

## Recombinant SETD2 (1392-2564) protein

**Catalog No:** 31399, 31799

**Expressed In:** Baculovirus

**Quantity:** 20, 1000 µg

**Concentration:** 0.2 µg/µl

**Source:** Human

**Buffer Contents:** Recombinant SETD2 (1392-2564) protein is supplied at a concentration of 0.2 µg/µl in 25 mM HEPES- NaOH pH 7.5, 300 mM NaCl, 5% glycerol, 0.04% Triton X-100, 0.2 mM TCEP.

**Background:** SETD2 (**SET Domain Containing 2**) is a histone methyltransferase that methylates histone H3 at lysine 36. H3K36 methylation is associated with active chromatin. SETD2 also contains a novel transcriptional activation domain and has been found associated with hyperphosphorylated RNA pol II.

**Recombinant SETD2 (1392-2564)** includes amino acids 1392-2564 that contains the SET domain of SETD2 that catalyzes methylation of lysine residues.

**Protein Details:** Recombinant SETD2 (1392-2564) (accession number NP\_054878.5) was expressed in Sf9 cells and contains an N-terminal FLAG tag with an observed molecular weight of 134 kDa. The recombinant protein is >70% pure by SDS-PAGE.

**Application Notes:** Recombinant SETD2 (1392-2564) is suitable for use in the study of enzyme kinetics, inhibitor screening, and selectivity profiling.

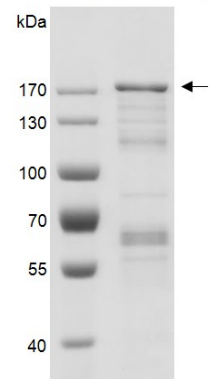
**Specific Activity:** H3K36 methyltransferase.

**HMT Assay Conditions:** 2 µg recombinant nucleosomes were incubated with SETD2 (1392-2564) in reaction buffer for 3 hr at RT. Western blot was used to detect reaction products (H3K36me2 Antibody, Cat# 39891). SETD2 (1392-2564) alone was used as negative control.

**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.

**SETD2 (1392-2564)**



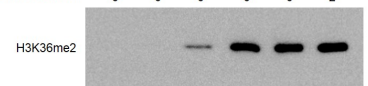
**Recombinant SETD2 (1392-2564) protein gel.**

SETD2 (1392-2564) protein was run on an 8% SDS-PAGE gel and stained with Coomassie Blue.

MW: 134 kDa.

Purity: > 70%

SETD2-SET	1	0	0.25	0.5	1	1	µg
Nucleosomes	0	2	2	2	2	0	
Histone Octamers	0	0	0	0	0	2	



**HMT Assay for Recombinant SETD2 (1392-2564) protein activity.**

2 µg recombinant nucleosomes were incubated with SETD2 (1392-2564) in reaction buffer for 3 hr at RT. Western blot was used to detect reaction products (H3K36me2 Antibody, Cat# 39891). SETD2 (1392-2564) alone was used as negative control.