

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.

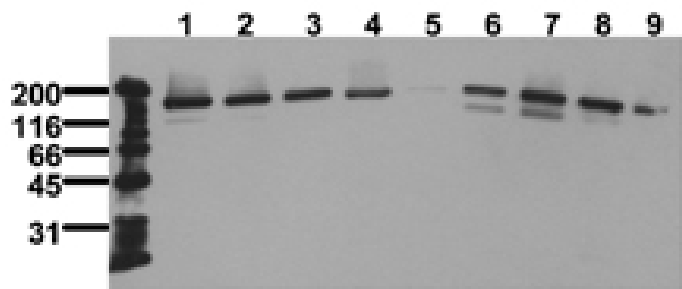
This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 1

[View Online »](#)

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 Notes	Detects 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.
Isotype	IgG1
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



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