

Data Sheet

Glutathione S-Transferase (GST) Fusion Protein

S. japonicum, Recombinant

Catalog #:100343

Lot #: 190405 Conc.: 1.31 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
110 mM NaCl, 2.2 mM KCl and 20%
Glycerol.

Stability: At least 6 months at -80°C .
Avoid freeze/thaw cycles.

References:

1. Strange R.C., *et al.*, *Mutation Research*, **482(1-2)**: 21-26 (2001).
2. KunLiang G., *et al.*, *Analytical Biochemistry*, **192(2)**: 262-267 (1991).

Description: S. japonicum Glutathione S-Transferase also known as GST, GenBank Accession No. AAB59203, a.a. 1-218(end), expressed in an *E. coli* expression system, MW=28 kDa.

Applications: Glutathione S-transferase (GST) can be used as a control for GST-fusion tag proteins. It is found as a fusion tag in the commonly used pGEX vectors.

Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:

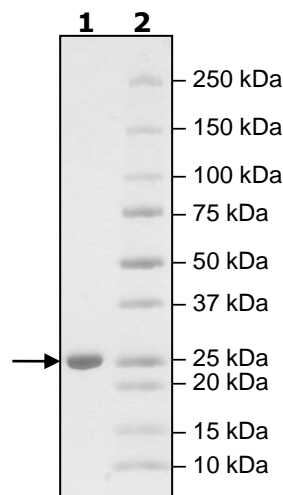
2 μg GST

Lane 2:

Protein Marker

MW: 28 kDa

Purity: $\geq 90\%$



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com