

RhoGDI-GST Protein (human recombinant)

Cat. # GDI01

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

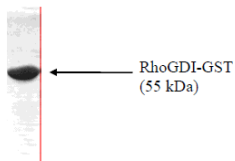
Material

The human Rho Guanine nucleotide Dissociation Inhibitor protein (RhoGDI) has been produced in a bacterial expression system. The protein contains a GST protein tag (~28 kDa) at its amino terminal. 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; 1.7 mM Tris pH 7.5, 0.3 mM MgCl₂, 0.3% sucrose and 0.1% dextran.

Purity

Purity is determined by scanning densitometry of proteins on SDS-PAGE gels. Samples are ~93% pure. The gel below shows 10 µg of RhoGDI-GST stained with coomassie blue.

RhoGDI-GST



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.