

Data Sheet

AHCY

Human, recombinant, N-terminal GST tag

Catalog #: 50260

Lot#: 110503DS **Conc.:** 1.1 mg/ml

Formulated in: 45 mM Tris-HCl, pH 8.0,
124 mM NaCl, 2.4 mM KCl, 10% 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:

Human AHCY (human S-adenosyl homocysteine hydrolase) (GenBank Accession No. NM_000687), (amino acids 2-end) with N-terminal GST tag, MW = 73 kDa, expressed in an *E. coli* expression system.

Application:

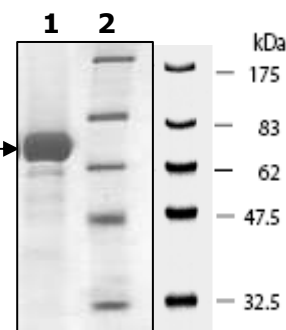
Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

10% SDS-PAGE Coomassie staining

Lane 1:
7 µg AHCY →

Lane 2:
Protein Marker
BioLabs (#P7708L)



MW: 73 kDa
Purity: >90%

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