

Aryl hydrocarbon receptor polyclonal antibody

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

Citations: 97

[View Online »](#)

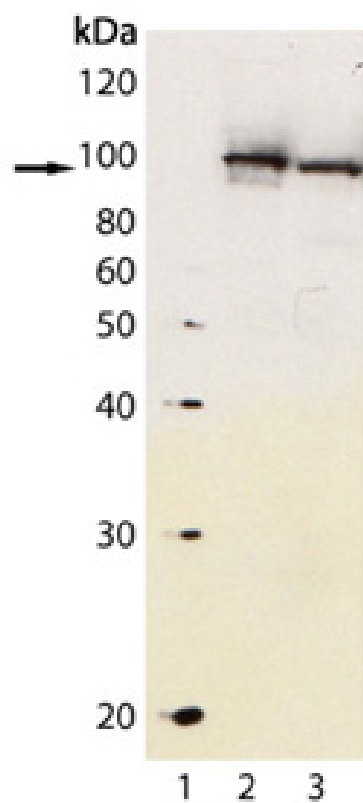
Ordering Information

[Order Online »](#)

BML-SA210-0025	25µg
BML-SA210-0100	100µg

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of Aryl hydrocarbon receptor, pAb (Prod. No. BML-SA210): Lane 1: MW marker; Lane 2: Mouse liver extract; Lane 3: Rat liver extract.

Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	AhR
Application	ChIP, IP, WB
Application Notes	Detects a band of ~95 kDa by Western blot.
Crossreactivity	Does not cross-react with Arnt.
Formulation	Liquid. In TBS containing 0.02% sodium azide.
Host	Rabbit
Immunogen	Recombinant protein corresponding to aa 1-402 of the N-terminal B-1 AhR (Aryl hydrocarbon receptor).
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1:2,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from rabbit serum.
Species Reactivity	Mouse
Technical Info / Product Notes	Cited samples: For an overview on cited samples please click here.
UniProt ID	P30561

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