

MUTYH Monoclonal Antibody [4D10]

[Catalog No. A-6179-100]

Description

Mouse monoclonal antibody raised against a partial recombinant MUTYH, clone 4D10.

Size

100 µg

Purification

Protein A purified

Immunogen

MUTYH (AAH03178, 436 a.a. ~ 536 a.a.) partial recombinant protein with GST tag. MW of the GST tag alone is 26 KDa.

Cross Reactivity

Human

Isotype

IgG1 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.

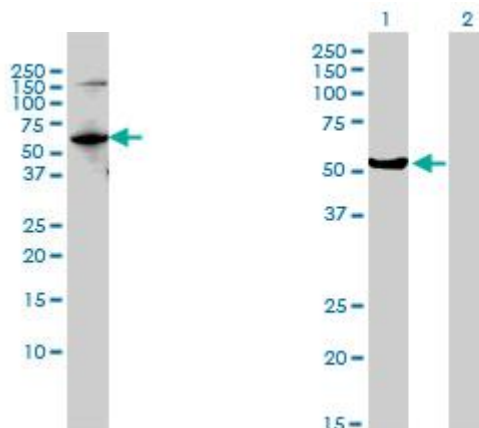
Applications & Suggested Dilutions*

Western Blot: 0.5-1 µg/ml

ELISA: 0.25-1 µg/ml

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

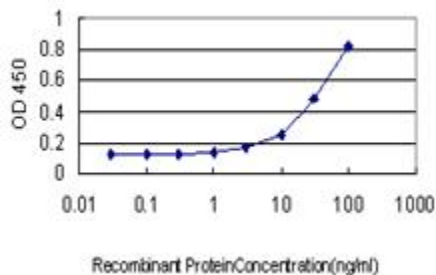
Data



MUTYH Monoclonal Antibody, clone 4D10. Western Blot analysis of MUTYH expression in HeLa NE.

Western Blot analysis of MUTYH expression in transfected 293T cell line by MUTYH Monoclonal Antibody, clone 4D10.

Lane 1: MUTYH transfected lysate (59.1 KDa).
Lane 2: Non-transfected lysate.



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

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