

GST Protein

Recombinant protein expressed in E. coli cells

Catalog # G52-30U

Lot # F664-2

Product Description

Recombinant artificial sequences GST (1-226) was expressed in E. coli cells and purified on glutathione column. The gene accession number is [U13850](#).

Gene Aliases

None

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

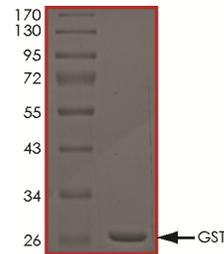
Scientific Background

GST or glutathione S-transferase family of enzymes comprises a long list of cytosolic, mitochondrial, and microsomal proteins that are used to create the so-called GST gene fusion system (1). These proteins are capable of multiple reactions with a multitude of substrates. GST tag has the size of 220 amino acids, which is relatively large compared to other tags like the myc- or the FLAG-tag that are quite small. GST can be used to purify and detect proteins of interest. In a GST gene fusion system, an expression vector consists of the gene sequence encoding the protein of interest and GST sequence alongside. After induction, the fusion proteins can be purified using affinity chromatography via its high affinity for glutathione (1).

References

1. Beckett, G.J. and Hayes, J.D. (1987). Glutathione S-transferase measurements and liver disease in man. J. Clin. Biochem. Nutr. 2, 1-24.

Purity



The purity of GST was determined to be **>95%** by densitometry. Approx. MW **25kDa**.

GST Protein

Recombinant protein expressed in E. coli cells

Catalog Number	G52-30U
Specific Lot Number	F664-2
Purity	>95%
Concentration	0.2 µg/µl
Stability	1yr At -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.