

## FANCG Monoclonal Antibody [2C8]

[Catalog No. A-6174-100]

### Description

Mouse monoclonal antibody raised against a full length recombinant FANCG, clone 2C8.

### Size

100 µg

### Purification

Protein A purified

### Immunogen

FANCG (AAH11623, 1 a.a. ~ 623 a.a) full length 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\*

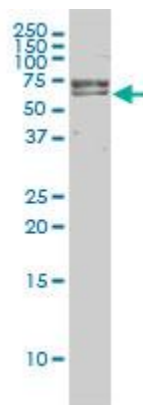
Western Blot: 0.5-1 µg/ml

ELISA: 0.25-1 µg/ml

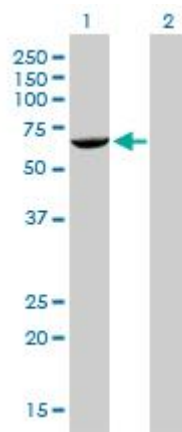
Immunofluorescence: 5 µg/ml

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

### Data



*FANCG Monoclonal Antibody, clone 2C8. Western Blot analysis of FANCG expression in HeLa NE.*



*Western Blot analysis of FANCG expression in transfected 293T cell line by FANCG Monoclonal Antibody, clone 2C8.*

*Lane 1: FANCG transfected lysate (68.6 KDa).  
Lane 2: Non-transfected lysate.*

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