

ADPGK Monoclonal Antibody [1E4]

[Catalog No. A-8142-100]

Description

Mouse monoclonal antibody raised against a partial recombinant ADPGK, clone 1E4.

Size

100 µg

Purification

Protein A purified

Immunogen

ADPGK (NP_112574, 425 a.a. ~ 497 a.a) partial recombinant protein with GST tag. MW of the GST tag alone is 26 KDa.

Cross Reactivity

Human, Mouse, Rat

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.

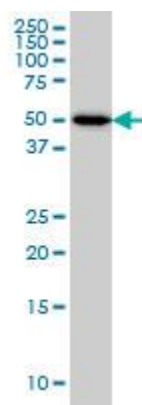
Applications & Suggested Dilutions*

Western Blot: 0.5-1 µg/ml

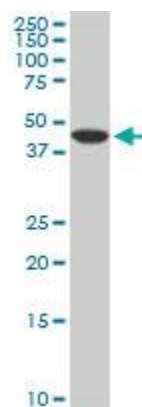
ELISA: 0.25-1 µg/ml

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

Data



ADPGK Monoclonal Antibody, clone 1E4. Western Blot analysis of ADPGK expression in HeLa.



ADPGK Monoclonal Antibody, clone 1E4. Western Blot analysis of ADPGK expression in PC-12.

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