

PDCD6 Monoclonal Antibody [2B4]

[Catalog No. A-6200-100]

Description

Mouse monoclonal antibody raised against a partial recombinant PDCD6, clone 2B4.

Size

100 µg

Purification

Protein A purified

Immunogen

PDCD6 (AAH12384, 92 a.a. ~ 191 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*

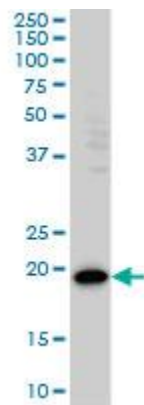
Western Blot: 0.5-1 µg/ml

ELISA: 0.25-1 µg/ml

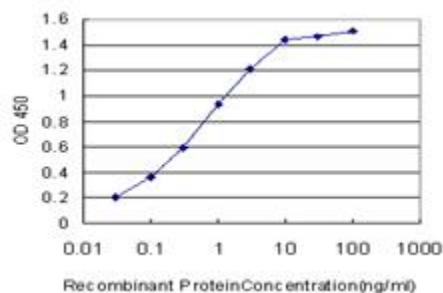
Immunofluorescence: 5 µg/ml

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

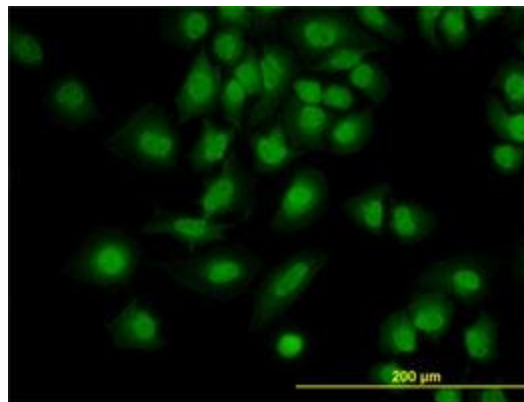
Data



PDCD6 Monoclonal Antibody, clone 2B4. Western Blot analysis of PDCD6 expression in HeLa.



Detection limit for recombinant GST tagged PDCD6 is approximately 0.03 ng/ml as a capture antibody.



Immunofluorescence of monoclonal antibody to PDCD6 on HeLa cell. [Antibody Concentration: 10 µg/ml]

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