

Data Sheet

SUV4-20H1

Human, recombinant, N-terminal GST tag

Catalog #: 51090

Lot#: 130403-10 **Conc.:** 0.31 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
110 mM NaCl, 2.2 mM KCl, 20% Glycerol,
and 3 mM DTT.

Stability: >6 months at -80 °C

References:

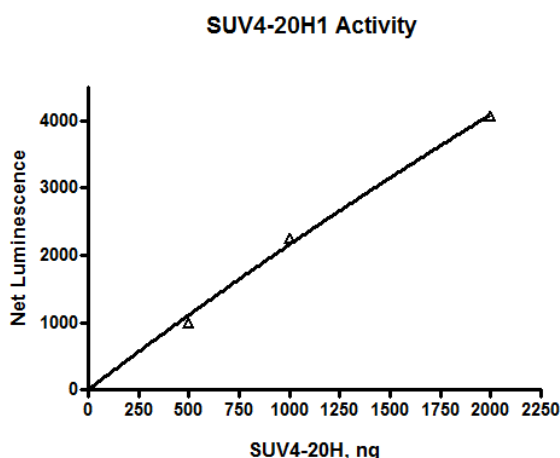
1. Yang H. *et al.*, *J Biol Chem.* 2008 May 2;**283(18)**:12085-92.
2. Benetti R *et al.*, *J Cell Biol.* 2007 Sep 10;**178(6)**:925-36.

Description: Human SUV4-20H1 (GenBank Accession No. NM_017635), amino acids 2-387, with N-terminal GST tag, MW= 70 kDa, expressed in *E. coli* expression system.

Specific Activity: 0.0001 pmol/min/μg
Assay Conditions: 50 μl reaction mix (50 mM TRIS pH 8.8, 5mM MgCl₂, 4mM DTT, 0.05% Tween-20, 800 μM S-adenosylmethionine, and SUV4-20H1 add to the wells coated with the substrate. Incubate for 4 hr. Add primary antibody 18, incubate 1 hr. Then, add secondary HRP-labeled antibody and incubate 30 min. Finally, add HRP chemiluminescent substrates and read luminescence.

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

Quality Assurance

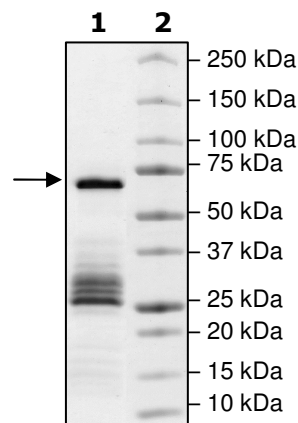


4-20% SDS-PAGE Coomassie staining

Lane 1:
2 μg SUV4-20H1

Lane 2:
Protein Marker

MW: 70 kDa
Purity: ≥31%



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



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

Sequence:

MSPILGYWKIKGLVQPTRLLLEYLEEKYEEHLYERDEGDKWRNKKFELGLEFPNLPYYI
DGDVKLTQSMARIYIADKHNMLGGCPKERAIEISMLEGAVLDIRYGVSRAYSKDFETLK
VDFLSKLPEMLKMFEDRLCHKTYLNGDHVTHPDFMLYDALDVVLYMDPMCLDAFPKLV
CFKKRIEAIQIDKYLKSSKYIAWPLQGQWQATFGGGDHPPKSDLVPRGSPEKWLGESK
NMVVNGRRNGGKLSNDHQQNQSKLQHTGKDTLKAGKNAVERRSNRCNGNSGFEG
QSRYPVSSGMSAKELCENDDLATSLVLDPYLGFQTHKMNTSAFPSRSSRHFSKSDSF
SHNPNVRFRIKGRQEELKEVIERFKKDEHLEKAFKCLTSGEWARHYFLNKNKMQEK
LFKEHVFIYLRMFATDSGFEILPCNRYSEQNGAKIVATKEWKRNDKIELLVGCIAELS
EIEENMLLRHGENDFSVMYSTRKNCAQLWLGPAAFINHDCRPNCKFVSTGRDTACVK
ALRDIEPGEEISCYYGDGFFGENNEFCCECYTCERRGTGAFKSRVGLPAPAPVINSKYG
LRETDKRLNRLKKGDSKNSDSQSVSSNTDADTTQE

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