

3-Nitrotyrosine ELISA Kit ab116691

★★★★★ 1 Abreviews 62 References 画像数 1

医薬用外毒物

製品の概要

製品名

3-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%
1	8			6.9%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
1	3			13%

サンプルの種類

Cell culture extracts, Tissue Extracts

アッセイタイプ

Sandwich (quantitative)

検出感度

8 ng/ml

検出範囲

8 ng/ml - 1000 ng/ml

ステップ

Multiple steps standard assay

製品の概要

ab116691 is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of 3-nitrotyrosine in cell and tissue lysates. The assay employs an antibody specific for 3-nitrotyrosine coated on a 96-well plate. Standards and samples are pipetted into the wells and 3-nitrotyrosine present in the sample is bound to the wells by the immobilized antibody. The wells are washed and a biotin labeled anti-3-nitrotyrosine detector antibody is added. After washing away unbound detector antibody, HRP-conjugated streptavidin specific for the biotin labeled detector antibody is pipetted into the wells. The wells are again washed, an HRP substrate solution (TMB) is added to the wells and color develops in proportion to the amount of 3-nitrotyrosine bound. The developing blue color is measured at 600 nm. Optionally the reaction can be stopped by adding hydrochloric acid which changes the color from blue to yellow and the intensity can be measured at 450 nm.

特記事項

Store all components at 4°C. This kit is stable for 6 months from receipt. After reconstitution the standard should be stored at -80°C. Unused microplate strips should be returned to the pouch containing the desiccant and resealed.

試験プラットフォーム

Microplate

製品の特性

保存方法

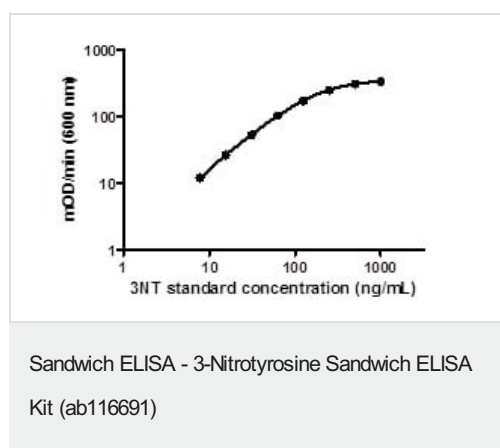
Store at +4°C. Please refer to protocols.

内容	1 x 96 tests
10X 3-nitrotyrosine Detector Antibody	1 x 1ml
10X Blocking Buffer	1 x 6ml
10X HRP Label	1 x 1ml
20X Buffer	1 x 20ml
3-nitrotyrosine BSA standard (Lyophilized)	1 vial
Extraction Buffer (ab260490)	1 x 15ml
Microplate 96 antibody coated wells in 12 strips	1 unit
HRP Development Solution	1 x 12ml

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



Example standard curve.

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