

AhR polyclonal antibody

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

Citations: 19

[View Online »](#)

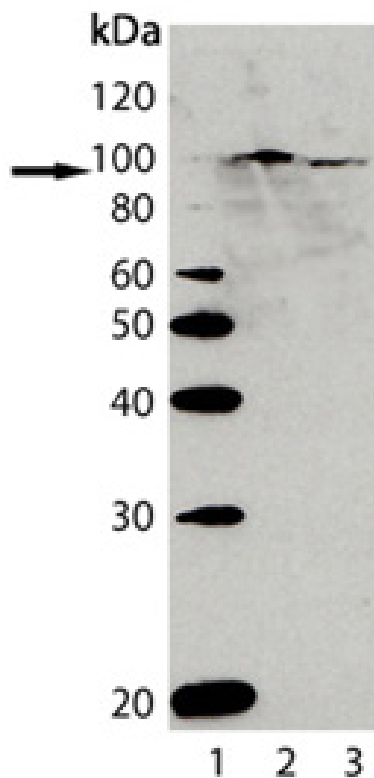
Ordering Information

[Order Online »](#)

BML-SA550-0025	25µl
BML-SA550-0100	100µl

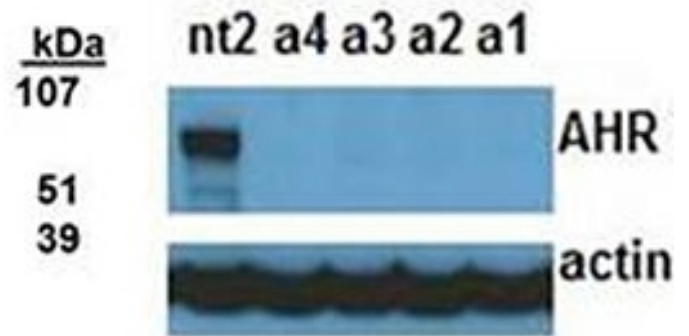
Manuals, SDS & CofA

[View Online »](#)



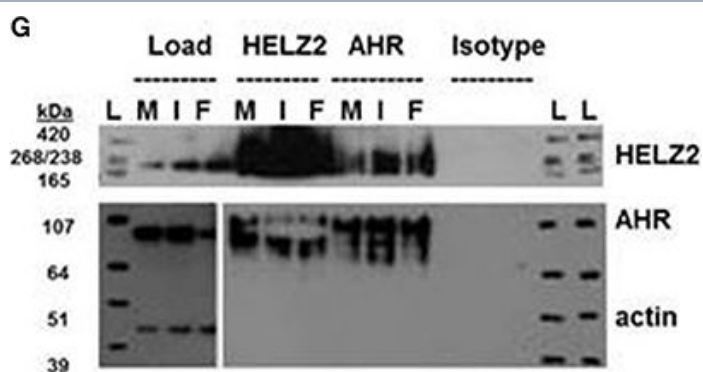
Western blot analysis of AhR, pAb (Prod. No. BML-SA550): Lane 1: MW marker; Lane 2: Rat liver lysate; Lane 3: Mouse liver lysate.

C



AHR is an IFN Effector Gene and Interacts with HELZ2. RNAi-mediated knockdown of AHR with four unique siRNAs leads to weak but consistent rescue of DENV from IFN in HeLa cells (A) and Huh7.5.1 hepatocytes (B). (C) Knockdown of AHR by the four unique siRNAs from (A) was validated at the protein level using western blot (C) with quantification (D) and at the mRNA level using qRT PCR (E). (F) Co-immunoprecipitation of AHR with HELZ2 (H) and AHR (A) antibodies and isotype control (I) was performed in Huh7.5.1 cells treated with IFN 1,600 IU/ml \times 24 h or mock, followed by immunoblot for AHR. Load controls (Ld) were run in parallel, and immunoblotted for AHR and actin (G). Co-immunoprecipitation was repeated in the setting of mock (M), IFN 1,600 IU/ml (I), or the AHR agonist FICZ 0.9 μ M (F) \times 24 h, using HELZ2 and AHR antibodies for IP, followed by HELZ2 and AHR immunoblot. Load controls were stained for HELZ2, AHR, and actin. Ladder is indicated by L. * $p < 0.05$ compared to control.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: HELZ2 Is an IFN Effector Mediating Suppression of Dengue Virus. *Front Microbiol* (2017)



AHR is an IFN Effector Gene and Interacts with HELZ2. RNAi-mediated knockdown of AHR with four unique siRNAs leads to weak but consistent rescue of DENV from IFN in HeLa cells (A) and Huh7.5.1 hepatocytes (B). (C) Knockdown of AHR by the four unique siRNAs from (A) was validated at the protein level using western blot (C) with quantification (D) and at the mRNA level using qRT PCR (E). (F) Co-immunoprecipitation of AHR with HELZ2 (H) and AHR (A) antibodies and isotype control (I) was performed in Huh7.5.1 cells treated with IFN 1,600 IU/ml \times 24 h or mock, followed by immunoblot for AHR. Load controls (Ld) were run in parallel, and immunoblotted for AHR and actin (G). Co-immunoprecipitation was repeated in the setting of mock (M), IFN 1,600 IU/ml (I), or the AHR agonist FICZ 0.9 μ M (F) \times 24 h, using HELZ2 and AHR antibodies for IP, followed by HELZ2 and AHR immunoblot. Load controls were stained for HELZ2, AHR, and actin. Ladder is indicated by L. * $p < 0.05$ compared to control.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: HELZ2 Is an IFN Effector Mediating Suppression of Dengue Virus. *Front Microbiol* (2017)

Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Aryl hydrocarbon receptor
Application	IP, WB
Application Notes	Detects a band of ~95kDa by Western blot.
Formulation	Liquid. Neat serum.
Host	Rabbit
Immunogen	Recombinant B-1 AhR (aa 1-402).
Source	Whole rabbit serum.
Species Reactivity	Mouse, Rat
UniProt ID	P30561 (mouse)
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)