

Dnmt1 monoclonal antibody (60B1220.1)

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

Citations: 20

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-3690	100µg
--------------	-------

Manuals, SDS & CofA

[View Online »](#)

MW (kDa)

205-

120-

84-

52-

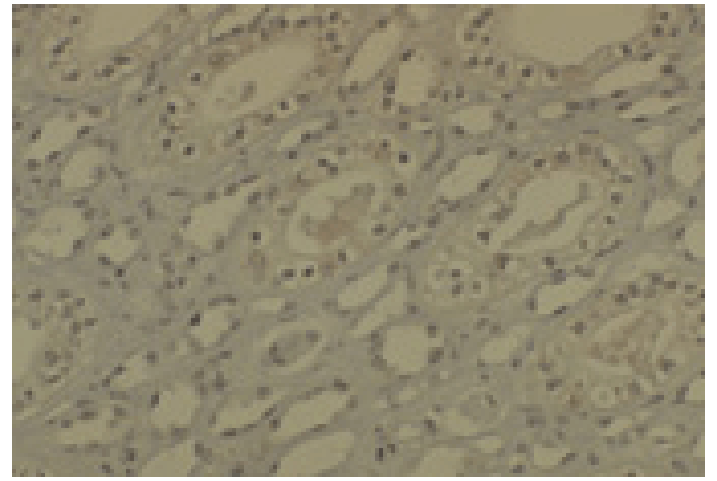


Figure 2: Immunohistochemistry of Dnmt1 using ALX-804-369 at 2 μ g/ml on formalin-fixed paraffin-embedded medullar kidney tissue sections.

Handling & Storage

Use/Stability	Stable for 6 months when stored at +4°C.
Handling	Avoid freeze/thaw cycles.
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-1
Application	ChIP, ICC, IHC (PS), IP, WB
Clone	60B1220.1
Formulation	Liquid. In PBS containing 0.05% BSA and 0.05% sodium azide.
Host	Mouse
Immunogen	Synthetic peptide corresponding to aa 637-650 of human Dnmt1 (DNA methyltransferase-1).
Isotype	IgG1
Purity Detail	Protein G-affinity purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (paraffin sections, 1-2µg/ml)Western Blot (0.1-0.5µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse, Zebrafish
Technical Info / Product Notes	Replacement for ADI-KAM-TF040

UniProt ID

P26358

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)