

PRODUCT INFORMATION



PDHK1 (human recombinant)

Item No. 42257

Overview and Properties

Synonyms: PDH Kinase 1, Pyruvate Dehydrogenase (acetyl-transferring) Kinase Isozyme 1, Pyruvate Dehydrogenase Kinase Isoform 1
Source: Active recombinant human PDHK1 purified from *E. coli*
Amino Acids: 30-436
Uniprot No.: Q15118
Molecular Weight: 46.3 kDa
Storage: -80°C (as supplied)
Stability: ≥1 year
Purity: ≥85% estimated by SDS-PAGE
Supplied in: 50 mM potassium phosphate, pH 7.5, with 250 mM potassium chloride, 800 mM lysine, 5% glycerol, and 20 mM DTT

Protein

Concentration: *batch specific* mg/ml

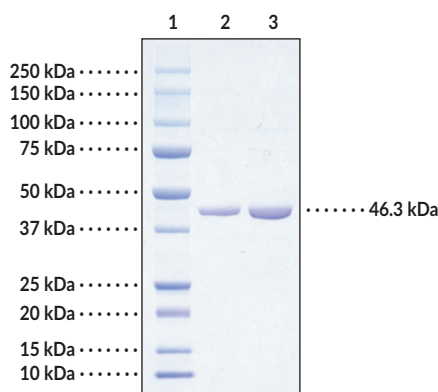
Activity: *batch specific* U/ml

Specific Activity: *batch specific* U/mg

Unit Definition: One unit is defined as the amount of enzyme required to produce 1 nmol of ADP per minute at 30°C in 40 mM Tris, pH 7.5, 50 mM MgCl₂, 1mM DTT, in the presence of 4 μM of PDHK1 substrate (peptide), and 1 mM ATP.

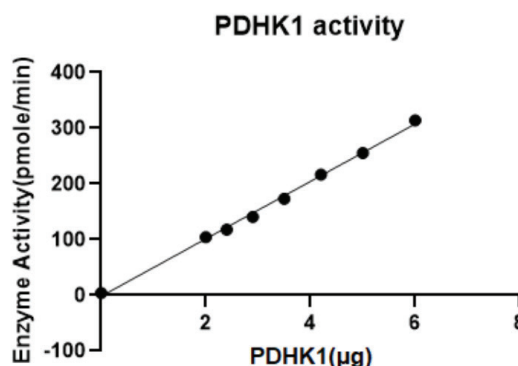
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Images



Lane 1: MW Markers
Lane 2: PDHK1 (2 μg)
Lane 3: PDHK1 (4 μg)

SDS-PAGE Analysis of PDHK1.



PDHK1 Activity Assay. PDHK1 activity was monitored by luminescence. In the presence of ATP, PDHK1 catalyzes the phosphorylation of a PDHK1 specific peptide substrate. The amount of ADP produced through the reaction was monitored using ADP-Glo Kinase Assay Kit (Promega).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 09/25/2025

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Pyruvate dehydrogenase kinase 1 (PDHK1) is a kinase and regulator of glycolysis and oxidative phosphorylation.¹⁻³ It is expressed primarily in the heart and to a lesser extent in skeletal muscle, liver, brain, placenta, lung, pancreas, and kidney and is localized to the mitochondria.¹ PDHK1 phosphorylates three serine residues on the E1 subunit of the pyruvate dehydrogenase complex, inhibiting the oxidative decarboxylation of pyruvate to acetyl-CoA and instead promoting its conversion to lactate.^{2,3} Pharmacologic inhibition of PDHK1 inhibits both proliferation and survival of T helper 17 (Th17) cells and suppresses Th17-mediated inflammation in rodent models of inflammatory bowel disease (IBD) and experimental autoimmune encephalomyelitis (EAE).² Levels of PDHK1 are increased in and associated with malignancy in lung cancers and pancreatic ductal adenocarcinomas, and pharmacologic inhibition of PDHK1 reverses the Warburg phenotype in various tumor models *in vitro* and *in vivo*.^{3,4} Cayman's PDHK1 (human, recombinant) protein can be used for activity assay, binding study, and Western blot (WB) applications.

References

1. Gudi, R., Bowker-Kinley, M.M., Kedishvili, N.Y., *et al.* Diversity of the pyruvate dehydrogenase kinase gene family in humans. *J. Biol. Chem.* **270**(48), 28989-28994 (1995).
2. Gerriets, V.A., Kishton, R.J., Nichols, A.G., *et al.* Metabolic programming and PDHK1 control CD4⁺ T cell subsets and inflammation. *J. Clin. Invest.* **125**(1), 194-207 (2015).
3. Xu, H., Ding, D., Han, X., *et al.* Discovery of ATP competitive PDHK1/2 dual inhibitors. *Bioorg. Med. Chem. Lett.* **122**, 130190 (2025).
4. Wu, P., Zhang, Z., Zhou, Y., *et al.* Novel dichloroacetophenone-based PDHK1 inhibitors as potent anticancer agents. *Drug Des. Devel. Ther.* **18**, 4661-4679 (2024).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM