

## XRCC5 Monoclonal Antibody [3D8]

[Catalog No. A-6194-100]

### Description

Mouse monoclonal antibody raised against a full length recombinant XRCC5, clone 3D8.

### Size

100 µg

### Purification

Protein A purified

### Immunogen

XRCC5 (AAH19027, 1 a.a. ~ 733 a.a) full length 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

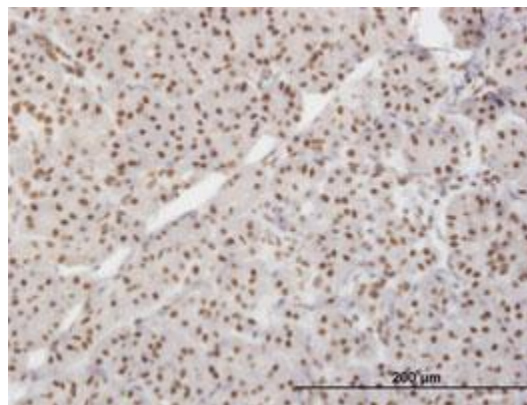
Immunohistochemistry (IHC): 2.5 µg/ml

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

### Data



*XRCC5 Monoclonal Antibody, clone 3D8. Western Blot analysis of XRCC5 expression in A-431.*



*Immunoperoxidase of monoclonal antibody to XRCC5 on formalin-fixed paraffin-embedded human pancreas. [Antibody Concentration: 3 µg/ml]*

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