

UBE3A Monoclonal Antibody [2F6]

[Catalog No. A-8115-100]

Description

Mouse monoclonal antibody raised against a partial recombinant UBE3A, clone 2F6.

Size

100 µg

Purification

Protein A purified

Immunogen

UBE3A (AAH09271, 51 a.a. ~ 151 a.a) partial recombinant protein with GST tag. MW of the GST tag alone is 26 KDa.

Cross Reactivity

Human

Isotype

IgG2a 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

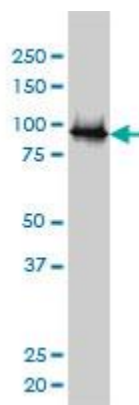
ELISA: 0.25-1 µg/ml

Immunohistochemistry (IHC): 2.5 µg/ml

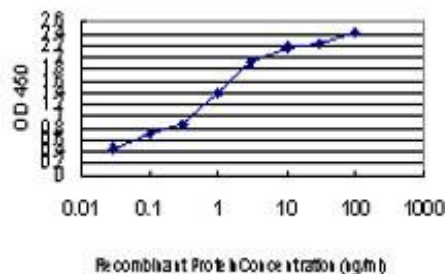
Immunofluorescence: 5 µg/ml

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

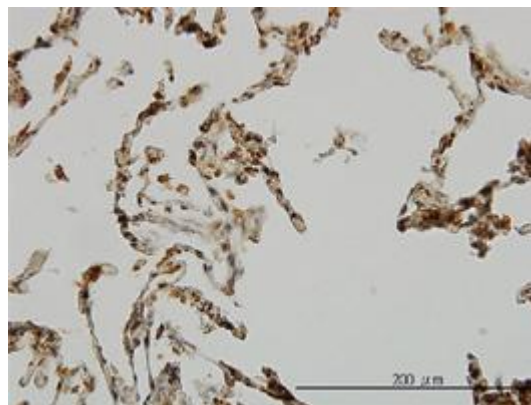
Data



UBE3A Monoclonal Antibody, clone 2F6. Western Blot analysis of UBE3A expression in Jurkat.



Detection limit for recombinant GST tagged UBE3A is approximately 0.03 ng/ml as a capture antibody.



Immunoperoxidase of monoclonal antibody to UBE3A on formalin-fixed paraffin-embedded human lung. [Antibody Concentration: 3 µg/ml]

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