

## Data Sheet

### IDH2 (R140Q)

Human, recombinant, C-terminal FLAG-tag

**Catalog #:** 71100

**Lot#:** 130718

**Conc.:** 0.12 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 80 µg/ml FLAG peptide, 0.04% Tween-20, and 20% glycerol.

**Stability:** >6 months at -70°C. Avoid freeze/thaw cycles.

#### 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.

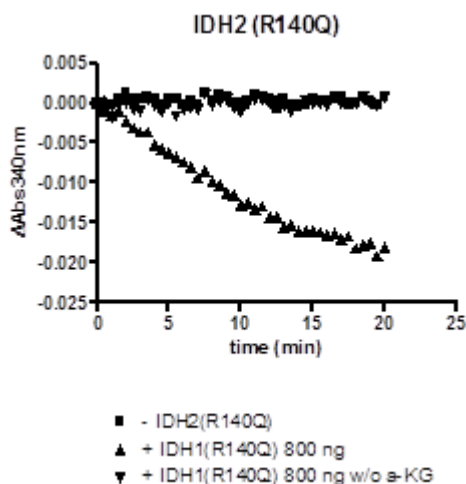
**Description:** Human isocitrate dehydrogenase 2, also known as IDH2, GenBank Accession No. NM\_002168, a.a. 1-452(end) with with Arg to Gln mutation on a.a. 140 and C-terminal FLAG-tag, MW = 52 kDa, expressed in a baculovirus infected Sf9 cell expression system. Note: This protein runs aberrantly by SDS-PAGE.

**Specific Activity:** 61 pmol/min/µg.

Assay conditions: IDH reductive activity was measured in 200 µl reaction containing 25 mM Tris (pH7.4), 150 mM NaCl, 10 mM MgCl<sub>2</sub>, 0.03% BSA, 1 mM α-Ketoglutarate, 10 µM NADPH and IDH. Depletion of NADPH was monitored continuously at Abs340 nm for 20 min. Molar extinction coefficient of NADPH is 6,200 M<sup>-1</sup>cm<sup>-1</sup>.

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

## Quality Assurance

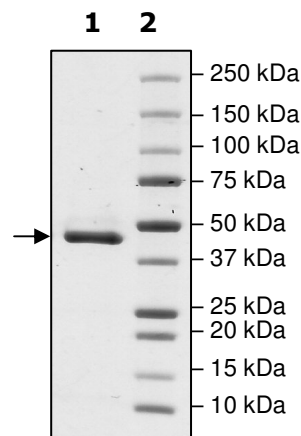


### 4-20% SDS-PAGE Coomassie staining

**Lane 1:**  
2.4 µg IDH2 (R140Q)

**Lane 2:**  
Protein Marker

**Purity:** ≥90%  
**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:**

MAGYLRVVRS�CRASGSRPAWAPAALTAPTSQEPPRRHYADKRIKVAKPVVEMDGDE  
MTRIIWQFIKEKLILPHVDIQLKYFDLGLPNRDQTDQVTIDSALATQKYSVAVKCATITP  
DEARVEEFKLLKMWKSPNGT**IQ**NILGGTVFREPIICKNIPRLVPGWTKPITIGRHAHGQ  
YKATDFVADRAGTFKMVFTPKDGSGVKEWEVYNFPAGGVGMGMYNTDESISGFAHSC  
FQYAIQKKWPLYMSTKNILKAYDGRFKDIFQEIFDKHYKTDFDKNKIWYEHRLIDDMVA  
QVLKSSGGFVWACKNYDGDVQSDILAQGFSLGLMTSVLVCPDGKTIEAAAHGTVTR  
HYREHQGRPTSTNPIASIFAWTRGLEHRGKLDGNQDLIRFAQMLEKVCVETVESGAM  
TKDLAGCIHGLSNVKLNEHFLNTTDFLDTIKSNLDRALGRQ**DYKDDDDK**

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)