

USP34 Monoclonal Antibody [2E2]

[Catalog No. A-8136-100]

Description

Mouse monoclonal antibody raised against a partial recombinant USP34, clone 2E2.

Size

100 µg

Purification

Protein A purified

Immunogen

USP34 (NP_055524, 3296 a.a. ~ 3396 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

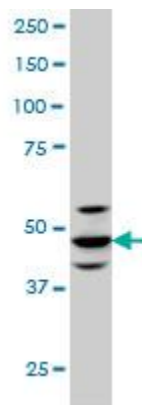
ELISA: 0.25-1 µg/ml

Immunohistochemistry (IHC): 5 µg/ml

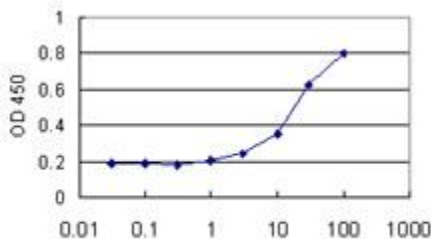
Immunofluorescence: 5 µg/ml

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

Data

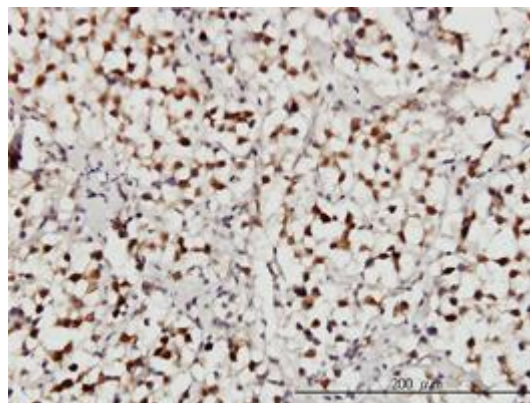


USP34 Monoclonal Antibody, clone 2E2. Western Blot analysis of USP34 expression in A-549.



Recombinant Protein Concentration (ng/ml)

Detection limit for recombinant GST tagged USP34 is approximately 1 ng/ml as a capture antibody.



Immunoperoxidase of monoclonal antibody to USP34 on formalin-fixed paraffin-embedded human seminoma. [Antibody Concentration: 3 µg/ml]

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