

His-Ran Protein
(Human recombinant)
Cat. # RN01

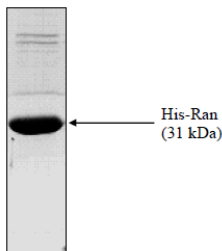
Upon arrival store at 4°C (desiccated)
See datasheet for storage after reconstitution

Material

The human Ran protein has been produced in a bacterial expression system. The protein contains a six amino acid histidine tag at its amino terminus. The protein is supplied as a lyophilized powder. When it is reconstituted in distilled water to 1 mg/ml, the protein is in the following buffer: 1.4 mM Tris pH 7.5, 0.28 mM MgCl₂, 0.28% sucrose, 0.07% dextran.

Purity

Purity is determined by scanning densitometry of proteins on SDS-PAGE gels. Samples are >95% pure. The gel below shows 10 µg of His-Ran stained with Coomassie blue. The protein is 31kD (but will run between 30-36 kDa on a 12% acrylamide gel).



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 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 protein.

Product Citations/Related Products

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