

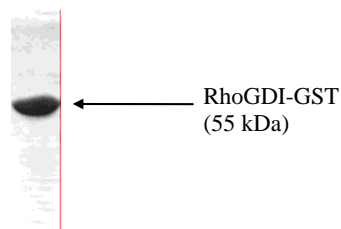
RhoGDI-GST Protein
(human recombinant)
Lot 115

Product Profile: GDI01

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.

Catalog Number	GDI01-A	1 x 25 µg
	GDI01-B	4 x 25 µg