

# mTOR/Raptor/MLST8 Complex, FLAG-Tag, His-Tag (Human) Recombinant

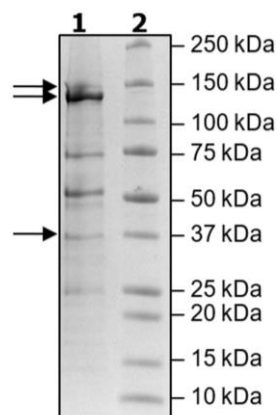
Catalog: 40300  
Lot: 220516

## Product Information

<b>Construct:</b>	mTOR (FLAG-1362-2549(end)) / Raptor (His-2-1335(end)) / MLST8 (His-2-326(end))
<b>Concentration:</b>	0.13 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, and 80 ng/μl FLAG peptide
<b>Expression System:</b>	Sf9
<b>Format:</b>	Aqueous buffer solution.
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>Genbank Accession:</b>	mTOR: NM_004958; Raptor: NM_020761; MLST8: NM_022372
<b>MW:</b>	mTOR: 137 kDa; Raptor: 150 kDa; MLST8: 37 kDa
<b>Purity:</b>	63%
<b>Specific Activity:</b>	6.8 pmol/min/μg
<b>Assay Conditions:</b>	Activity was measured using ADP-Glo™ (Promega) reagents with 100 μM ATP, 0.1 mg/ml 4EBP1 and 5 mM MnCl <sub>2</sub> in 25 μl using kinase buffer (40 mM Tris HCl, pH 7.4, 20 mM MgCl <sub>2</sub> , 0.1 mg/ml BSA, and 2 mM DTT). Reaction was incubated for 45 minutes at 30°C.
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

4-20% SDS-PAGE Coomassie Staining



mTOR/Raptor/MLST8 Complex Activity

