# Nrf2 polyclonal antibody

Nuclear factor erythroid 2-related factor 2 (Nrf2) is a transcription factor, which regulates the expression of many detoxification and antioxidant enzymes. Nrf2 can potentially play a significant role in adaptive responses to oxidative stress. Nrf2 belongs to the Cap N Collar subfamily of basic leucine zipper (CNC-bZIP) transcription factors. The NF-E2 related family comprises of Nrf1 (TCFII), Nrf2, and Nrf3 and these proteins heterodimerize with Maf proteins to bind to Maf recognition element (MARE). Under normal conditions, Nrf2 is inactive and localized in the cytoplasm through its association with Keap1. When cells are exposed to electrophiles or oxidative stress generating agents cytoplasm localization of Nrf2 is inactivated and Nrf2 translocates into the nucleus and binds to specific DNA sequence elements known as NF-E2 binding site, Maf recognition element (MARE), stress response element or anti-oxidant response element (ARE). Studies have investigated the role of Nrf2 in urinary bladder carcinogenesis, pulmonary fibrosis, and smoke-induced emphysema. A study reported Nrf2-mediated ARE transcription to be a likely target of preventative therapy in neurodegenerative disorders such as Huntington's disease.

This antibody is covered by our Worry-Free Guarantee.

Citations: 9

View Online »

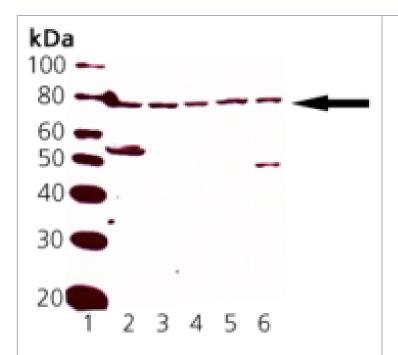
### **Ordering Information**

**Order Online** »

ADI-KAP-TF125-D	50µg
ADI-KAP-TF125-F	200μg

Manuals, SDS & CofA

**View Online** »



Western blot analysis of Nrf2: Lane 1: MW Marker, Lane 2: Rat Liver Microsomes, Lane 3: Mouse Liver Microsomes, Lane 4: Human Liver Microsomes, Lane 5:

Canine Liver Microsomes, Lane 6: HeLa Cell Lysate.

### **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 Nuclear factor erythroid 2-related factor 2

**Application** WB

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

azide.

Gene/Protein Identifier NP\_006155 (RefSeq)

**Host** Rabbit

Immunogen Synthetic peptide derived from the sequence of human

Nrf2 (residues 40-60) with a cysteine residue added and

coupled to KLH.

Purity Detail Protein A affinity purified.

Source Purified from rabbit serum.

Species Reactivity Dog, Guinea pig, Hamster, Human, Mouse, Rat

UniProt ID Q16236

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 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg
Phone: +32 3 466 0420
infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

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