

SET7/9 (human), (recombinant)

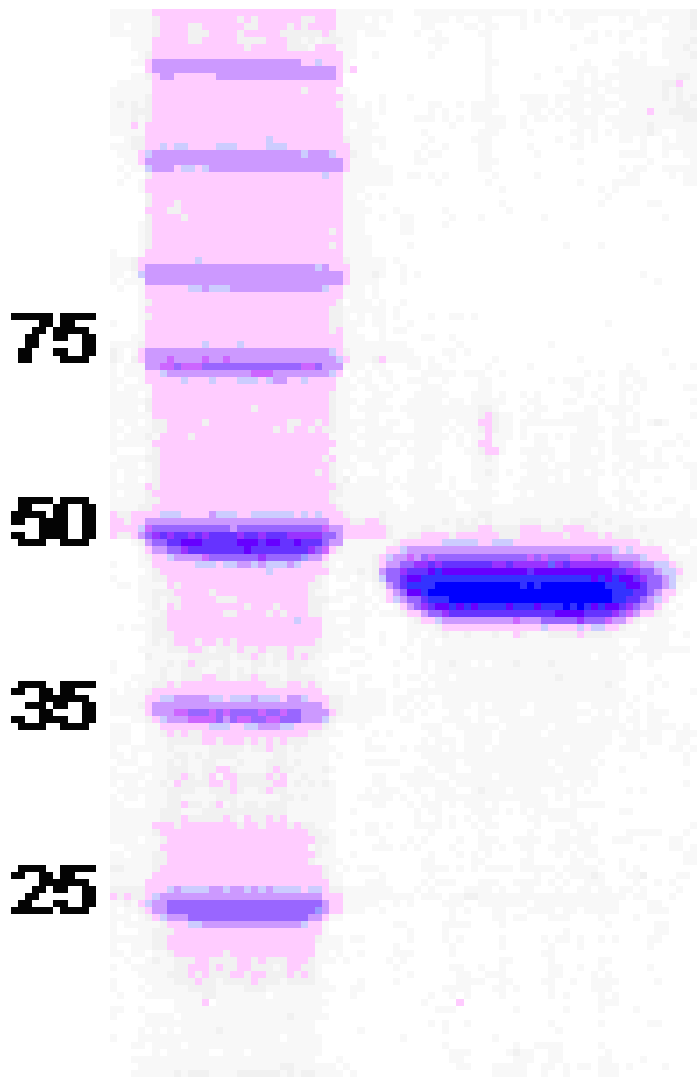
Histone methyltransferase that transfers methyl groups to Lys⁴ of histone H3, in complex with S-adenosyl-L-methionine.

Citations: 6

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



10% SDS-PAGE of human SET7/9 (Prod. No. ALX-201-178).

Handling & Storage

Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Histone H3-K4 methyltransferase
Concentration	1mg/ml
Endotoxin Content	<1.0EU/μg protein (LAL test)
Formulation	Liquid. In 50mM TRIS-HCl, pH 7.5, containing 0.2M sodium chloride, 5mM DTT and 20% glycerol.
MW	~40.7kDa.
Purity	≥95% (SDS-PAGE)
Source	Produced in <i>E. coli</i> .
UniProt ID	Q8WTS6

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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