

SAE1 Monoclonal Antibody [1G4-1G5]

[Catalog No. A-8152-100]

Description

Mouse monoclonal antibody raised against a full length recombinant SAE1, clone 1G4-1G5.

Size

100 µg

Purification

Protein A purified

Immunogen

SAE1 (AAH00344, 1 a.a. ~ 347 a.a) full length recombinant protein with GST tag. MW of the GST tag alone is 26 KDa.

Cross Reactivity

Human

Isotype

IgG1 Lambda

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

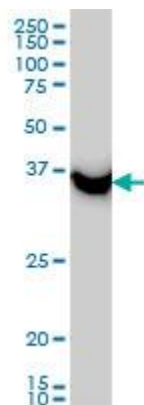
ELISA: 0.25-1 µg/ml

Immunohistochemistry (IHC): 5 µg/ml

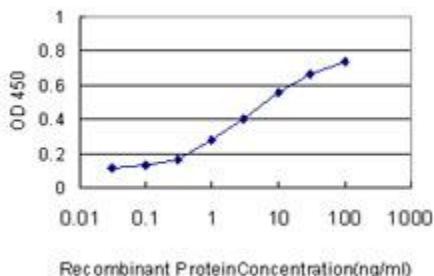
Immunofluorescence: 5 µg/ml

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

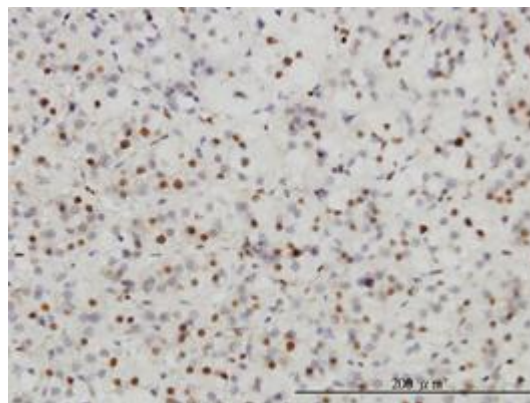
Data



SAE1 Monoclonal Antibody, clone 1G4-1G5. Western Blot analysis of SAE1 expression in HeLa.



Detection limit for recombinant GST tagged SAE1 is approximately 0.1 ng/ml as a capture antibody.



Immunoperoxidase of monoclonal antibody to SAE1 on formalin-fixed paraffin-embedded human adrenal gland. [Antibody Concentration: 3 µg/ml]

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