

## Data Sheet

### IDH2

Human, recombinant, C-terminal Flag-tag

**Catalog #:** 71074

Lot#: 110106-2      Conc.: 0.07 mg/ml

**Formulated in:** 45 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 mM KCl, 90 µg/ml FLAG peptide, 10% glycerol, and 3 mM DTT.

**Stability:** >6 months at -80 °C

#### References:

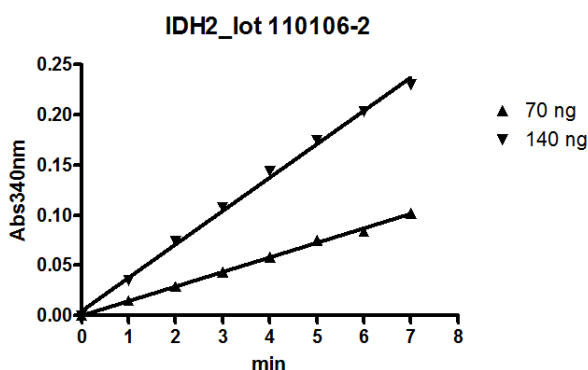
1. Schaap, F.G., *et al.*, *Adv Anat Pathol* 2013;**20(1)**:32-38.
2. Grassian, A.R., *et al.*, *J Biol Chem* 2012;**287(50)**:42180-42194.

**Description:** Human isocitrate dehydrogenase 2, also known as IDH2, (GenBank Accession No. NM\_002168), a.a. 1-452(end) with C-terminal Flag-tag, MW = 52 kDa, expressed in a baculovirus infected Sf9 cell expression system.

**Specific Activity:** 14,200 pmol/min/µg.  
 Assay conditions: The reaction mixture contained 50 mM Tris, pH 7.4, 5 mM MgCl<sub>2</sub>, 200 µM DL-isocitrate, 200 µM NADP<sup>+</sup>, and various amount of IDH2 in a volume of 200 µL. Reactions were initiated with the addition of IDH2, and the increase in NADPH absorbance at 340 nm was continuously monitored for 7 min. Molar extinction coefficient of NADPH is 6,220 M<sup>-1</sup>cm<sup>-1</sup>.

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

## Quality Assurance

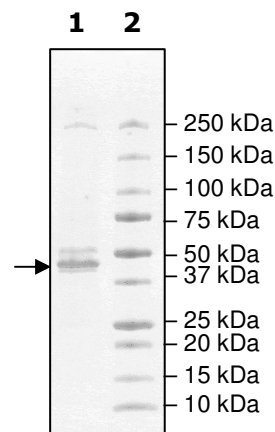


### 10% SDS-PAGE Coomassie staining

**Lane 1:**  
1.4 µg IDH2

**Lane 2:**  
Protein Marker

**Purity:** ≥59%  
**MW:** 52 kDa



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)



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

**Sequence:**

MAGYLRVVRSLCRASGSRPAWAPAALTAPTSQEQPRRHYADKRIKVAKPVVEMDGDE  
MTRIIWQFIKEKLILPHVDIQLKYFDLGLPNRDQTDDQVTIDSALATQKYSVAVKCATITP  
DEARVEEFKLLKMWKSPNGTIRNILGGTVFREPIICKNIPRLVPGWTKPITIGRHAHGDQ  
YKATDFVADRAGTFKMFVTPKDGSGVKEWEVYNFPAGGVGMGMYNNTDESISGFAHSC  
FQYAIQKKWPLYMSTKNTILKAYDGRFKDIFQEIFDKHYKTDFDKNKIWEHRLIDDMVA  
QVLKSSGGFVWACKNYDGDVQSDILAQQFGSLGLMTSVLVCPDGKTI EAEAAHGTVTR  
HYREHQGRPTSTNPIASIFAWTRGLEHRGKLDGNQDLIRFAQMLEKVCVETVESGAM  
TKDLAGCIHGLSNVKLNEHFLNTTDFLDTIKSNLDRALGRQDYKDDDDK

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)