



## XPA (Xeroderma pigmentosum group A protein) Monoclonal Antibody

Cat No: G01068 - 100 µL

### General Data

<b>Shipping:</b>	wet ice
<b>Stability:</b>	Store at -20degrees; shelf life 3 year maximum after production Avoid repeat freezing and thawing cycles. Does not contain any preservative
<b>Formulation:</b>	Liquid Ascite, does not contain any preservative therefore avoid repeat freezing and thawing cycles
<b>Host:</b>	Mouse
<b>Antigen:</b>	synthetic peptide
<b>Clone:</b>	1E11
<b>Isotype:</b>	IgG1, κ



<b>Application(s):</b>	ELISA
	Western Blot
	Immunoprecipitation
	Immunocytochemistry
	Recommended dilutions: 1/500-1/5000

**Specificity:** human XPA.

### Scientific Literature

Riedl T, Hanaoka F, Egly JM. The comings and goings of nucleotide excision repair factors on damaged DNA. EMBO J. 2003 Oct 1;22(19):5293-303.

Alekseev S, Ayadi M, Brino L, Egly JM, Larsen AK, Coin F. A Small Molecule Screen Identifies an Inhibitor of DNA

FP/20/20

For research laboratory use only – Not for human diagnostic use.

Buyers agree to purchase the material subject to Purchasing Terms that can be found on our website. Seek appropriate training to safely handle this product under normal conditions. Use the recommended personal protective equipment to prevent chemical exposures.

Bertin Bioreagent does not make any other warranty or representation whatsoever whether expressed or implied, with respect to these products. In no event will Bertin Bioreagent be liable for incidental, consequential or punitive damages.

#### Contact Bertin Bioreagent

Parc d'Activités du Pas du Lac  
10 bis avenue Ampère  
78180 Montigny le Bretonneux - France  
Tel.: +33 (0)139 306 036

<https://www.bertin-bioreagent.com/pa206/contact-us>  
<https://www.bertin-bioreagent.com>