

MAP2K7 Monoclonal Antibody [2G5]

[Catalog No. A-7158-100]

Description

Mouse monoclonal antibody raised against a partial recombinant MAP2K7, clone 2G5.

Size

100 µg

Purification

Protein A purified

Immunogen

MAP2K7 (NP_660186, 1 a.a. ~ 100 a.a) partial recombinant protein with GST tag. MW of the GST tag alone is 26 KDa.

Cross Reactivity

Human

Isotype

IgG1 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*

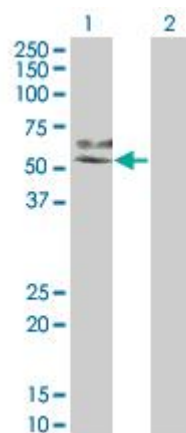
Western Blot: 0.5-1 µg/ml

ELISA: 0.25-1 µg/ml

Immunohistochemistry (IHC): 2.5 µg/ml

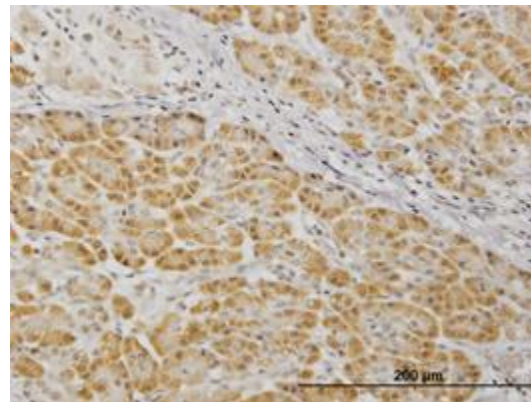
**The end user is responsible for determining optimal working dilutions.*

Data



Western Blot analysis of MAP2K7 expression in transfected 293T cell line by MAP2K7 Monoclonal Antibody, clone 2G5.

Lane 1: MAP2K7 transfected lysate (47.485 KDa).
Lane 2: Non-transfected lysate.



Immunoperoxidase of monoclonal antibody to MAP2K7 on formalin-fixed paraffin-embedded in human pancreas.
[Antibody Concentration: 1.2 µg/ml]

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