

# Dnmt3a (mouse) monoclonal antibody (64B1446)

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

Citations: 11

[View Online »](#)

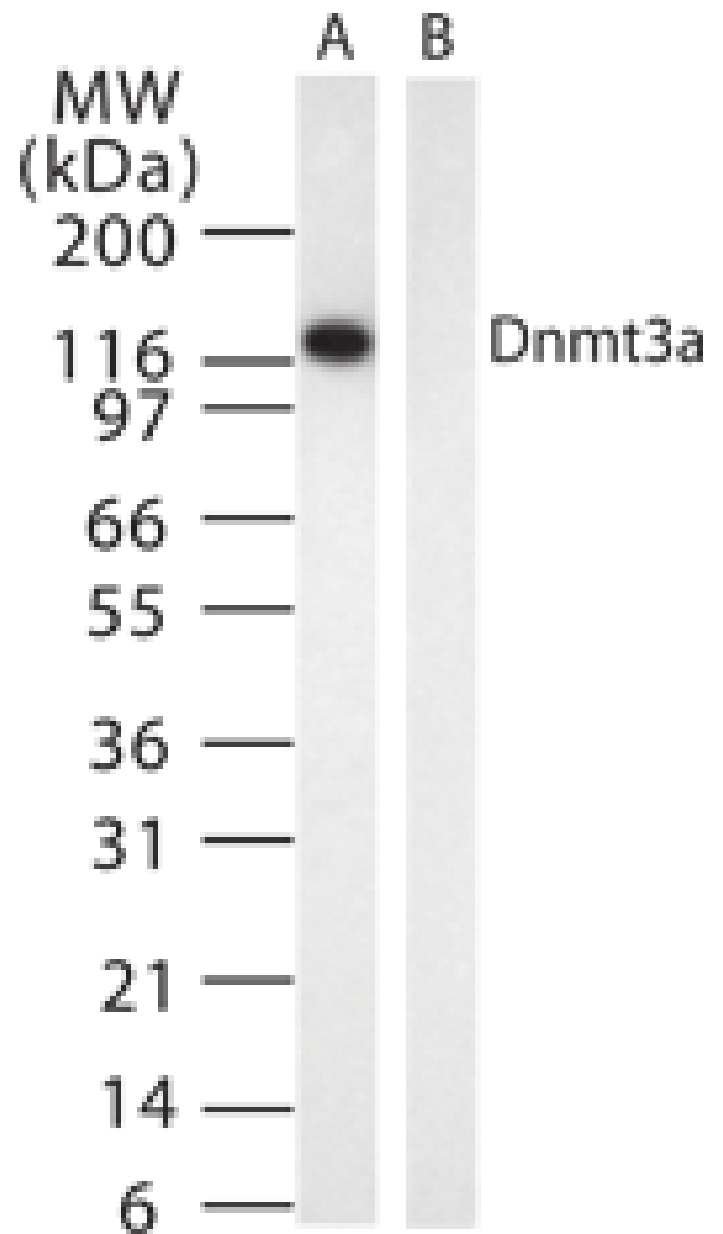
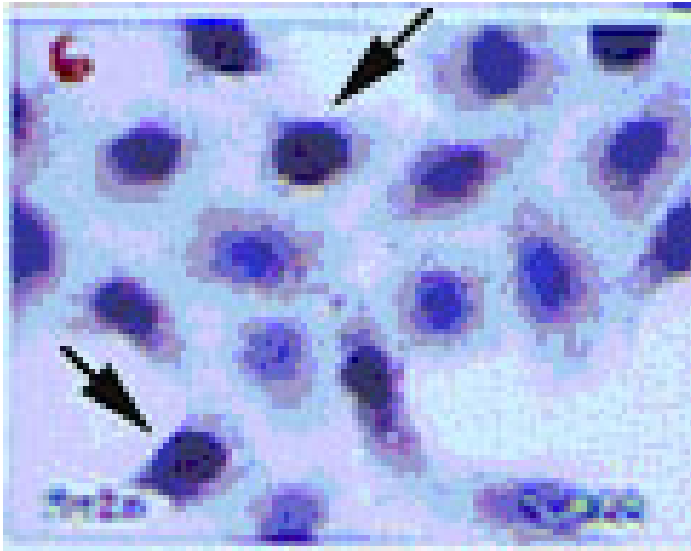
## Ordering Information

[Order Online »](#)

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

Manuals, SDS & CofA

[View Online »](#)



**Figure:** Western blot analysis of Dnmt3a using ALX-804-370 at 2 $\mu$ g/ml against 10 $\mu$ 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 Dilutions/Conditions	Immunocytochemistry (5µg/ml)Immunohistochemistry (1:400)Western Blot (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

Worry-free Guarantee

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

Last modified: May 29, 2024



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)