

p47Phox polyclonal antibody

Originally recognized for their role in the respiratory burst of phagocytes, the NOX family of NADPH oxidases is now known to consist of at least 13 enzymes and enzyme subunits that catalyze the production of superoxide and other downstream reactive oxygen species (ROS) in nearly all cells and tissues. NOX enzymes contribute to a multitude of physiological functions including host defense, cell signaling, and differentiation, with dysregulation linked to pathological manifestations including immunosuppression, hypothyroidism, cardiovascular disease, and neurodegeneration. p47Phox, also known as neutrophil cytosolic factor 1 (NCF1), is considered the organizing subunit of NADPH oxidase enzyme complexes containing the prototypic NOX2 enzyme. Cell stimulation results in phosphorylation of p47Phox, exposing a proline rich region which mediates translocation and association with membrane-associated NOX2/p22 complexes. Translocation of p47Phox is accompanied by the p40Phox subunit and p67Phox activator subunit, which along with GTP-bound Rac completes formation of the active NOX2 oxidase complex. Defects in p47Phox are associated with chronic granulomatous disease (CGD), characterized by an inability of host macrophages and phagocytes to kill ingested microbes, leading to life-threatening infections.

This antibody is covered by our [Worry-Free Guarantee](#).

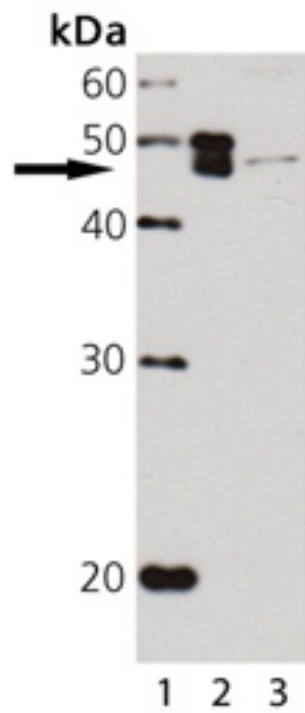
Ordering Information

[Order Online »](#)

ADI-OSA-550-0050	50µg
ADI-OSA-550-0200	200µg

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of p47phox: Lane 1: MWM, Lane 2: p47Phox (human), (recombinant) (His-tag) (Prod No. ADI-OSP-550), Lane 3: RAW 264.7 Cell Lysate.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Neutrophil cytosolic factor 1, NCF1

Application WB

Application Notes Detects a band of ~50kDa by Western blot.

Formulation Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.

Host Rabbit

Immunogen Recombinant human p47Phox.

Purity Detail Protein A-affinity purified.

Recommendation Dilutions/Conditions Western Blot (1:1,000, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Source Purified from rabbit serum.

Species Reactivity Human, Mouse

UniProt ID P14598

Worry-free Guarantee This antibody is covered by our [Worry-Free Guarantee](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)