

# HIF-1 $\alpha$ monoclonal antibody (H $\alpha$ 111a)

Hypoxia inducible factor-1 (HIF-1) is a transcriptional complex which regulates systemic, local and intracellular homeostatic responses elicited by hypoxia. HIF-1 is a heterodimer of HIF-1 $\alpha$  and HIF-1 $\beta$ . HIF-1 $\beta$  is constitutively expressed and serves as a dimerization partner for several other transcription factors, while HIF-1 $\alpha$  and the closely related HIF-2 are unique to HIF-1 and their expression is tightly regulated by cellular oxygen concentration. HIF-1 regulates a diverse group of genes, including erythropoietin, VEGF, glucose transporters, heme oxygenase, and NOS.

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

Citations: 4

[View Online »](#)

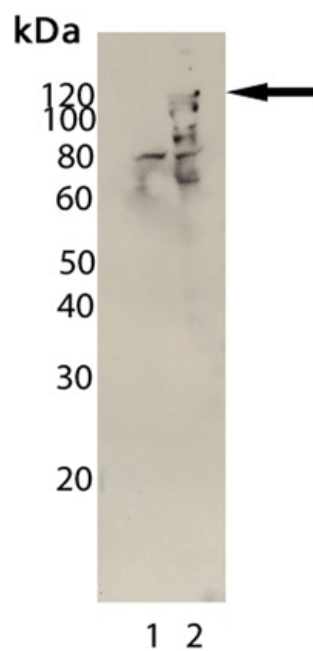
## Ordering Information

[Order Online »](#)

ADI-OSA-602-E	100 $\mu$ g
---------------	-------------

## Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of HIF-1 $\alpha$ , mAb (H $\alpha$ 111a) (Prod. No. ADI-OSA-602): Lane 1: HeLa Cell Lysate (Prod. No. ADI-LYC-HL100), Lane 2: HeLa+CoCl<sub>2</sub> Cell Lysate.

## Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Hypoxia-inducible factor 1 $\alpha$ , ARNT-interacting protein

**Application** IHC, WB

**Application Notes** Detects a band of ~120kDa by Western blot.

**Clone** Halpha111a

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

**Host** Mouse

**Immunogen** Recombinant human HIF-1 $\alpha$ .

**Isotype** IgG2a

**Purity Detail** Protein G affinity purified.

**Recommendation Dilutions/Conditions** Western Blot (4 $\mu$ g/ml) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

**Source** Purified from hybridoma tissue culture supernatant.

**Species Reactivity** Human

**Technical Info / Product Notes** Recommended by the Human Protein Atlas Organization for IHC (Ensembl No. ENSG00000100644).

**UniProt ID** Q16665



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

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

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

France  
Phone: +33 1 42 44 55 55  
[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)

**Worry Free Guarantee** This antibody is covered by our **Worry Free Guarantee**