

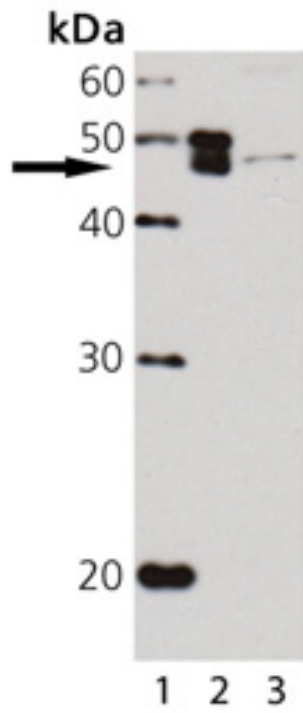
p47Phox polyclonal antibody

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](#).

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

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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