

Data Sheet

FDH

P. putida, recombinant, N-terminal His tag
Catalog #: 71076
 Lot #: 121211 Conc.: 1.44 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
 110 mM NaCl, 2.2 mM KCl, 200 mM
 imidazole, and 20% glycerol.

Stability: >6 months at -80 °C

References:

1. Ito, K., *et al.*, *J Bacteriol.*
 1994;**176**:2483-2491.
2. Tanaka, K., *et al.*, *J Mol Biol.*
 2002;**324**:519-533.

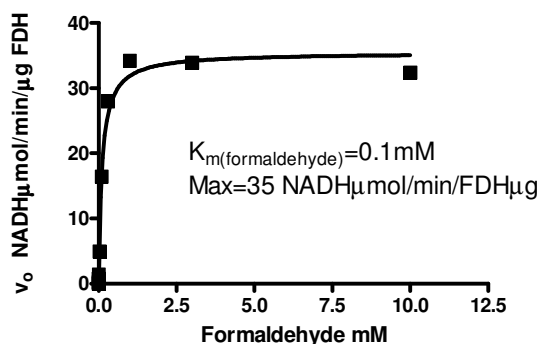
Description: *Pseudomonas putida*
 formaldehyde dehydrogenase (FDH)
 amino-acids 2 – 399 (end) (GenBank
 Accession No. D21201) with N-terminal
 His tag, MW = 42.9 kDa, expressed in
 an *E. coli* expression system.

Specific Activity: 35 $\mu\text{mol}/\text{min}/\mu\text{g}$
 Assay conditions: The reaction was
 monitored by measuring NADH
 production by following absorbance at
 340nm at room temperature for 5min.
 Buffer: pH 7.3, 20 mM Tris, 50 mM
 NaCl, 20 μM Fe^{+2} , 500 mM α -
 ketoglutarate, 2 mM ascorbic acid,
 0.01% Tween-20, and 1 mM NAD+.

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

Quality Assurance

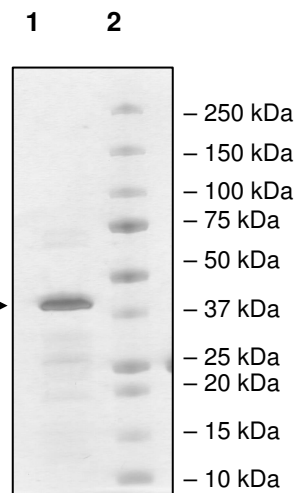
**Kinetics of Formaldehyde Dehydrogenase
 for Formaldehyde**



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

Lane 1:
 2.88 μg FDH
Lane 2:
 Protein Marker

MW: 42.9 kDa →
Purity: ≥74%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Sequence:

MHHHHHHHSGNRGVVYLGSGKVEVQKIDYPKMQDPRGKKIEHGVILKVVSTNICGSDQ
HMVRGRRTTAQVGLVLGHEITGEVIEKGRDVENLQIGDLVSVPFNVACGRRCRSCKEMH
TGVCLTVNPARAGGAYGYVDMGDWTGGQAEYLLVPYADFNLLKLPDRDKAMEKIRD
LTCLSDILPTGYHGAVTAGVGPSTVYVAGAGPVGLAAAASARLLGAAVVIVGDLNP
ARLAHAKAQGF EIADLSLDTPLHEQIAALLGEPEVDCAVDAVGFEARGHGHEGAKHE
APATVLNSLMQVTRVAGKIGIPGLYVTEDPGAVDAAKIGSLSIRFGLGWAKSHSFHT
GQTPVMKYNRALMQAIMWDRINIAEVVGVQVISLDDAPRGYGEFDAGVPPKFFVIDPHK
TFSAA

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com