

DAXX Monoclonal Antibody [4C2]

[Catalog No. A-6217-100]

Description

Mouse monoclonal antibody raised against a partial recombinant DAXX, clone 4C2.

Size

100 µg

Purification

Protein A purified

Immunogen

DAXX (NP_001341, 561 a.a. ~ 661 a.a) partial recombinant protein with GST tag. MW of the GST tag alone is 26 KDa.

Cross Reactivity

Human

Isotype

IgG2b Kappa

Formulation

In 1x PBS, pH 7.2

Storage

Aliquot and store at -20°C or lower. Avoid repeated freezing and thawing. Multiple freeze/thaw cycles may result in decreased performance. Non-hazardous. No MSDS necessary.

Handling Recommendations

For maximum recovery of the products, centrifuge the vial prior to opening the cap.

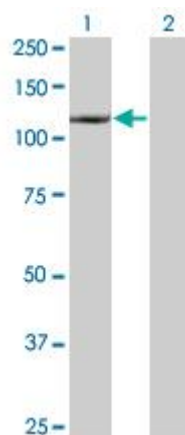
Applications & Suggested Dilutions*

Western Blot: 0.5-1 µg/ml

ELISA: 0.25-1 µg/ml

**The end user is responsible for determining optimal working dilutions.*

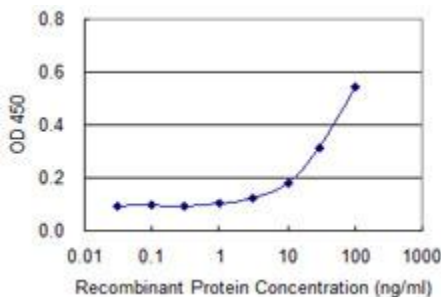
Data



Western Blot analysis of DAXX expression in transfected 293T cell line by DAXX Monoclonal Antibody, clone 4C2.

Lane 1: DAXX transfected lysate (81.373 KDa).

Lane 2: Non-transfected lysate.



Detection limit for recombinant GST tagged DAXX is approximately 1 ng/ml as a capture antibody.

This product is for research purposes only. Not intended for use in diagnostic procedures.