

Aryl hydrocarbon receptor monoclonal antibody (RPT9)

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

Citations: 1

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	AhR
Application	ChIP, Flow Cytometry, Gel Shift, ICC, IF, IHC (FS), IP, WB
Application Notes	Detects a band of ~95kDa by Western blot. Cited Applications include: ELISA
Clone	RPT9
Formulation	Liquid. Contains 0.05% sodium azide.
Host	Mouse
Immunogen	Synthetic peptide corresponding to aa 12-17 (R12KRRKP17) and aa 22-31 (V22KPIPAEGIK31) of mouse AhR (aryl hydrocarbon receptor).
Isotype	IgG3
Purity Detail	Protein A-affinity purified.
Recommendation Dilutions/Conditions	Chromatin Immunoprecipitation (1:100)Flow Cytometry (1:50)Immunofluorescence (1:10-1:100)Western Blot (1:500 – 1:2,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse, Rat
UniProt ID	P30561

Last modified: May 29, 2024



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)