

## Data Sheet

### SUV4-20h2

Human, recombinant, N-terminal GST tag

**Catalog #:** 51060

Lot#: 80828

Conc.: 0.66 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0,  
110 mM NaCl, 2.2 mM KCl, 0.05% tween-  
20, 3 mM DTT, 16 mM Glutathione and  
20% Glycerol**Stability:** >6 months at -80 °C**References:**

1. Tryndyak, V.P et al., Cancer Biol. Ther. **5** (1), 65-70 (2006)
2. Wan, D et al., Proc. Natl. Acad. Sci. U.S.A. **101** (44), 15724-15729 (2004)

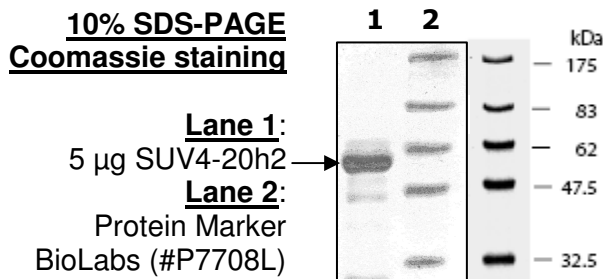
**Description:**

Human SUV4-20h2 (GenBank Accession No. NM\_032701), amino acids 2-280, with N-terminal GST tag, MW= 58 kDa, expressed in an E.coli expression system.

**Application:**

Useful for Western blot, ELISA, and SDS-PAGE analysis, antibody production, and for protein arrays.

### Quality Assurance

**MW:** 58 kDa**Purity:** ≥50%

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)