Nitric oxide synthase (universal) monoclonal antibody (NOS-3F7-B11-B5)

This antibody is covered by our Worry-Free Guarantee.

Citations: 3

View Online »

Ordering Information

Order Online »

ALX-804-029-R200

200µl

Manuals, SDS & CofA

View Online »

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name NOS (universal)

Application IHC (FS), WB

Application NotesDetects bands of ~130kDa (representing iNOS in samples

first induced with interferon (IFN) and lipopolysaccarides (LPS)), ~155kDa (in tissues expressing nNOS) and ~140kDa (in tissues expressing eNOS) by Western blot.

Clone NOS-3F7-B11-B5

Formulation Liquid. Ascites diluted in PBS containing 0.05% sodium

azide.

Host Mouse

Immunogen Purified neuronal nitric oxide synthase (nNOS; NOS I)

from bovine brain.

lsotype lgM

Recommendation Dilutions/Conditions Immunohistochemistry (frozen sections, 1:100)Western

Blot (1:50-1:500)Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be

determined individually for each application.

Species Reactivity Bovine, Mouse, Rat

Specificity Recognizes iNOS (NOS II), eNOS (NOS III) and nNOS

(NOS I).

UniProt ID E1BKI3

Last modified: May 29, 2024



info-

eu@enzolifesciences.com