

Data Sheet

GLS1, His-tag

Human, Recombinant, N-terminal His-tag
Catalog #: 71102
Lot #: 200423-A **Conc.:** 0.20 mg/ml

Formulated in: 40 mM Tris-HCl pH 8.0,
 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-
 20, 20% glycerol.

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

References:

1. Van den Heuvel, A.P., *et al.*, *Cancer Biol. Ther.* 2012;**13(12)**:1185-1194.
2. Thangavelu, K., *et al.*, *Proc. Natl. Acad. Sci. USA* 2012;**109(2)**:7705-7710.

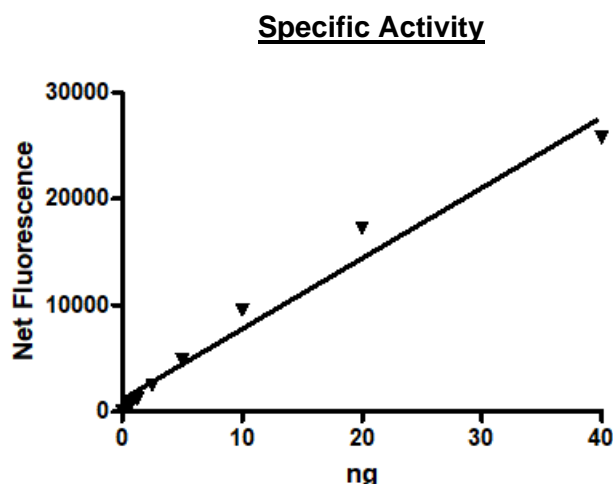
Description: Human glutaminase transcript variant 1 (GLS1), also known as GAC, GAM and KGA, GenBank Accession No. NM_014905, a.a. 17-669(end), with N-terminal His-tag, expressed in an *E. coli* cell expression system. MW=73 kDa.

Specific Activity: ≥25 nmol/min/μg

Assay conditions: Assay was performed according to the GLS1 Inhibitor Screening Assay Kit (BPS Bioscience, #79596) with GLS1 titrated from 40- 0 ng.

Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

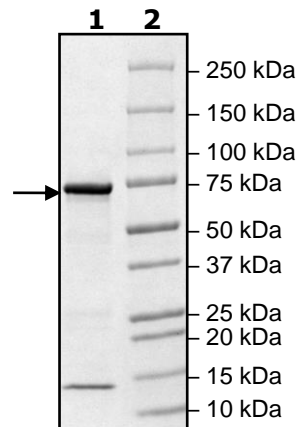


**4-20% SDS-PAGE
 Coomassie staining**

Lane 1:
 2 μg GLS1
Lane 2:
 Protein Marker

MW: 73 kDa

Purity: 80%



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