

HIF-1 α monoclonal antibody (H1 α 67)

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

Citations: 1

[View Online »](#)

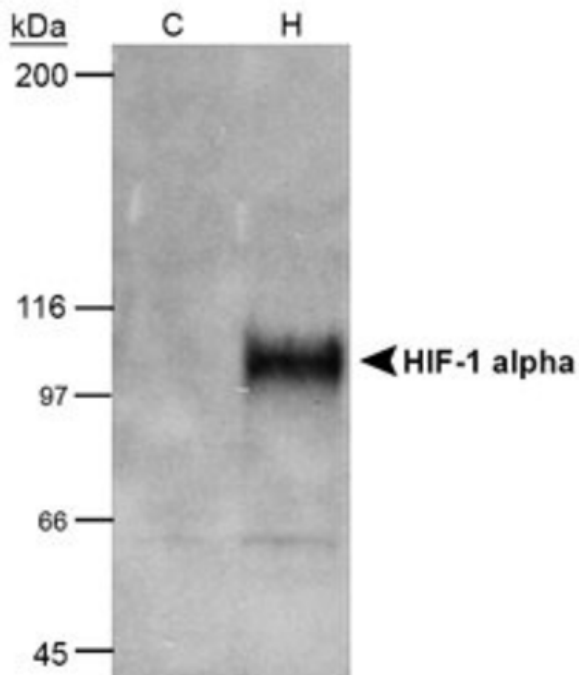
Ordering Information

[Order Online »](#)

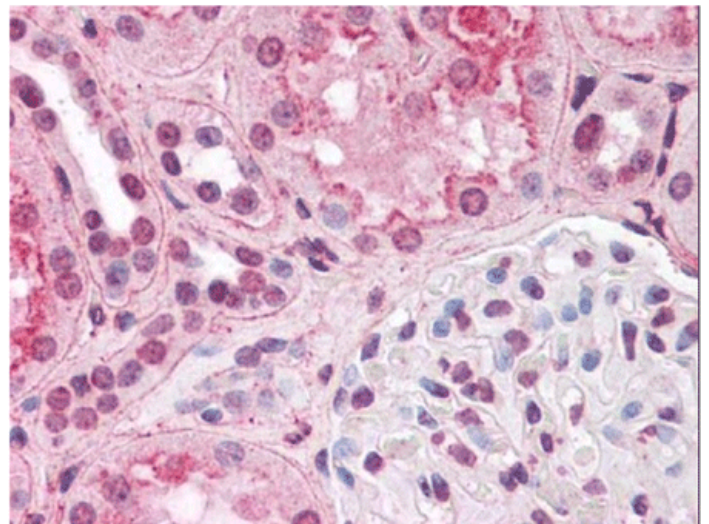
BML-SA287-0100	100 μ l
----------------	-------------

Manuals, SDS & CofA

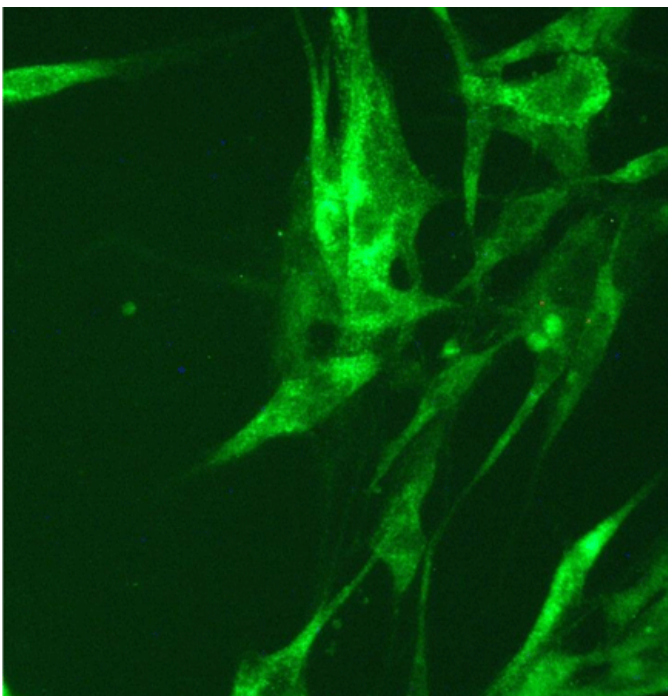
[View Online »](#)



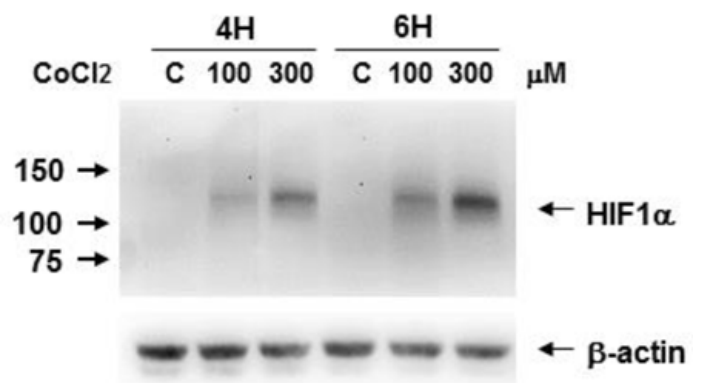
Western Blot: HIF-1 alpha, mAb (H1alpha67) [Prod.no. BML-SA287] – Analysis of 50ug cobalt chloride (CoCl₂) induced COS-7 nuclear extracts.



Immunohistochemistry: HIF-1 alpha, mAb (H1alpha67) [Prod. no. BML-SA287] – Staining of HIF1 alpha in human kidney. Renal tubular epithelium showed moderate membranous, cytoplasmic and nuclear staining, and glomeruli showed faint to moderate nuclear staining.



Immunocytochemistry/Immunofluorescence: HIF-1 alpha, mAb (H1alpha67) [Prod. no. BML-SA287] – Expression of HIF-1 in interstitial pig cells under 12% of O₂ condition.



Western Blot: HIF-1 alpha, mAb (H1alpha67) [Prod.no BML-SA287] – HIF-1 alpha induction by CoCl₂ on Caki-1 cell lysate.

Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Hypoxia-inducible factor 1 α , ARNT-interacting protein
Application	ELISA, Flow Cytometry, ICC, IF, IHC, IHC (FS), IHC (PS), IP, WB
Application Notes	Detects a band of ~120kDa by Western blot.
Clone	H1 α 67
Formulation	Liquid. In PBS with 1% BSA containing 0.05% sodium azide.
Host	Mouse
Immunogen	Recombinant human hypoxia-inducible factor-1 α (HIF-1 α) (aa 432-528).
Isotype	IgG
Purity Detail	Purified IgG.
Species Reactivity	Bovine, Ferret, Human, Monkey, Mouse, Porcine, Rabbit, Rat, Sheep
UniProt ID	Q16665
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee .



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

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

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

France
Phone: +33 472 440 655
[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)