

## Recombinant SMYD4 protein

**Catalog No:** 31408, 31808

**Expressed In:** Baculovirus

**Quantity:** 20, 1000 µg

**Concentration:** 0.6 µg/µl

**Source:** Human

**Buffer Contents:** Recombinant SMYD4 protein is supplied at a concentration of 0.6 µg/µl in 25 mM HEPES pH 7.5, 300 mM NaCl, 5% glycerol, 0.04% Triton X-100, 0.2 mM TCEP.

**Background:** SET and MYND domain-containing protein 4 (SMYD4) is a member of the SMYD family of histone methyltransferases that may function as a potential tumor suppressor in breast cancer cells.

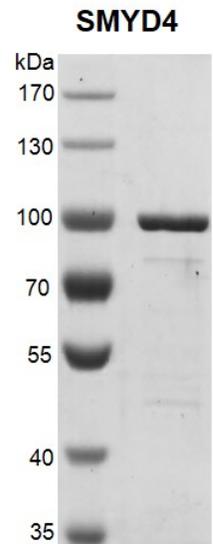
**Protein Details:** Recombinant SMYD4 (accession number NP\_443160.2) was expressed in Sf9 cells and contains an N-terminal FLAG tag with an observed molecular weight of 90.9 kDa. The recombinant protein is >90% pure by SDS-PAGE.

**Application Notes:** Recombinant SMYD4 is suitable for use in the study of enzyme kinetics, inhibitor screening, and selectivity profiling.

**Specific Activity:** Not published.

**Storage and Guarantee:** Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



**Recombinant SMYD4 protein gel.**  
Recombinant SMYD4 protein was run on an 8% SDS-PAGE gel and stained with Coomassie Blue.