NaCl, Tris-HCl, bovine serum albumin (BSA), bovine gamma globulins, Tween 40, diethylenetriaminepentaacetic acid (DTPA), < 0.1 % NaN₃ and an inert red dye. If this assay buffer cannot be used, it is recommended to use a Tris-HCl buffer containing 20 µmol/L EDTA or DTPA to keep the fluorescence background low.

The labelled protein as such is practically non-fluorescent. After binding assay DELFIA Enhancement Solution dissociates Tb ions from labelled protein into solution. For detection of Tb-labelled proteins first add 200 µL of Enhancement Solution per well and shake **slowly** for 5 minutes. Then add 50 µL of DELFIA Enhancer to each well and shake **slowly** for 5 minutes before measuring with the time-resolved fluorometer (1420 VICTOR or 1234 DELFIA Research Fluorometer).

WARRANTY

Purchase of the product gives the purchaser the right to use this material in his own research, development, and investigational work. The product is not to be injected into humans or used for diagnostic procedures. Wallac Oy reserves the right to discontinue or refuse orders to any customer who plans to use these products for any other purposes.

Wallac Oy does not warrant or guarantee that the product is merchantable or satisfactory for any particular purpose, nor free from any claim of foreign or domestic patent infringement by a third party, and there are no warranties, expressed or implied, to such effect. Wallac Oy will not be liable for any incidental, consequential or contingent damages involving their use including damages to the property or personal injuries.

All information supplied with the product and technical assistance given is believed to be accurate, but it remains the responsibility of the investigator to confirm all technical aspects of the application. We appreciate receiving any additions, corrections, or updates to information supplied to the customer.

REFERENCES

1. Hemmilä, I., Mukkala, V-M., Latva, M. and Kiilholma, P. (1993): Di- and tetracarboxylate derivates of pyridines, bipyridines and terpyridines as luminogenic reagents for time-resolved fluorometric determination of terbium and dysprosium. J. Biochem. Biophys. Methods, **26**, 283-290.

PATENTS

The reagents are covered by the following patents on both the chemical structure and the dissociation enhancement principle:

- Patent 1. Mikola, H., Mukkala, V-M. and Hemmilä, I. (1987): Eur. Patent No. 139,675.
Mikola, H., Mukkala, V-M. and Hemmilä, I. (1989): US Patent No. 4,808,541.
- Patent 2. Hemmilä, I. and Dakubu S. (1982): Eur. Patent No. 64,484.
Hemmilä, I. and Dakubu S. (1982): US Patent No. 4,565,790.