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 »

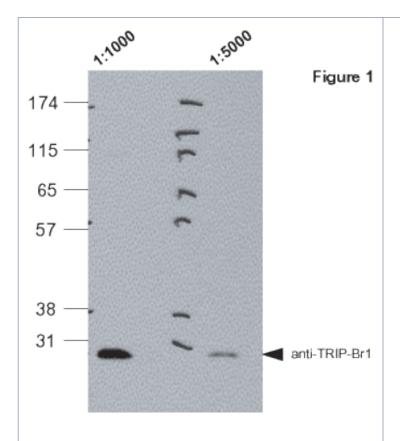


Figure 1: Western blot analysis of pig kidney lysates (LLC-PK1) using MAb to TRIP-Br1 (TB1-27/2) (Prod. No. ALX-804-645) at 1:1'000 and 1:5'000.

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

Last modified: May 29, 2024



info-

eu@enzolifesciences.com