

Nitric Oxide Synthase Activity Assay Kit (Fluorometric) ab211084

★★★★★ [1 Abreviews](#) [4 References](#) [4 图像](#)

概述

产品名称	Nitric Oxide Synthase Activity Assay试剂盒(Fluorometric)
检测方法	Fluorescent
样品类型	Cell Lysate, Tissue Lysate
检测类型	Enzyme activity (quantitative)
灵敏度	0.75 μ U
种属反应性	与反应: Mammals, Other species
产品概述	Nitric Oxide Synthase Activity Assay Kit (Fluorometric) (ab211084) provides an accurate and convenient method to assay nitric oxide synthase (NOS) activity in cell and tissue lysates. In this assay, nitric oxide (NO) generated by NOS undergoes a series of reactions and reacts with the fluorescent probe to generate a stable signal at Ex/Em = 360/450 nm, which is directly proportional to NOS activity. The assay is simple, sensitive and high-throughput adaptable and can detect as low as 0.75 μ U of NOS activity.
说明	<p>This product is manufactured by BioVision, an Abcam company and was previously called K206 Nitric Oxide Synthase (NOS) Activity Assay Kit (Fluorometric). K206-100 is the same size as the 100 test size of ab211084.</p> <p>Nitric oxide synthases (EC 1.14.13.39) (NOS) are a family of enzymes that catalyze the production of nitric oxide (NO) from L-arginine. Nitric oxide (NO) plays an important role in neurotransmission, vascular regulation, immune response and apoptosis. In presence of NADPH, FAD, FMN, (6R)-5,6,7,8-tetrahydrobiopterin, calmodulin and heme, NOS catalyzes a five-electron oxidation of the guanidino nitrogen of L-arginine with molecular oxygen to generate NO and L-citrulline. There are three isoforms of NOS: endothelial (eNOS), neuronal (nNOS), and inducible (iNOS). nNOS accounts for the production of NO in central nervous system, where NO participates in cell communication and information storage. eNOS produces NO in blood vessels and is involved in regulation of vascular function. In contrast to other isoforms, iNOS is expressed de novo under oxidative stress conditions and produces large amounts of NO as a part of body's defense mechanism.</p>
平台	Microplate reader

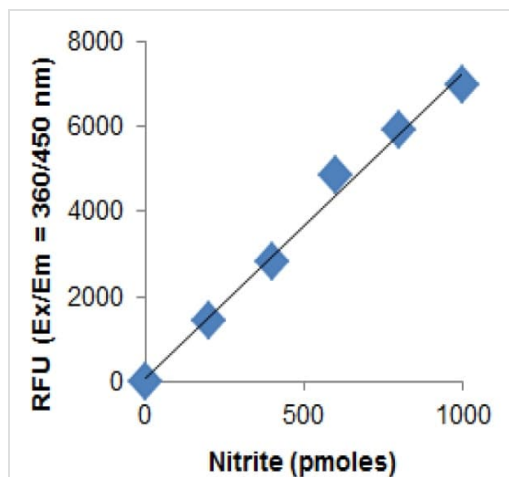
性能

存放说明 Store at -80°C. Please refer to protocols.

组件	100 tests
25X NOS Cofactor 2	1 x 100µl
DAN Probe	1 x 1ml
Enhancer II	1 vial
Nitrate Reductase II	1 vial
Nitrite Standard	1 vial
NOS Assay Buffer	1 x 25ml
NOS Cofactor I	1 vial
NOS Dilution Buffer	1 x 1.5ml
NOS Positive Control	1 x 4µl
NOS Substrate	1 x 500µl
Sodium Hydroxide	1 x 1ml

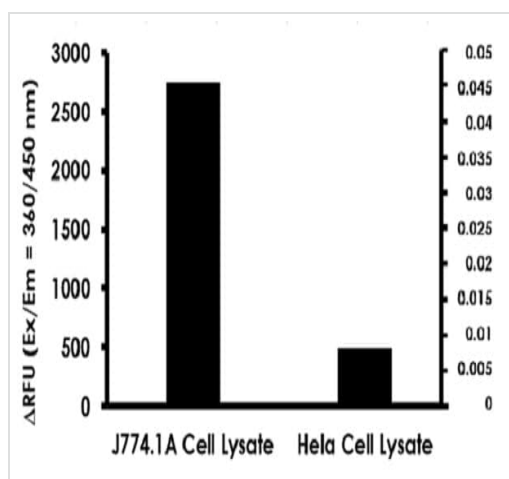
功能	Produces nitric oxide (NO) which is a messenger molecule with diverse functions throughout the body. In the brain and peripheral nervous system, NO displays many properties of a neurotransmitter. Probably has nitrosylase activity and mediates cysteine S-nitrosylation of cytoplasmic target proteins such SRR.
组织特异性	Isoform 1 is ubiquitously expressed: detected in skeletal muscle and brain, also in testis, lung and kidney, and at low levels in heart, adrenal gland and retina. Not detected in the platelets. Isoform 3 is expressed only in testis. Isoform 4 is detected in testis, skeletal muscle, lung, and kidney, at low levels in the brain, but not in the heart and adrenal gland.
序列相似性	Belongs to the NOS family. Contains 1 FAD-binding FR-type domain. Contains 1 flavodoxin-like domain. Contains 1 PDZ (DHR) domain.
结构域	The PDZ domain in the N-terminal part of the neuronal isoform participates in protein-protein interaction, and is responsible for targeting nNos to synaptic membranes in muscles. Mediates interaction with VAC14.
翻译后修饰	Ubiquitinated; mediated by STUB1/CHIP in the presence of Hsp70 and Hsp40 (in vitro).
细胞定位	Cell membrane > sarcolemma. Cell projection > dendritic spine. In skeletal muscle, it is localized beneath the sarcolemma of fast-twitch muscle fiber by associating with the dystrophin glycoprotein complex. In neurons, enriched in dendritic spines.

图片



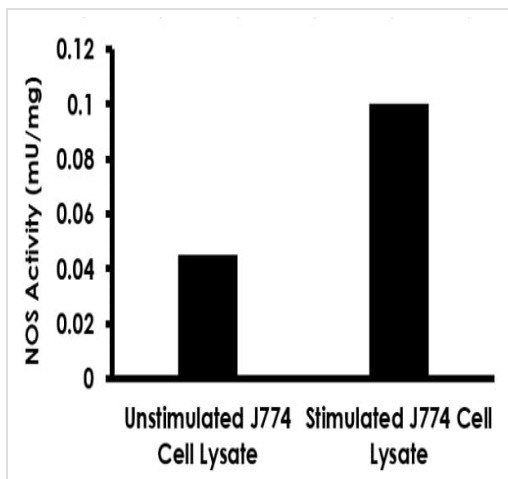
Nitric Oxide Synthase Activity Assay Kit (Fluorometric) (ab211084).
Typical standard calibration curve.

Nitric Oxide Synthase Activity Assay Kit
(Fluorometric) (ab211084)



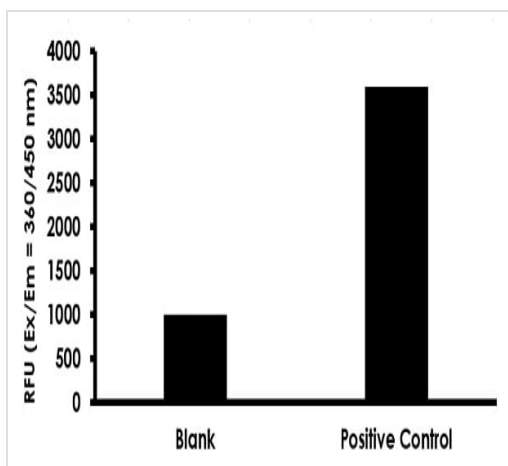
Nitric Oxide Synthase Activity Assay Kit (Fluorometric) (ab211084).
Detection of endogenous NOS activity in cell lysates from J774.1A mouse monocytes (135 μ g) and HeLa cells (157 μ g).

Nitric Oxide Synthase Activity Assay Kit
(Fluorometric) (ab211084)



Nitric Oxide Synthase Activity Assay Kit
(Fluorometric) (ab211084)

Nitric Oxide Synthase Activity Assay Kit (Fluorometric) (ab211084). Detection of endogenous NOS activity in J774.1A cell lysates. J774 cells were stimulated with 200 ng/mL LPS and 100 ng/mL murine IFN-gamma. Unstimulated control included. Cell lysates were prepared (225 µg) and assayed following the kit protocol.



Nitric Oxide Synthase Activity Assay Kit
(Fluorometric) (ab211084)

Nitric Oxide Synthase Activity Assay Kit (Fluorometric) (ab211084). Measurement of NOS Positive Control activity (5 µL) compared to blank (assay buffer).

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.cn/abpromise> or contact our technical team.

Terms and conditions

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