

MRP4 monoclonal antibody (M4I-80)

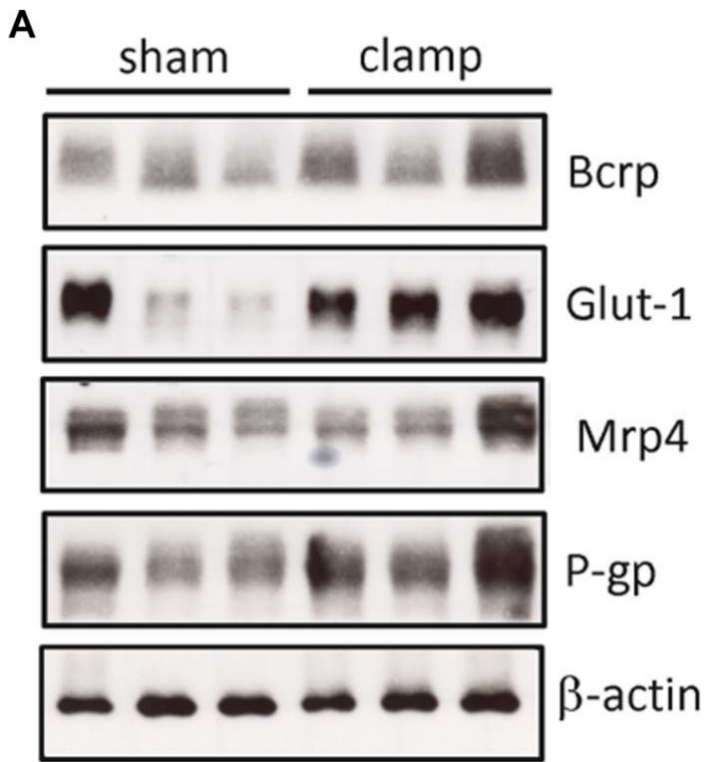
This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 8

[View Online »](#)

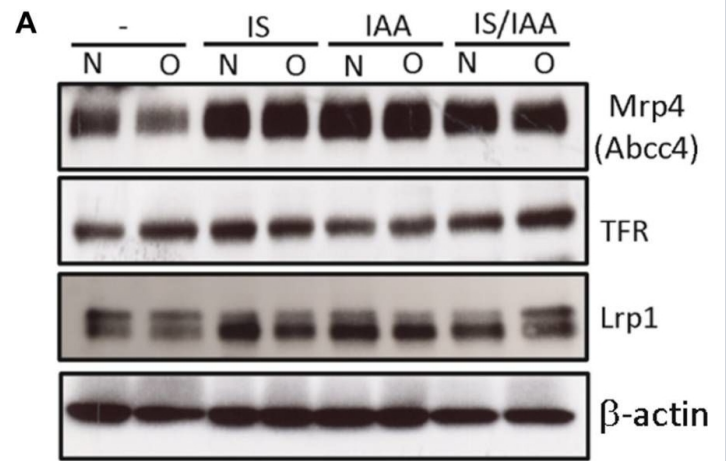
Manuals, SDS & CofA

[View Online »](#)



Expression of selected transporters and tight junction proteins in brain capillaries of mice exposed to bilateral renal ischemia/reperfusion injury at the protein level. Brain capillaries were isolated from frozen brains of mice exposed to bilateral renal ischemia/reperfusion injury followed by protein extraction and Western blot (A). The protein content of Bcrp (Abcg2, ATP Binding Cassette Subfamily G Member 2) (B), Glut1 (Slc2a1, Solute Carrier Family 2 Member 1) (C), Mrp4 (Abcc4, ATP Binding Cassette Subfamily C Member 4) (D), and P-gp (Mdr1, Abcb1b ATP Binding Cassette Subfamily B Member1) (E) was normalized to β -actin and is shown as average of $n = 3$ with standard deviations. * $p < 0.05$, *** $p < 0.001$.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Kidney Ischemia/Reperfusion Injury Induces Changes in the Drug Transporter Expression at the Blood-Brain Barrier in vivo and in vitro. *Front Physiol* (2020)



Effects of ischemic HK-2 and uremic toxins on efflux pumps and cellular receptor protein level in brain microvascular endothelial cells. BMECs treated as described in Figure 1 were lysed and used for protein extraction. The protein content of Mrp4 (Abcc4, ATP Binding Cassette Subfamily C Member 4), Tfr (Transferrin Receptor), and Lrp1 (LDL Receptor Related Protein 1) was measured by Western blot (A). The protein content was normalized to β -actin and to the untreated control. The densitometric analysis is shown as an average of four independent experiments with standard deviations (B–D). One-way ANOVA with Dunnett's or Sidak's multiple comparison test was used to compare each treatment to the untreated control. * $p < 0.05$.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: Kidney Ischemia/Reperfusion Injury Induces Changes in the Drug Transporter Expression at the Blood-Brain Barrier in vivo and in vitro. *Front Physiol* (2020)

Handling & Storage

Short Term Storage +4°C

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Multidrug resistance-associated protein 4, ABCC4
Application	ICC, IHC (FS), WB
Application Notes	Immunocytochemistry: cytospin preparations
Clone	M4I-80
Formulation	Liquid. In serum free culture supernatant containing 0.7% BSA and 0.1% sodium azide.
Host	Rat
Immunogen	Recombinant human MRP4 (multidrug resistance-associated protein 4) (aa 372-431).
Isotype	IgG2a
Recommendation Dilutions/Conditions	Immunocytochemistry: Use 1:20-1:50 dilution on acetone fixed cytospin preparations. Immunohistochemistry (frozen sections): Use 1:20 dilution on acetone fixed frozen sections followed by incubation with biotin-labelled rabbit anti-rat IgG and HRP-labelled streptavidin. Western Blot: Use 1:20-1:50 dilution and HRP-anti-rat-IgG. Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse
UniProt ID	O15439

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)