

## **XRCC3 Monoclonal Antibody [1H7]**

[Catalog No. A-6191-100]

---

### **Description**

Mouse monoclonal antibody raised against a partial recombinant XRCC3, clone 1H7.

### **Size**

100 µg

### **Purification**

Protein A purified

### **Immunogen**

XRCC3 (AAH11725, 31 a.a. ~ 131 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\***

ELISA: 0.25-1 µg/ml

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

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