Rictor monoclonal antibody (1G11)

The mTOR (mammalian target of rapamycin) pathway coordinates nutrient signals with growth factor dependent signaling. mTOR is organized in two independent protein complexes, mTORC1 and mTORC2. The mTORC1 complex containing mTOR, GβL/mLST8 and Raptor is sensitive to rapamycin, while the mTORC2 complex containing mTOR, GβL/mLST8, SIN1, and Rictor is rapamycin insensitive. The mTOR/Rictor complex phosphorylates numerous substrates associated with regulation of the actin cytoskeleton, and is also responsible for phosphorylation of Akt/PKB at Ser473, a critical step in Akt activation.

This antibody is covered by our Worry-Free Guarantee.

Citations: 1

View Online »

Ordering Information

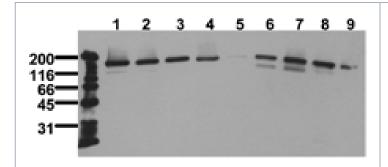
Order Online »

ADI-905-766-100

100µg

Manuals, SDS & CofA

View Online »



Western blot analysis of serum starved cancer cells probed with Rictor mAb (1G11); HeLa (1), HepG2 (2), HEK293 (3), SH-SY5Y (4), MDCK (5), PC12 (6), CTM93 (7) Neuro 2A (8), NIH-3T3 (9).

Handling & Storage

Use/Stability Stable at -80°C up to 1 year, at 4°C up to 3 months.

Handling Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -80°C.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application WB

Application NotesDetects a band of ~190kDa by Western blot.

Clone 1G11

Formulation Lyophilized from 1ml of 2x PBS containing 0.09% sodium

azide, PEG, and sucrose.

Host Mouse

Immunogen Synthetic peptide corresponding to a portion of Rictor.

lgG1

Purity Detail Thiophilic adsorption and size exclusion chromatography

purified.

Recommendation Dilutions/Conditions Western Blot (0.5µg/ml, ECL)Suggested

dilutions/conditions may not be available for all applications. Optimal conditions must be determined

individually for each application.

Reconstitution Reconstitute with 1ml water (15 minutes at room

temperature).

Species Reactivity Human, Mouse, Rat

UniProt ID Q6R327

Last modified: May 29, 2024



info-

eu@enzolifesciences.com