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

This antibody is covered by our Worry-Free Guarantee.

Citations: 2

View Online »

**Ordering Information** 

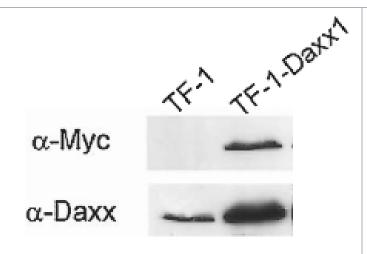
Order Online »

ALX-804-404-C100

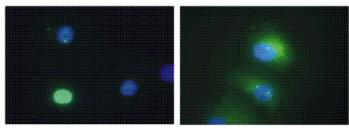
100µg

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:* Cotransfection 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 Worry-Free Guarantee

.

### Last modified: May 29, 2024



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

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