

# Daxx (human) monoclonal antibody (DAXX-01)

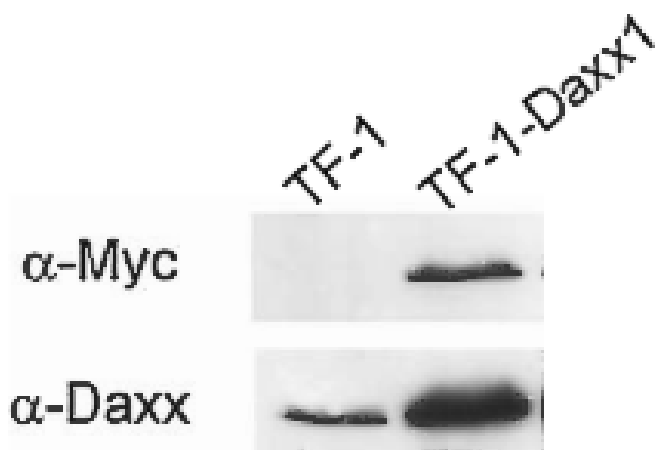
This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 2

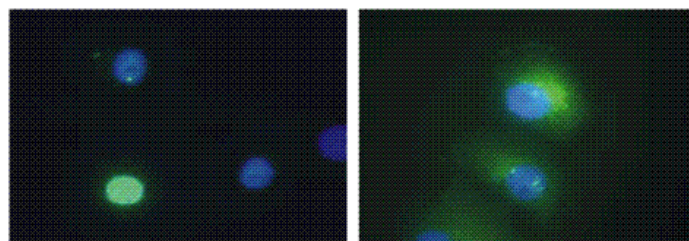
[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



**Figure:** Detection of Daxx in cell lysate of Myc-epitope tagged Daxx overexpressing TF-1 cells. Lane 1 shows endogenous Daxx detected by Daxx-01.



**Figure:** Immunofluorescence microscopy of HeLa cells transfected with pCDNA3-MycDaxx and stained with anti-Daxx (green) MAb or anti-Myc epitope MAb (blue). **Left:** nuclear localization of Myc-Daxx. **Right:** Co-transfection with pCDNA3-ASK1HA, leading to translocation of Daxx from nucleus to the cytoplasm.

## Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name	Death domain-associated protein 6
Application	ICC, WB
Clone	DAXX-01
Formulation	Liquid. In PBS, pH 7.4, containing 15mM sodium azide.
Host	Mouse
Immunogen	Recombinant human Daxx (aa 558-740).
Isotype	IgG1
Purity Detail	Protein A-affinity purified.
Species Reactivity	Human
UniProt ID	Q9UER7
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>

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)