

## DNMT2 Monoclonal Antibody [1E12]

[Catalog No. A-1101-100]

### Background

This gene is specifically involved in the methylation of aspartic acid transfer RNA at the cytosine position 38. Although it is considered part of the family of DNA cytosine methyltransferases, DNMT2 does not methylate CpG in genomic DNA because it lacks the large N-terminal regulatory domain found in other eukaryotic methyltransferases.

### Description

Mouse monoclonal antibody raised against a full length recombinant DNMT2, clone 1E12.

### Size

100 µg

### Purification

Protein A purified

### Immunogen

DNMT2 (AAH47733, 1 a.a. ~ 392 a.a) full length recombinant protein with GST tag. MW of the GST tag alone is 26 KDa.

### Cross Reactivity

Human

### Isotype

IgG1 Lambda

### 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\*

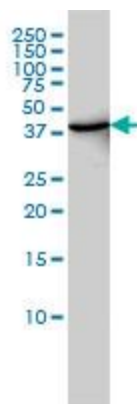
Western Blot: 0.5-1 µg/ml

ELISA: 0.25-1 µg/ml

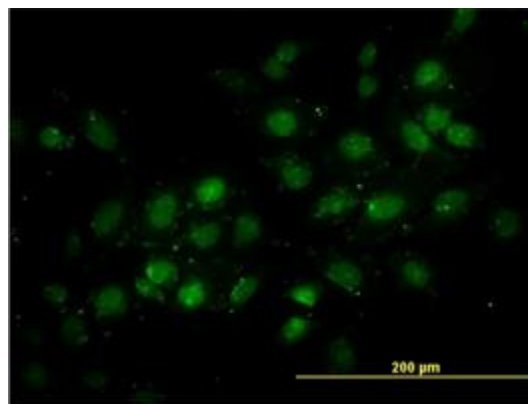
Immunofluorescence: 5 µg/ml

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

### Data



DNMT2 Monoclonal Antibody, clone 1E12. Western Blot analysis of DNMT2 expression in HeLa NE.



Immunofluorescence of monoclonal antibody to DNMT2 on HeLa cell. [Antibody Concentration: 40 µg/ml]

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