

TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

XPF Monoclonal Antibody (clone 3A11)

Catalog #: 4423-AMC-100

Size: 100 µl

Background: XPF (5' endonuclease) /ERCC1 (Excision Repair Cross Complementing 1) forms a heterodimer complex and catalyzes 5' incision to the DNA damage during nucleotide excision repair (NER). ERCC1/XPF is also required for repair of DNA interstrand crosslinks (ICLs) (1) and involved in repair of double strand breaks (2).

Physical State: Protein G affinity purified mouse monoclonal antibody raised against full length recombinant human XPF/ERCC1 complex. Provided at 1.0 µg/µl in phosphate buffered saline buffer with 0.1% sodium azide.

Specificity: Recognizes human XPF.

IgG class: IgG₁

Storage: Store at 4°C for at least 1 month. For long time storage, aliquote and store at -20 °C to avoid repeated freeze-thawing.

Applications: Suitable for Western blotting and immunocytochemistry (IC). For Western blot analysis, a starting dilution of 1:500 – 1:1,000 is recommended, whereas for IC, a starting dilution of 1:100 is recommended. Empirical determination of antibody concentration is required for optimal results.

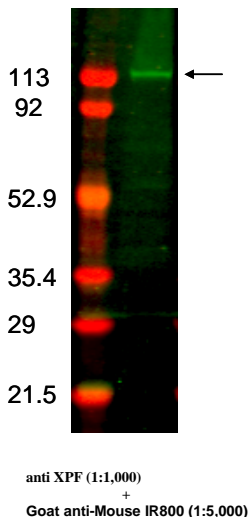


Fig. 1. Western blot analysis of A375 cells using Trevigen's anti-XPF antibody. Cells were lysed in Trevigen Lysis Reagent supplement with Protease Inhibitor Cocktail (Thermo Scientific) at the concentration 1×10^7 cells/ml. 10 µl of each lysate were loaded per well of 4-20% Tris-Glycine gel. Proteins were transferred onto an Immobilon FL membrane and XPF was detected with anti-XPF antibody (Cat# 4423-AMC-100) followed by IR800-conjugated secondary anti-body (Licor). Membranes were scanned using an Odyssey Infrared Imaging System (Licor).

References:

1. Niedernhofer, L.J. et al. (2004) *Mol Cell Biol* 24, 5776–87.
2. Ahmad, A. et al. (2008) *Mol Cell Biol* 28, 5082–92.

Related Products:

Catalog #	Description	Size
4422-AMC-100	Anti-XPF monoclonal (1E4)	100 µl
4424-AMC-100	Ant-ERCC1 monoclonal (3G7)	100 µl
6370-MC-100	Anti-human/murine-Cytochrome C	100 µg
6380-MC-100	Anti-human/murine-Holocytochrome C	100 µg
2291-MC-100	Anti-human-Bcl-2 mAb (clone YTH-8C8)	100 µg
2281-MC-100	Anti-human-Bax mAb (clone YTH-6A7)	100 µg
6361-PC-100	Anti-human/mouse-PBR polyclonal	100 µl
4335-MC-100	Anti-PAR polymer mAb (10HA)	100 µl
4336-BPC-100	Anti- PAR polymer polyclonal	100 µl
4338-MC-50	Anti-human/murine-PARP mAb (clone C2-10)	50 µg

Anti-XPF**Monoclonal Antibody****Catalog #: 4423-AMC-100****Storage: -20 °C****TREVIGEN®****1-800-873-8443**