

PRODUCT INFORMATION



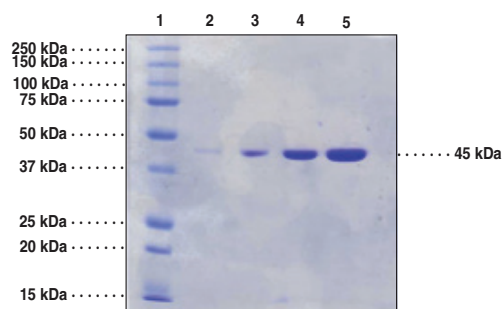
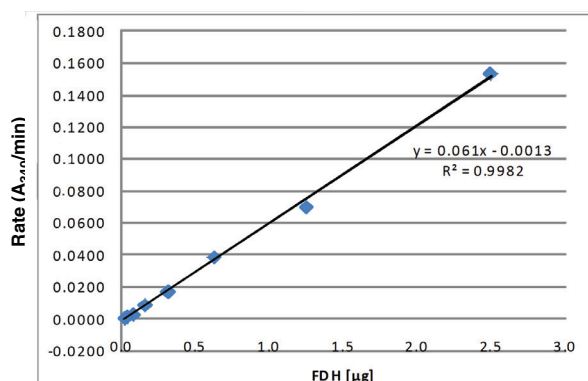
Formaldehyde Dehydrogenase (*P. putida* recombinant)

Item No. 10352

Overview and Properties

Synonym: FDH
Source: Recombinant N-terminal His-tagged protein expressed in *E. coli*
Amino Acids: S2-A399 (full-length)
Uniprot No.: P46154
Molecular Weight: 44.6 kDa
Storage: -80°C (as supplied)
Stability: ≥1 year
Purity: *batch specific* (≥95% estimated by SDS-PAGE)
Supplied in: *batch specific*
Protein
Concentration: *batch specific* mg/ml
Activity: *batch specific* U/ml. One unit is defined as the amount of enzyme required to produce 1 μmole of NADH per minute at 25°C in 50 mM potassium phosphate, pH 7.5, containing 200 μM formaldehyde.
Specific Activity: *batch specific* μmoles/min/mg

Images



Lane 1: MW Markers
Lane 2: Purified FDH (1 μg)
Lane 3: Purified FDH (2 μg)
Lane 4: Purified FDH (5 μg)
Lane 5: Purified FDH (10 μg)

Representative gel image shown; actual purity may vary between each batch but protein will be ≥95% pure.

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, 11/10/2021

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

Formaldehyde dehydrogenase (FDH) is a zinc-containing metalloenzyme that catalyzes the oxidation of formaldehyde to formate. *P. putida* FDH is an NAD-dependent enzyme that is a member of the class III alcohol dehydrogenase family.¹

Reference

1. Ito, K., Takahashi, M., Yoshimoto, T., *et al.* Cloning and high-level expression of the glutathione-independent formaldehyde dehydrogenase gene from *Pseudomonas putida*. *J. Bacteriol.* **176(9)**, 2483-2491 (1994).

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