

### Nitrotyrosine ELISA Kit ab210603

9 References [画像数 1](#)

#### 製品の概要

製品名 Nitrotyrosine ELISA Kit

検出方法 Colorimetric

再現性 Intra-Assay (同時再現性)

サンプル	N	平均値	SD	CV%
3 Samples	30			< 10%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
3 Samples	30			< 15%

サンプルの種類 Urine, Serum, Plasma, Other biological fluids, Cell Lysate

アッセイタイプ Competitive

検出感度 = 50 nM

検出範囲 62.5 nM - 8000 nM

全工程の試験時間 2h 00m

ステップ Multiple steps standard assay

製品の概要 Abcam's Nitrotyrosine *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit (ab210603) is designed for the competitive measurement of Free Nitrotyrosine protein in plasma, serum, cell lysates, urine, and other sample matrices. The ELISA utilizes an Nitrotyrosine-coated plate and an HRP-conjugated antibody for detection which allows for an assay range of 62.5 to 8000 nM Free Nitrotyrosine, with a sensitivity of 50 nM. The other highlights of this kit are a quick incubation time of 60 minutes, stable reagents, and an easy to use protocol.

特記事項 Nitrotyrosine has been identified as a marker of inflammation and NO production. Nitrotyrosine is formed in presence of the active metabolite NO. Various pathways including the formation of peroxynitrite lead to nitrotyrosine production. Since nitrotyrosine is a stable end product of peroxynitrite oxidation, assessment of its plasma concentration may be useful as a marker of NO-dependent damage in vivo. Since NO<sub>x</sub> is only an indicator for enhanced NO production, protein associated nitrotyrosine might be a more suitable marker for damage induced by reactive nitrogen intermediates derived from NO. Furthermore, most proteins have a longer half-life in the circulation than NO<sub>x</sub> levels. The presence of nitrotyrosine has been detected in various

inflammatory processes including atherosclerotic plaques, celiac disease, rheumatoid arthritis, chronic renal failure and septic shock. In normal plasma low, undetectable, levels of nitrotyrosine are present.

Nitrosylation of the amino acid tyrosine occurs both for free tyrosine and for protein bound tyrosine.

#### 試験プラットフォーム

Microplate

#### 製品の特性

#### 保存方法

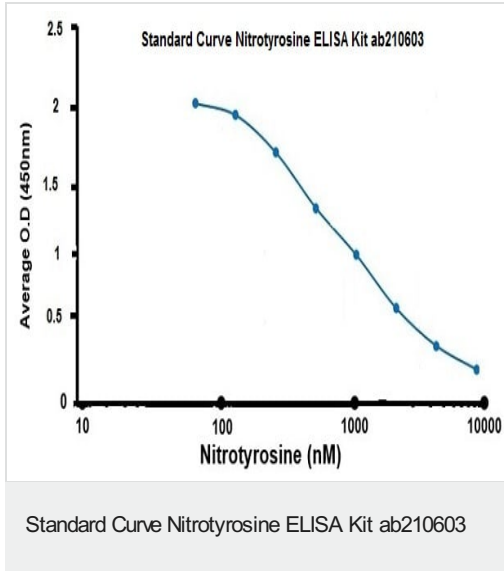
Store at +4°C. Please refer to protocols.

内容	ラベル	1 x 96 tests
Anti-Nitrotyrosine:HRP Conjugated Detection Antibody		1 x 75µl
Nitrosylated BSA Coated Plate		1 x 96 tests
Nitrotyrosine Antibody Diluent	Blue	1 x 13ml
Nitrotyrosine Standard		1 x 110µl
Plate Cover		2 units
Sample and Standard Diluent	Red	1 x 50ml
Stop Solution		1 x 13ml
TMB Substrate		1 x 13ml
Wash Buffer Concentrate (10X)		1 x 50ml

#### 関連性

The cellular production of highly reactive nitrogen species derived from nitric oxide, such as peroxynitrite, nitrogen dioxide and nitryl chloride, leads to the nitration of tyrosine residues in tissue proteins. The extent of protein nitrotyrosine formation provides an index of the production of reactive nitrogen species and potential cell damage over a period of time. Nitrotyrosine can be measured by amino-acid analysis of protein hydrolysates and detected, estimated semi-quantitatively and located in cells and tissues by immunocytochemical techniques using antibodies directed against the nitrotyrosine hapten.

#### 画像



Standard Curve Nitrotyrosine ELISA Kit ab210603

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

### Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.co.jp/abpromise> or contact our technical team.

### Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors