

Raptor monoclonal antibody (10E10)

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.

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

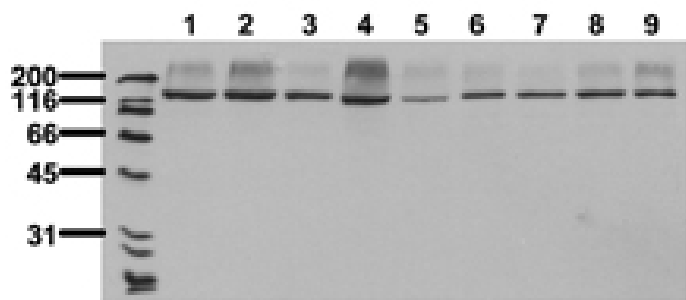
Ordering Information

[Order Online »](#)

ADI-905-765-100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of serum starved cancer cells, probed with Raptor mAb (10E10); HeLa (1); HepG2 (2); HEK293 (3); SH-SY5Y (4); MDCK (5); PC12 (6); CMT-93 (7); Neuro 2A (8); and 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 ~160kDa by Western blot.
Clone	10E10
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 Raptor.
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	Dog, Human, Mouse, Rat
UniProt ID	Q8N122

Worry-free Guarantee



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)

This antibody is covered by our **Worry-Free Guarantee**.