

Recombinant GLU protein

Catalog No: 81298, 81998

Expressed In: *E. coli*

Quantity: 50, 1000 µg

Concentration: 1.5 µg/µl

Source: Human

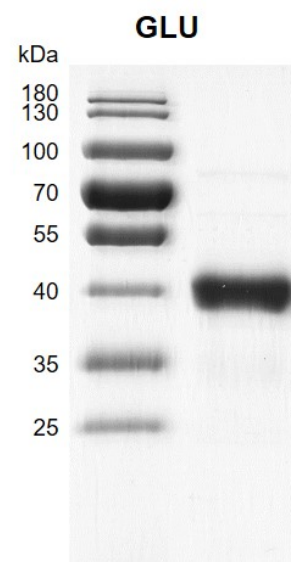
Buffer Contents: Recombinant GLU protein is supplied in 25 mM Tris pH 8.0, 300 mM NaCl, 10% glycerol, 0.5 mM TCEP.

Background: GLU (Glutamine synthetase), also called as GLU, is an enzymatic that catalyzes the synthesis of glutamine from glutamate and ammonia in an ATP-dependent reaction. This protein plays a role in ammonia and glutamate detoxification, acid-base homeostasis, cell signaling, and cell proliferation. Glutamine is an abundant amino acid, and is important to the biosynthesis of several amino acids, pyrimidines, and purines. Mutations in this gene are associated with congenital glutamine deficiency, and overexpression of this gene was observed in some primary liver cancer samples. There are six pseudogenes of this gene found on chromosomes 2, 5, 9, 11, and 12. Alternative splicing results in multiple transcript variants.

Protein Details: GLU protein was expressed in *E. coli* cells (accession number NP_001028216.1) with a C-terminal 6×His-Tag. The molecular weight of the protein is 43.1 kDa.

Application Notes: This product was manufactured as described in Protein Details. Where possible, Active Motif has developed functional or activity assays for recombinant proteins. Additional characterization such as enzyme kinetic activity assays, inhibitor screening or other biological activity assays may not have been performed for every product. All available data for a given product is shown on the lot-specific Technical Data Sheet.

Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of arrival.



Recombinant GLU protein gel
12.5% SDS-PAGE with Coomassie blue staining

MW: 43.1 kDa

Purity: >92%