

N17 Rac1-GST Mutant Protein
(Human recombinant)
Cat. # R17G01

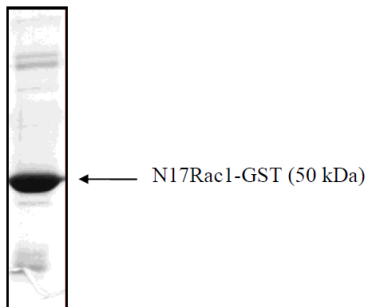
Upon arrival store at 4°C (desiccated)
See datasheet for storage after reconstitution

Material

The dominant negative form of the human Rac1 protein has been produced in a bacterial expression system. The protein has a threonine to asparagine substitution at amino acid 17 creating a dominant negative mutant protein. The protein contains a GST protein tag (~28 kDa) at its amino terminus. The protein is supplied as a lyophilized white powder. The approximate Molecular weight of N17 Rac1-GST is 50 kDa. When it is reconstituted in distilled water to 1 mg/ml, the protein will be in the following buffer: 20 mM Tris pH 7.5, 0.4 mM MgCl₂, 0.4% sucrose and 0.1% dextran.

Purity

Purity is determined by scanning densitometry of proteins on SDS-PAGE gels. Samples are ~80% pure. The gel below shows 10 µg of N17 Rac1-GST stained with Coomassie Blue.



Storage

Stable at 4°C for 1 year as a lyophilized powder. Once reconstituted as a 1 mg/ml solution the protein can be stored at -70°C for 6-8 months. For frozen storage we recommend that the protein be supplemented with DTT to 1 mM final concentration, aliquoted into "experiment sized" amounts and snap frozen in liquid nitrogen. We do not recommend freeze thawing of the proteins.

Product Citations/Related Products

For the latest citations and related products please visit www.cytoskeleton.com.