

## CAND1 Monoclonal Antibody [5D7]

[Catalog No. A-8135-100]

### Description

Mouse monoclonal antibody raised against a partial recombinant CAND1, clone 5D7.

### Size

100 µg

### Purification

Protein A purified

### Immunogen

CAND1 (NP\_060918, 1 a.a. ~ 101 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

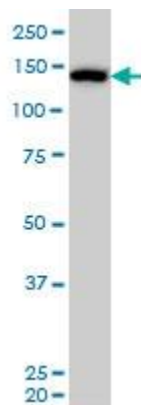
ELISA: 0.25-1 µg/ml

Immunohistochemistry (IHC): 5 µg/ml

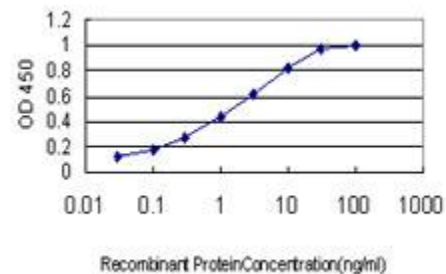
Immunofluorescence: 5 µg/ml

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

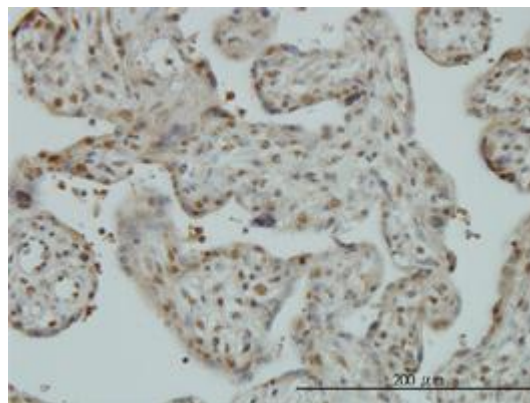
### Data



CAND1 Monoclonal Antibody, clone 5D7. Western Blot analysis of CAND1 expression in HeLa NE.



Detection limit for recombinant GST tagged CAND1 is approximately 0.03 ng/ml as a capture antibody.



Immunoperoxidase of monoclonal antibody to CAND1 on formalin-fixed paraffin-embedded human placenta. [Antibody Concentration: 3 µg/ml]

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