

TSNAX Monoclonal Antibody [4D5]

[Catalog No. A-6159-200]

Description

Mouse monoclonal antibody raised against a partial recombinant TSNAX, clone 4D5.

Size

200 µl

Purification

Protein A purified

Immunogen

TSNAX (-, 191 a.a. ~ 291 a.a) partial recombinant protein with GST tag.

Cross Reactivity

Human

Isotype

IgM Kappa

Formulation

In ascites fluid

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.

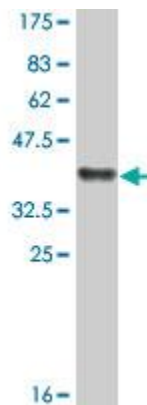
Applications & Suggested Dilutions*

Western Blot: 1:500 - 1:1000 dilution

ELISA: 1:1000 - 1:2000 dilution

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

Data



Western Blot detection against Immunogen (37.11 KDa).

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