

Dnmt3a (mouse) monoclonal antibody (64B1446)

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

Citations: 11

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-370-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

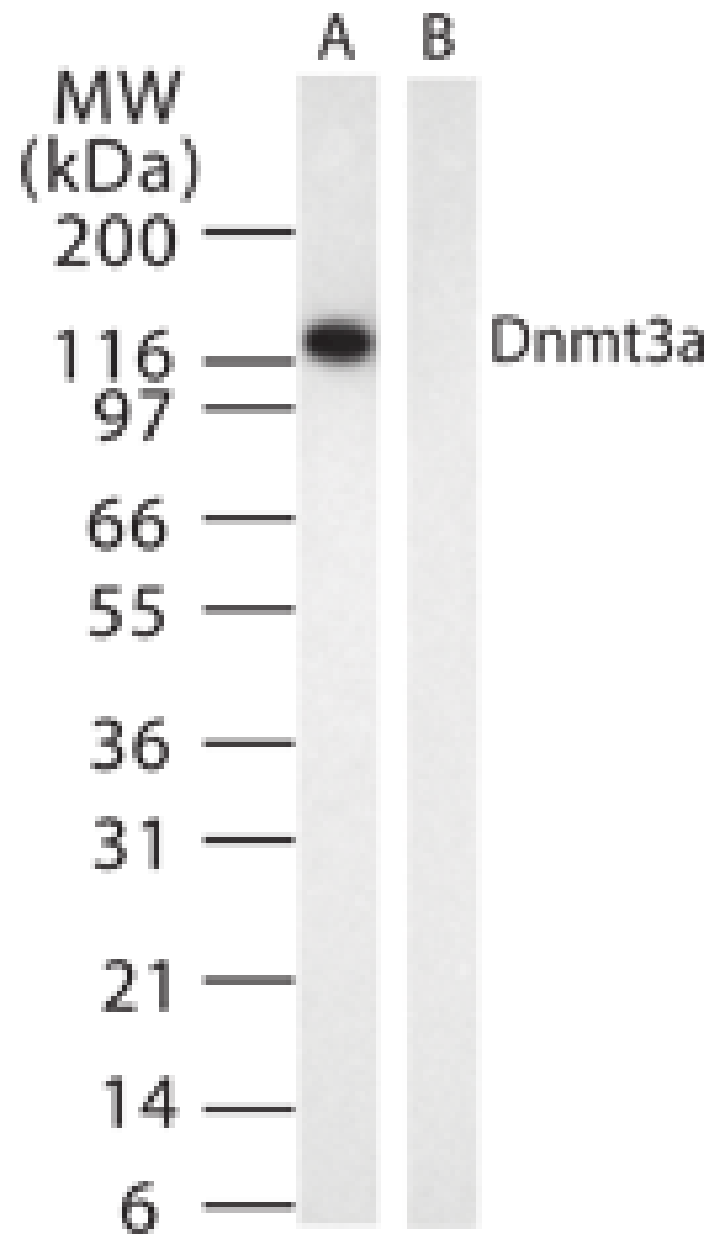
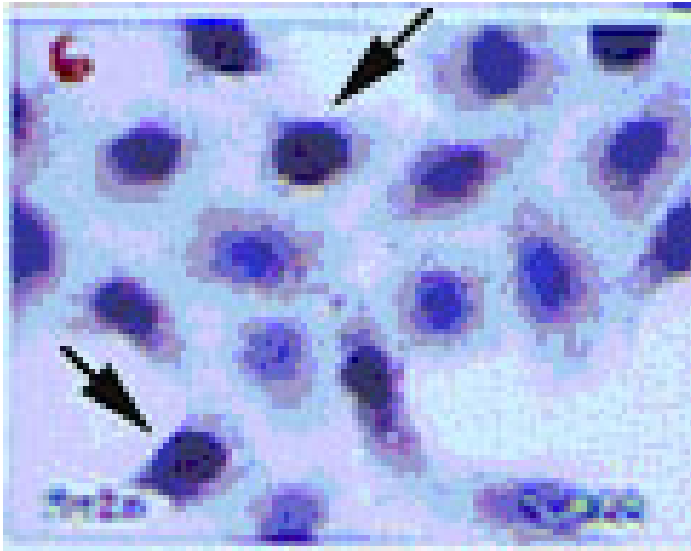


Figure: Western blot analysis of Dnmt3a using ALX-804-370 at 2 μ g/ml against 10 μ g of 293 cell lysate transfected with either mouse Dnmt3a (lane A) or mouse Dnmt3b (lane B).

Handling & Storage

Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	DNA methyltransferase-3a
Application	ChIP, ICC, IHC, IP, WB
Clone	64B1446
Crossreactivity	Weakly cross-reacts with human Dnmt3a.
Formulation	Liquid. In PBS containing 0.05% sodium azide
Host	Mouse
Immunogen	Recombinant mouse Dnmt3a (DNA methyltransferase-3a).
Isotype	IgG1
Purity Detail	Protein G-affinity purified.
Recommendation	Immunocytochemistry (5µg/ml)Immunohistochemistry (1:400)Western Blot
Dilutions/Conditions	(2µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Mouse
Specificity	Recognizes C-terminal epitope aa 705-908 of mouse Dnmt3a.
UniProt ID	O88508



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

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

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

France
Phone: +33 1 42 44 55 55
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

This antibody is covered by our **Worry Free Guarantee**