

# SET07 monoclonal antibody

## (43AT551.86.76)

SET07 is a histone methyltransferase that methylates Lys20 of histone H4, a specific tag for epigenetic transcriptional repression. The nuclear SET07 protein, which associates with silent chromatin on euchromatic arms but shows no association with constitutive heterochromatin, prefers nucleosomes as substrate compared to free histones. SET07 may play a role in maintaining silent chromatin by preventing neighboring acetylation of the H4 tail.

This antibody is covered by our [Worry-Free Guarantee](#).

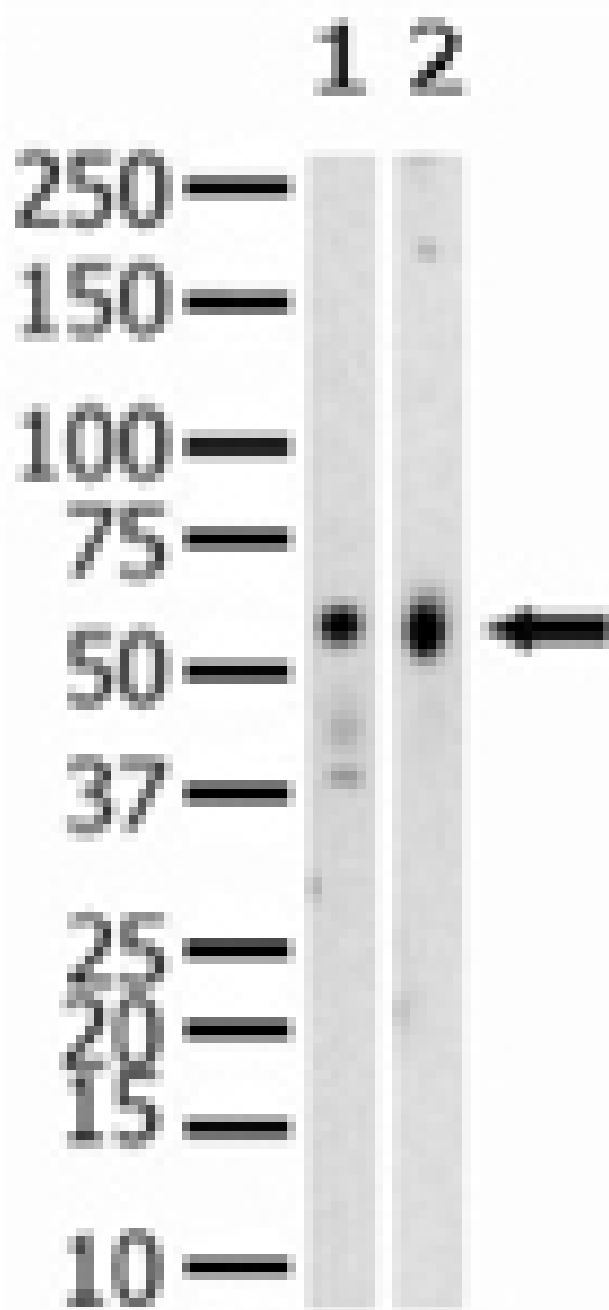
### Ordering Information

[Order Online »](#)

ADI-905-694-100	100µg
-----------------	-------

### Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of SET07: Lane 1 – SAOS-2 Lane 2 – NIH/3T3 cell lysates.

## Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Application ELISA, WB

Clone 43AT551.86.76

Formulation Liquid. In PBS containing 0.09% sodium azide.

Host Mouse

Immunogen Recombinant human SET07 (N-terminal fragment).

Isotype IgG1

Purity Detail Peptide affinity purified.

**Recommendation** ELISA (1:1,000)Western Blot (1:100-1:500)Suggested dilutions/conditions may not be  
**Dilutions/Conditions** available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Human, Mouse

UniProt ID Q8WTS6

**Worry-free Guarantee** This antibody is covered by our [Worry-Free Guarantee](#).



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)