

## Data Sheet

### IDH1

Human recombinant, with C-terminal FLAG-tag

**Catalog #:** 71075

**Lot#:** 120131      **Conc.:** 0.55 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0,  
110 mM NaCl, 2.2 mM KCl, 80 ng/μl Flag  
peptide, 20% glycerol and 3 mM DTT.

**Stability:** >6 months at -80°C.

#### References:

1. Dang, L., *et al.* (2010). *Trends Mol. Med.* **16(9)**:387-97.
2. Paschka, P., *et al.* (2010). *J. Clin. Oncol.* **28(22)**:3636-43
3. Fu, Y., *et al.* (2010). *Biochem. Biophys. Res. Commun.* **397(2)**:127-30.

#### Description:

Human Isocitrate Dehydrogenase (IDH1), also known as oxalosuccinate decarboxylase, GenBank Accession No. NM\_005896, full length [a.a. 1-414(end)] with C-terminal FLAG-tag, MW= 47.6 kDa, expressed in Sf9 cells via a baculovirus expression system.

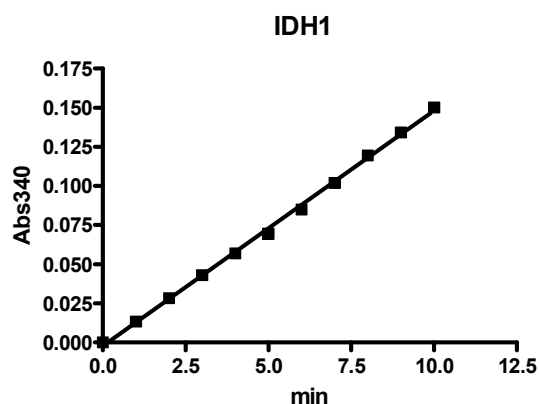
**Specific Activity:** 7000 pmol/min/μg.

**Assay conditions:** Reaction was conducted in 50 mM Tris (pH7.4), 10 mM MgCl<sub>2</sub>, 200 mM DL-isocitrate, and 200 mM NADP. Incubate for 30 minutes at room temperature. Read at Abs340 using Tecan Infinite 1000 (5 flash/10 ms settle time).

#### Application:

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

### Quality Assurance



#### 10% SDS-PAGE Coomassie staining

**Lane 1:**

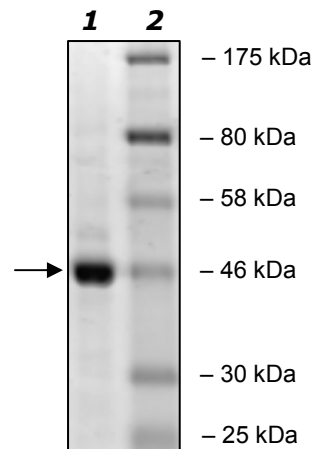
3 μg IDH1

**Lane 2:**

Protein Marker  
BioLabs (#P7708L)

**MW:** 48 kDa

**Purity:** ≥94%



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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)