

iNOS Polyclonal Antibody

(Catalog No. A-0475)

Background

Nitric oxide (NO) is a diffusible free radical gas that acts as a biological messenger between cells. Due to its rapid reaction with different substances, it has an exceedingly short half-life of only a few seconds. NO is synthesized in many tissues and affects a variety of physiological functions. Production of NO is mediated by members of the nitric oxide synthase (NOS) family. NOS catalyzes the oxidization of L-arginine to produce L-citrulline and NO. Three known isoforms of these enzymes exist: nNOS, type I (neuronal); eNOS, type III (endothelial cell); and iNOS, type II (inducible). All of these isoforms contain calmodulin, nicotinamide adenine dinucleotide phosphate (NADPH), flavin adenine dinucleotide (FAD), and flavin mononucleotide (FMN) binding domains. Inducible nitric oxide synthase (iNOS), which produce large amounts of NO, is expressed in a broad range of cell types including macrophages, hepatocytes, synoviocytes, and smooth muscle cells. Inflammatory mediators such as cytokines and lipopolysaccharides (LPS) cause an increase in iNOS mRNA, protein, and activity levels. Protein kinase C-stimulating agents also display the same effect on iNOS activity.

Description

Rabbit polyclonal antibody to iNOS

Formulation

Liquid. In PBS containing 50% glycerol and 0.09% sodium azide.

Immunogen

Synthetic peptide corresponding to aa 1131-1144 of mouse iNOS.

Specificity

Mouse, Rat, Bovine

Purification

Protein A-affinity purified.

Storage

Store at -20°C. Avoid multiple freeze/thaw cycles.

Application

IP, WB (1:1000, ECL)

Ordering Information

Product

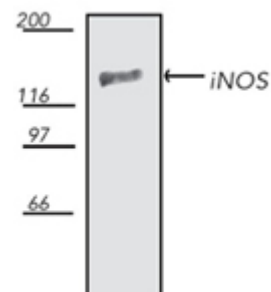
iNOS Polyclonal Antibody

Size

200 µg

Cat. No.

A-0475-200



▲ **Fig. 1.** Western blot analysis: rat brain extract, probed with iNOS Polyclonal Antibody

This product is for research purposes only. Not intended for use in diagnostic procedures.