

TRIP-Br1 monoclonal antibody (TB1-27/2)

TRIP-Br1 (SEI-1) functions as physiological integrator of transcriptional regulatory signals at E2F-responsive promoters to ensure proper execution of E2F-dependent cell cycle progression. TRIP-Br1 is defined as an oncogene.

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

Citations: 12

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-645-C050	50µg
ALX-804-645-C100	100µg

Manuals, SDS & CofA

[View Online »](#)

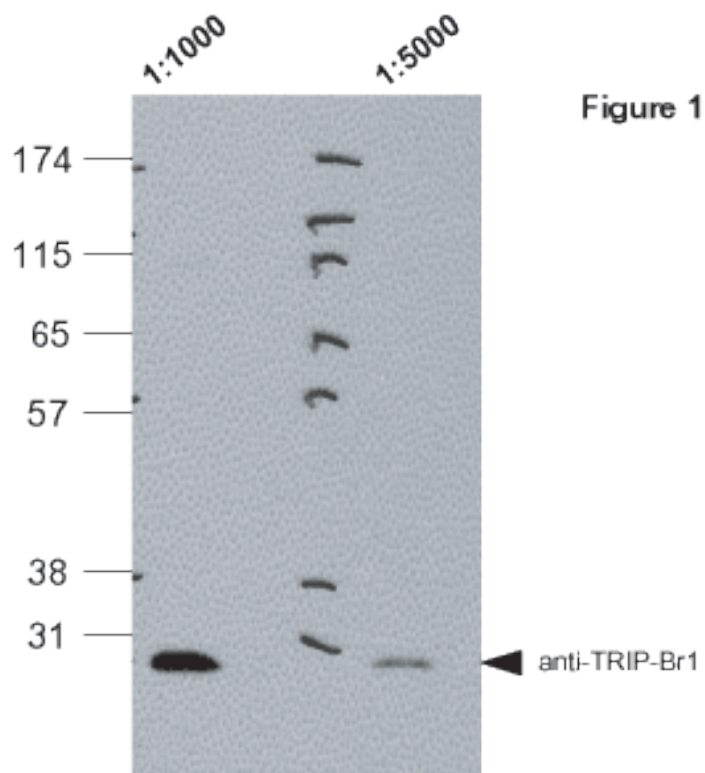


Figure 1: Western blot analysis of pig kidney lysates (LLC-PK1) using MAb to TRIP-Br1 (TB1-27/2) (Prod. No. ALX-804-645) at 1:1'000 and 1:5'000.

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	Transcriptional regulator interacting with the PHD-bromodomain 1, SEI-1, SERTA domain-containing protein 1, CDK4-binding protein p34SEI1
Application	IHC, IP, WB
Clone	TB1-27/2
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Mouse
Immunogen	Synthetic peptide corresponding to aa 27-42 (D27PGHAAVAQAPPAVAS42) of human TRIP-Br1 [SEI-1].
Isotype	IgG2a
Recommendation Dilutions/Conditions	Immunohistochemistry (1:200)Western Blot (1:1,000-1:5,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from hybridoma tissue culture supernatant.
Species Reactivity	Human, Monkey, Mouse, Porcine
UniProt ID	Q9UHV2

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