

Data Sheet

GYG/GYS2 Complex

Human, Recombinant

Catalog #: 100317**Lot #:** 190501**Conc.:** 0.15 mg/ml

Formulated in: 40 mM Tris-HCL, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% Glycerol and 250 mM Imidazole.

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

References:

1. Kamura, T., *et al.*, *Science*. 1999; **284(5414)**: 657-661.
2. Stebbins, C.E., *et al.*, *Science*. 1999; **284(5413)**: 455-461.

Description: GYG/GYS2 complex of human **GYG** (GenBank Accession No. NM_004130) a.a. 1-350(end) with N-terminal His-tag and MW= 42 kDa, and human **GYS2** (GenBank Accession No. NM_021957), a.a. 1-703(end) with C-terminal Strep-tag and MW= 82 kDa, individually expressed in an Sf9 infected baculovirus expression system.

Quality Assurance

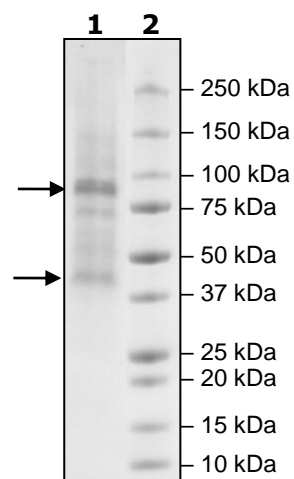
4-20% SDS-PAGE Coomassie staining

Lane 1:
3.6 μg GYG + GYS2

Lane 2:
Protein Marker

MW: GYG 42 kDa
GYS2 82 kDa

Purity: $\geq 69\%$



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