

Data Sheet

AHCY

Recombinant, N-terminal His tag

Catalog #: 70017**Lot#:** 120117 **Conc.:** 4.0 mg/ml

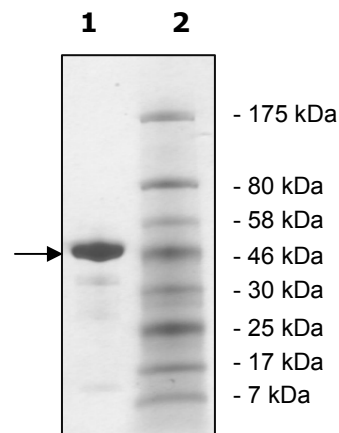
Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 200 mM imidazole, and 20% glycerol.**Stability:** >6 months at -80°C**References:**

1. Leal, J.F. *et al.*, *Carcinogenesis* **29** (11), 2089-2095 (2008).
2. Hermes, M. *et al.*, *Cell. Physiol. Biochem.* **22** (1-4), 223-236 (2008).

Description:Recombinant *S. solfataricus* S-adenosyl homocysteine hydrolase, or AHCY, (GenBank Accession No. Z50174) amino acids 2 – 417 (end) with N-terminal His tag, MW= 47 kDa, expressed in an *E. coli* expression system.**Application:**

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

Quality Assurance

**4-20% SDS-PAGE
Coomassie staining****Lane 1:**
2 µg AHCY**Lane 2:**
Protein Marker
BioLabs (#P7708L)**MW:** 47 kDa**Purity:** ≥86%

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.comPlease visit our website at: www.bpsbioscience.com



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

Protein Sequence:

MHHHHHHSYKIKDLSLASEGKKQIEWAERHMPTLMEIRKRFKAEKPLKGINISAVLHV
TKETAALVKTLKIGGANVALAGSNPLSTQDDVAAALVEEGISVFAWKGENETEYYSNI
ESIVKIHEPNIVMDDGADLHAYIHEKVSSKLDIYGGTEETTTGVIRLKAMEKDGVLKYPL
VAVNNAYTKYLFDNRYGTGQSAIDGILRATNILIAGKIAVVAGYGWVGRGIANRLRGM
GARVIVTEVDPIRALEAVMDGFDVMPIAEASKVGDIFVTATGNTKAIRVEHMLNMKDG
AILS NAGHFNVEVDVKGLKETAVKVRNIRPYVDEYTL PNGKRVYLLADGRLVNLAAAE
GHPSEVMDMSFANQALAVEYLVKNRGKLEKKVYNMPMELDYEVARIK LKSMGIQIDE
LTEEQKEYLEQWKS GT

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