

## Recombinant PKM2 protein

**Catalog No:** 81302, 81602

**Expressed In:** *E. coli*

**Quantity:** 50, 1000 µg

**Concentration:** 4 µg/µl

**Source:** Human

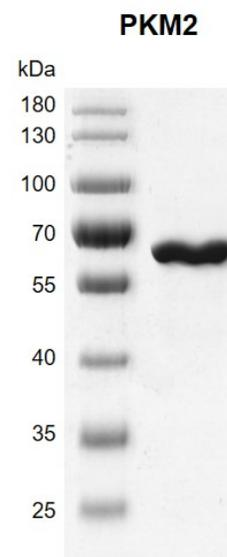
**Buffer Contents:** Recombinant PKM2 protein is supplied in 25 mM Tris-HCl pH 8.0, 300 mM NaCl, 10% glycerol and 0.5 mM TCEP.

**Background:** PKM (Pyruvate Kinase M2) is a pyruvate kinase that catalyzes the transfer of a phosphoryl group from phosphoenolpyruvate to ADP, generating ATP and pyruvate. This protein has been shown to interact with thyroid hormone and may mediate cellular metabolic effects induced by thyroid hormones. This protein has been found to bind Opa protein, a bacterial outer membrane protein involved in gonococcal adherence to and invasion of human cells, suggesting a role of this protein in bacterial pathogenesis. Several alternatively spliced transcript variants encoding a few distinct isoforms have been reported.

**Protein Details:** Full length human PKM2 protein was expressed in *E. coli* cells (accession number NP\_872270.1) with a C-terminal 6×His-Tag. The molecular weight of the protein is 60.1 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 PKM2 protein gel**  
10% SDS-PAGE with Coomassie blue staining

MW: 60.1 kDa

Purity: >95%