

HIF-1 alpha

Western Blot

<https://www.abcam.com/protocols/general-western-blot-protocol>

Target Overview in WB Application

HIF-1 alpha is stabilized only at O₂ concentration below 5%. Under normoxic conditions HIF-1 alpha has a short half-life and may be degraded within 5-8 minutes in both nuclear and cytoplasmic compartments. Therefore, **proper sample preparation** is critical to aiding WB success. If care hasn't been taken with sample preparation no bands may be seen on your blot.

Here are a few tips to help ensure the best results in WB:

Sample Preparation	<ul style="list-style-type: none"> • Add adequate protease inhibitors (or phosphatase inhibitors for proteins modified by phosphorylation) to avoid target protein degradation.
	<ul style="list-style-type: none"> • Ultrasonicate samples to enrich more target proteins.
	<ul style="list-style-type: none"> • Keep samples on ice during the whole WB process.
	<ul style="list-style-type: none"> • Perform a Bradford assay, a Lowry assay or a bicinchoninic acid (BCA) assay to determine the protein concentration.
Electrophoresis	<ul style="list-style-type: none"> • For large proteins (the MW of target protein >100 kDa), be sure to run samples in 8% or lower separating gel.
	<ul style="list-style-type: none"> • Load at least 50µg total protein per lane.
	<ul style="list-style-type: none"> • We strongly recommend the use of a positive control lysate when setting up a new experiment; this will give you immediate confidence in the protocol.
Transferring	<ul style="list-style-type: none"> • It is preferred to add SDS to a final concentration of 0.1% in the transfer buffer for large proteins.
	<ul style="list-style-type: none"> • Wash PVDF membrane to remove methanol completely.
	<ul style="list-style-type: none"> • To determine if the transfer is successful by visualization of proteins in membranes using Ponceau S.

You should pay attention to these notes to maximize the signal:

- ✓ **Hypoxic chambers** may be used to incubate samples overnight at low oxygen pressure to induce HIF-1 alpha levels.
- ✓ Cells should be lysed as quickly (within 2 mins) as possible if removed from hypoxia.
- ✓ Use **positive control samples** such as nuclear lysates of DFO or CoCl₂ treated cells.
- ✓ **Overnight incubation** at 4°C with the primary antibody can also help.
- ✓ The observed band size of HIF-1 alpha is not exactly as predicted 93 kDa in WB due to the different forms of HIF-1 alpha as blow:
 - **40-80 kDa** - degraded HIF-1 alpha
 - **110-130 kDa** - post translationally modified HIF-1 alpha
 - **~200 kDa** - heterodimer with HIF-1 beta

Protein Function	<p>Functions as a master transcriptional regulator of the adaptive response to hypoxia. Under such conditions, it can activate over 40 genes whose protein products increase oxygen delivery or facilitate metabolic adaptation to hypoxia.</p> <p style="text-align: right;">SwissProt: Q16665</p>				
Expression	<ul style="list-style-type: none"> - Most cell lines and tissues (under hypoxic conditions) - Highest levels in kidney and heart - Overexpressed in majority of human cancers <p>Under normoxic conditions, HIF-1 alpha is largely undetectable. Hypoxia needs to be induced in most cells and normal tissues.</p>				
Location	<p>Cytoplasmic in normoxic conditions. Nuclear translocation in response to hypoxia.</p> <p style="text-align: right;">PMID: 9822602</p>				
Isoforms	<ul style="list-style-type: none"> - Isoform 1: 93 kDa (predicted) - Isoform 2: 83 kDa (predicted) - Isoform 3: 96 kDa (predicted) <p>The observed band size of HIF-1 alpha may not be the same as predicted MWs in WB due to the different forms of HIF-1 alpha.</p>				
Modifications	<p>Snitrosylation/Phosphorylation/Sumoylation/ Acetylation/Polyubiquitnation/Hydroxyalation</p> <p>The observed band size of HIF-1 alpha may not be the same as predicted MWs in WB due to these modifications.</p>				
Positive Controls	<p>WB: Hypoxic samples such as HeLa-DFO treated whole cell lysate ab116322.</p> <p>For a stronger signal, HeLa-DFO treated nuclear extracts are recommended ab180880. The cell fractionation kit can also be purchased separately ab109719.</p>				
Negative Controls	<p>Most normal tissues or cells, other than kidney or heart.</p>				
Treatments	<p>CoCl₂ or DFO to induce hypoxia.</p> <p style="text-align: right;">PMID: 3217626</p> <p>Protease inhibitors like MG132 can stabilize HIF-1 alpha.</p> <p style="text-align: right;">PMID: 19347037</p>				
References	Journal	abID	Application	Species	PMID
	Nature Communications (2019)	ab1	WB	Mouse	31391533
	Cell (2019)	ab179483	WB	Mouse	31708126
	Science Advances (2019)	ab2185	WB	Mouse	31281892