

## 3-Nitrotyrosine Competitive ELISA Kit

Catalog Number EEL008 (96 tests)

Rev 2.1

For safety and biohazard guidelines, read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

### Product description

The 3-Nitrotyrosine (3-NT) Immunoassay Kit is a solid-phase Competitive Enzyme-Linked Immunosorbent Assay (ELISA). This assay is designed to detect and quantify the level of 3-NT in serum, plasma, and other biological fluids independent of species.

3-Nitrotyrosine is a product of tyrosine nitration mediated by reactive nitrogen species (ex. peroxynitrite anion and nitrogen dioxide). Nitrotyrosine is identified as an indicator of cell damage, inflammation, and NO (nitric oxide) production. In many disease states, oxidative stress increases the production of superoxide ( $O_2^-$ ) and free radical oxidant peroxynitrite ( $ONOO^-$ ).  $ONOO^-$  is capable of oxidizing several lipoproteins and nitrating tyrosine residues in various proteins.

### Contents and storage

An unopened kit can be stored at 2-8°C for 12 months. After opening, store the items separately according to the conditions specified in the table below.

**Note:** All reagent bottle caps must be tightened to prevent evaporation and microbial pollution. The volume of reagents in partial shipments is slightly more than the volume marked on the label, please use accurate measuring equipment instead of directly pouring into the vial(s).

Components	Quantity (96 tests)	Storage
3-NT Antigen Coated Microplate	8 wells x 12 strips	-20°C, 12 months
3-NT Standard	2 vials	
3-NT Biotinylated Detection Antibody (100X)	120 µL	
HRP Conjugate (100X)	120 µL	-20°C (Protect from light), 12 months
Standard & Sample Diluent	2×20 mL	2-8°C, 12 months
Biotinylated Detection Antibody Diluent	14 mL	
HRP Conjugate Diluent	14 mL	
Wash Buffer Concentrate (25X)	30 mL	
Substrate Reagent	10 mL	2-8°C (Protect from light), 12 months
Stop Solution; contains 1 M H <sub>2</sub> SO <sub>4</sub> , <b>CAUSTIC</b>	10 mL	2-8°C, 12 months
Plate Sealer	5	

## Required materials

- Distilled or deionized water
- Microtiter plate reader with software capable of measurement at or near 450 nm
- Plate washer—automated or manual (squirt bottle, manifold dispenser, or equivalent)
- Calibrated adjustable precision pipettes and glass or plastic tubes for diluting solutions
- Incubator capable of maintaining 37°C.

## Procedural guidelines

---

**IMPORTANT!** Reagents are lot-specific. Do not mix or interchange different reagent lots from various kit lots.

---

Allow reagents to reach room temperature before use. Mix to redissolve any precipitated salts.

### Sample preparation guidelines

**Serum:** Allow samples to clot for 1 hour at room temperature or overnight at 2-8°C before centrifugation for 20 min at 1000×g at 2-8°C. Collect the supernatant to carry out the assay.

**Plasma:** Collect plasma using EDTA or heparin (EDTA-Na<sub>2</sub> is most recommended) as an anticoagulant. Centrifuge samples for 15 min at 1000×g at 2-8°C within 30 min of collection. Collect the supernatant to carry out the assay.

**Other biological fluids:** Centrifuge samples for 20 min at 1000×g at 2-8°C. Collect the supernatant to carry out the assay.

Note:

- Collect samples in pyrogen/endotoxin-free tubes.
- Samples should be assayed within 7 days when stored at 2-8°C, otherwise samples must be aliquoted and stored at -20°C (≤ 1 month) or -80°C (≤ 3 months). Avoid multiple freeze-thaw cycles of frozen samples. Thaw completely and mix well (do not vortex) prior to analysis.
- Avoid the use of hemolyzed or lipemic sera.
- If large amounts of particulate matter are present in the sample, centrifuge, or filter sample prior to analysis.

### Prepare samples

Sample concentrations should be within the range of the standard curve. Because conditions may vary, the optimal dilution for each application should be determined (It is recommended to carry out the preliminary test referring to the expected values of samples on page 8).

Use all prepared samples within 2 hours of dilution. It is not recommended to conduct experiments after 2 hours.

## **Prepare 1X Wash Buffer**

1. Dilute 30 mL of Wash Solution Concentrate (25X) with 720 mL of deionized or distilled water. Label as 1X Wash Buffer.
2. Store the concentrate and 1X Wash Buffer at 2-8°C. Use the diluted buffer within 3 months.

Note: if crystals have formed in the concentrate, warm it in a 40°C-water bath and mix it gently until the crystals have completely dissolved.

## **Prepare 1X Biotinylated Detection Antibody Solution**

**Note: The working solution should be prepared just before use**

1. Calculate the required amount before the experiment (50 µL/well). In preparation, slightly more than calculated should be prepared.
2. Centrifuge the Concentrated Biotinylated Detection Ab at 800×g at 2-8°C for 1 min.
3. Dilute the Concentrated Biotinylated Detection Ab (100X) to 1X working solution with Biotinylated Detection Ab Diluent.

## **Prepare 1X HRP Conjugate Solution**

**Note: The working solution should be prepared just before use**

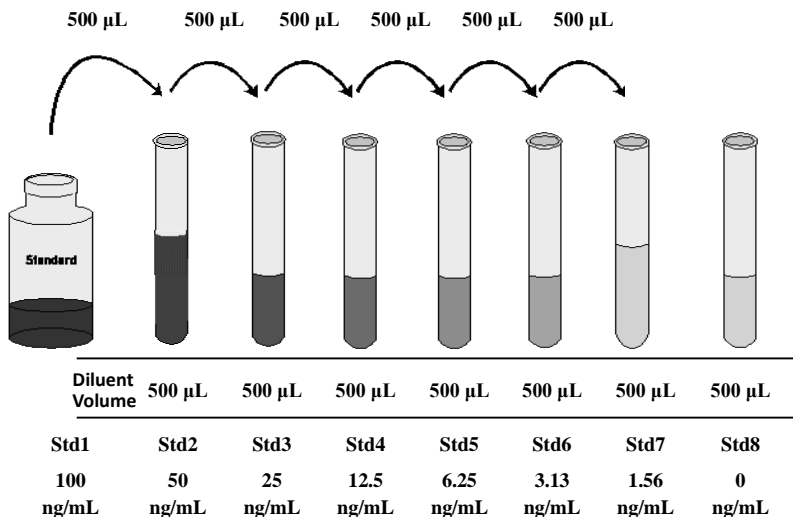
1. Calculate the required amount before the experiment (100 µL/well). In preparation, slightly more than calculated should be prepared.
2. Centrifuge the Concentrated HRP Conjugate at 800×g at 2-8°C for 1 min.
3. Dilute the Concentrated HRP Conjugate (100X) to 1X working solution with HRP Conjugate Diluent.

## **Prepare diluted standards**

Note: Use glass or plastic tubes for diluting standards.

1. Centrifuge the standard at 10,000×g at 2-8°C for 1 min to ensure the contents are at the bottom of vial.
2. Add 1 mL of Standard & Sample Diluent, let it stand for 10 min and invert it gently several times. Once fully dissolved, mix it thoroughly with a pipette. This reconstitution produces a working solution of 100 ng/mL.
3. Take 7 tubes, add 500 µL of Standard & Sample Diluent to each tube. Pipette 500 µL of the 100 ng/mL working solution to the first tube and mix up to produce a 50 ng/mL working solution. Pipette 500 µL of the solution from the former tube into the latter one according to this step. The last tube is regarded as a blank, don't pipette solution into it from the former tube. The recommended dilution gradient is as follows: 100、50、25、12.5、6.25、3.13、1.56、0 ng/mL.

4. The illustration of diluted standards below is for reference. Mix thoroughly between dilution gradients.



5. Use the diluted standards within 2 hours of preparation. If multiple standard tests are to be carried out, the redissolved standard solution (standard with the highest concentration) can be divided into 2-3 vials and frozen to  $-20^{\circ}\text{C}$ , to be used within half a month. Avoid multiple freeze-thaw cycles.

## Perform ELISA (Total assay time: 2 hours)

**IMPORTANT!** Perform a standard curve with each assay.

Allow all components to reach room temperature before use. Mix all liquid reagents prior to use.

Determine the number of 8-well strips required for the assay. Insert the strips in the frames for use. Re-bag any unused strips and frames, and store desiccated at  $-20^{\circ}\text{C}$  for future use. The silica pack in the bag keeps the plate dry.

### 1. Add antigen and biotinylated detection antibody

Note: solutions should be added to the bottom of the ELISA plate well, avoid touching the inside wall and causing foaming as much as possible.



- For the standard curve, add 50  $\mu\text{L}$  of standards to the appropriate wells. For samples, add 50  $\mu\text{L}$  of pretreated samples to the wells.
- Immediately add 50  $\mu\text{L}$  of **Biotinylated Detection Ab working solution** to each well
- Cover the plate with plate sealer and incubate for 45 min at 37  $^{\circ}\text{C}$ .
- Thoroughly aspirate the solution and wash wells 3 times with 350  $\mu\text{L}$  of 1X Wash Buffer. Decant the solution from each well, add 350  $\mu\text{L}$  of **wash buffer** to each well. Soak for 1 min and aspirate or decant the solution from each well and pat it dry against clean absorbent paper. Proceed immediately to the next step, making sure the wells do not dry out.



### 2. Add HRP conjugate

- Add 100  $\mu\text{L}$  **HRP Conjugate Working Solution** into each well.
- Cover the plate with plate sealer and incubate for 30 min at 37  $^{\circ}\text{C}$ .
- Thoroughly aspirate the solution and repeat the wash process for 5 times as conducted in step 1



### 3. Add substrate

- Add 90  $\mu\text{L}$  **Substrate Reagent** to each well.
- Cover the plate with plate sealer and incubate for about 15 min at 37  $^{\circ}\text{C}$ . Protect the plate from light.

**Note:** The reaction time can be shortened or extended according to the actual color change, but not more than 30 min.



### 4. Add stop solution

- Add 50  $\mu\text{L}$  **Stop Solution** to each well. This step should be done in the same order as the substrate solution. Tap the side of the plate gently to mix.
- The solution in the wells will change from blue to yellow.

● Ag



Biotinylated  
detection Ab



Streptavidin-  
HRP

## Read the plate and generate the standard curve

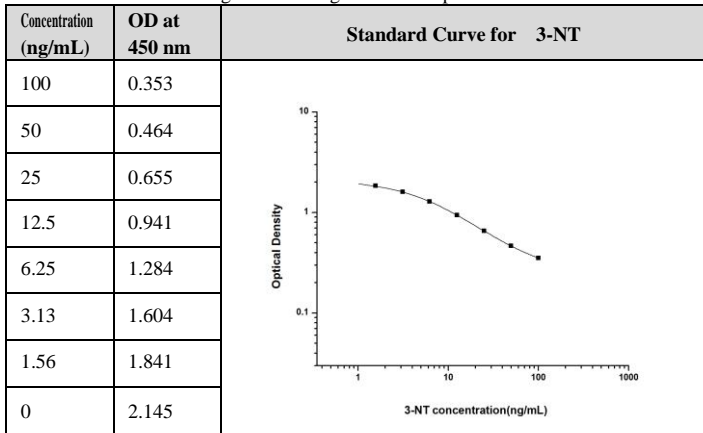
1. Preheat the Microplate Reader for about 15 min before OD measurement.
2. Read the absorbance at 450 nm. Read the plate within 10 minutes after adding the Stop Solution.
3. Use curve-fitting software to generate the standard curve. A four-parameter algorithm provides the best standard curve fit.
4. Read the concentrations for unknown samples and controls from the standard curve. Multiply value(s) obtained for sample(s) by the appropriate factor to correct for the sample dilution.

Note: If the OD of the sample under the lowest limit of the standard curve, you should re-test it with an appropriate dilution.

## Performance characteristics

### ■ Standard curve (example)

As the OD values of the standard curve may vary according to the conditions of the actual assay performance (e.g. operator, pipetting technique, washing technique or temperature effects), the operator should establish a standard curve for each test. Typical standard curve and data over the range of 0–100 ng/mL 3-NT is provided below for reference only.



### ■ Inter-assay Precision

Three Human serum samples with low, medium, and high level 3-NT were tested 10 ten times in duplicate to determine precision between assays.

Parameters	Sample 1	Sample 2	Sample 3
Mean (ng/mL)	5.1	13.1	36.6
Standard deviation	0.3	0.6	1.9
%CV	5.88	4.58	5.19

CV = Coefficient of Variation

### ■ Intra-assay Precision

Three Human serum samples with low, medium, and high level 3-NT were assayed in replicates of 20 to determine precision within an assay.

Parameters	Sample 1	Sample 2	Sample 3
Mean (ng/mL)	5	12.5	38.3
Standard deviation	0.3	0.7	1.8
%CV	6	5.6	4.7

CV = Coefficient of Variation

### ■ Expected values

Sixteen random Human serum/plasma samples were tested in the assay.

Sample Type	3-NT Range (ng/mL)	3-NT Average (ng/mL)
Serum (n=16)	4.1-13.2	5.9
Plasma (n=16)	3.7-9.8	4.3



### ■ Recovery

The recovery of 3-NT spiked at three different levels in Human samples throughout the range of the assay was evaluated in various matrices.

Sample Type	Range (%)	Average Recovery (%)
Serum (n=8)	97-108	102
EDTA plasma (n=8)	91-108	99
Cell culture media (n=8)	93-105	100

### ■ Linearity of dilution

Human Samples were spiked with high concentrations of 3-NT and diluted with Standard & Sample Diluent to produce samples with values within the range of the assay.

		Serum (n=5)	EDTA plasma (n=5)	Cell culture media(n=5)
1:2	Range (%)	87-99	92-105	97-112
	Average (%)	93	100	103
1:4	Range (%)	84-96	91-101	98-113
	Average (%)	90	96	105
1:8	Range (%)	88-99	93-107	101-116
	Average (%)	93	99	106
1:16	Range (%)	88-100	94-105	97-111
	Average (%)	94	100	103

## ▪ Specificity

This assay has been shown to detect 3-NT only from human, mouse, rat, monkey, rabbit and porcine samples. Please contact technical support for cross-reactivity information on other species.

## ▪ Sensitivity

The analytical sensitivity of the assay is 0.94 ng/mL 3-NT. This was determined by adding two standard deviations to the mean O.D. obtained when the zero standard was assayed 20 times and calculating the corresponding concentration.

## 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.thermofisher.com/us/en/home/global/terms-and-conditions.html](http://www.thermofisher.com/us/en/home/global/terms-and-conditions.html). If you have any questions, please contact Life Technologies at [www.thermofisher.com/support](http://www.thermofisher.com/support).

The information in this guide is subject to change without notice.

## DISCLAIMER

TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

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.

Corporate entity: Bender MedSystems GmbH | Campus Vienna Biocenter 2 | 1030 Vienna, Austria

© 2025 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.