

HIF-1 α monoclonal antibody (H α 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 α and HIF-1 β . HIF-1 β is constitutively expressed and serves as a dimerization partner for several other transcription factors, while HIF-1 α 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 »](#)

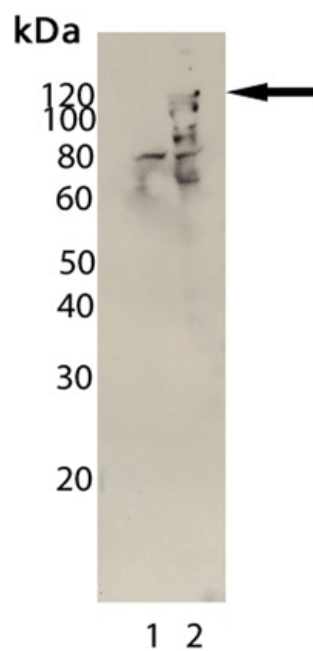
Ordering Information

[Order Online »](#)

ADI-OSA-602-E	100 μ g
---------------	-------------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of HIF-1 α , mAb (H α 111a) (Prod. No. ADI-OSA-602): Lane 1: HeLa Cell Lysate (Prod. No. ADI-LYC-HL100), Lane 2: HeLa+CoCl₂ Cell Lysate.

Long Term Storage -20°C

Regulatory Status RUO - Research Use Only

Alternative Name	Hypoxia-inducible factor 1α, ARNT-interacting protein
-------------------------	---

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

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

Immunogen	Recombinant human HIF-1α.
------------------	---------------------------

Purity Detail Protein G affinity purified.

Recommendation Dilutions/Conditions	Western Blot (4µg/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
1	1
2	2
3	3
4	4
5	5
6	6
7	7
8	8
9	9
10	10
11	11
12	12
13	13
14	14
15	15
16	16
17	17
18	18
19	19
20	20
21	21
22	22
23	23
24	24
25	25
26	26
27	27
28	28
29	29
30	30
31	31
32	32
33	33
34	34
35	35
36	36
37	37
38	38
39	39
40	40
41	41
42	42
43	43
44	44
45	45
46	46
47	47
48	48
49	49
50	50
51	51
52	52
53	53
54	54
55	55
56	56
57	57
58	58
59	59
60	60
61	61
62	62
63	63
64	64
65	65
66	66
67	67
68	68
69	69
70	70
71	71
72	72
73	73
74	74
75	75
76	76
77	77
78	78
79	79
80	80
81	81
82	82
83	83
84	84
85	85
86	86
87	87
88	88
89	89
90	90
91	91
92	92
93	93
94	94
95	95
96	96
97	97
98	98
99	99
100	100

Technical Info / Product Notes

Recommended by the Human Protein Atlas Organization for IHC (Ensembl No. ENSG00000100644).

UniProt ID Q16665

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-be@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