

TTK Monoclonal Antibody [3G7]

[Catalog No. A-7146-100]

Description

Mouse monoclonal antibody raised against a partial recombinant TTK, clone 3G7.

Size

100 µg

Purification

Protein A purified

Immunogen

TTK (AAH00633, 1 a.a. ~ 176 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.

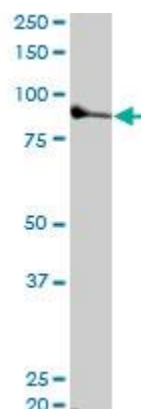
Applications & Suggested Dilutions*

Western Blot: 0.5-1 µg/ml

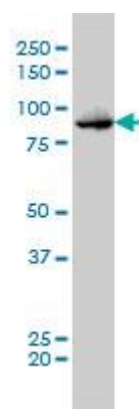
ELISA: 0.25-1 µg/ml

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

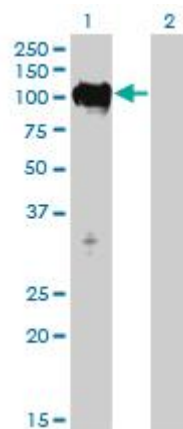
Data



TTK Monoclonal Antibody, clone 3G7. Western Blot analysis of TTK expression in HeLa.



TTK Monoclonal Antibody, clone 3G7. Western Blot analysis of TTK expression in K-562.



Western Blot analysis of TTK expression in transfected 293T cell line by TTK Monoclonal Antibody, clone 3G7.

Lane 1: TTK transfected lysate (97.1 KDa).

Lane 2: Non-transfected lysate.

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