

Product Information

Description:	Recombinant human full length RPTOR (regulatory-associated protein of mTOR). This construct contains an N-terminal Strep-tag. This protein was affinity purified.
Background:	mTOR (mammalian target of rapamycin) is a serine/threonine kinase protein of the PIK3 (phosphatidylinositol 3- kinase)-related family. It can form two different complexes, which differ in the accessory protein present and have different biological functions. mTORC1 (mTOR complex 1) includes the protein RPTOR, while mTORC2 is composed of RICTOR (RPTOR independent companion of mTOR complex 2). MLST8 (target of rapamycin complex subunit LST8) is common to both complexes and seems to stabilize the active site of mTOR and participate in its activation. mTORC1 is activated in response to growth factors and responds to activation of the PI3K-AKT (protein kinase B) pathway. It is involved in cell growth via phosphorylation of proteins like 4E-BP1 (eIF-4E-binding protein). Dysfunction of mTORC1 can result in cancer, and strategies targeting mTORC1 may result in major advances in cancer therapy.
Species:	Human
Construct:	RPTOR (Strep-2-1335(end))
Concentration:	0.28 mg/ml
Expression System:	Sf9
Purity:	88%
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
MW:	150 kDa
Genbank Accession:	NM_020761
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
Applications:	Useful for SDS-PAGE.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining

