



## DNA Damage Antibodies

Catalog #	Description	Clone	Size	Specificity	Subtype	IHC	WB	IP	E	ICC
<b>4335-MC-100-AC</b>	Anti-PAR w/Control	10HA	100 µl	U	IgG3	●	●	●	●	●
<b>4335-MC-100</b>	Anti-PAR	10HA	100 µl	U	IgG <sub>3</sub>	●	●	●	●	●
<b>4335-AMC-050</b>	Anti-PAR	10HA (AP)	50 µl	U	IgG3	●	●	●	●	●
<b>4336-APC-050</b>	Anti-PAR	poly. (AP)	50 µl	U	n/a		●		●	
<b>4336-BPC-100</b>	Anti-PAR	polyclonal	100 µl	U	n/a		●		●	
<b>4338-MC-50</b>	Anti-PARP	C2-10	50 µl	H,M,R,S,Ha	IgG <sub>1</sub>	●	●			●
<b>4354-MC-050</b>	Anti-8-oxo-dG	2E2	50 µl	U,FB	IgG <sub>2a</sub>	●			●	●
<b>4360-MC-100</b>	Anti-BPDE	8E11	100 µl	U	IgG <sub>1/k</sub>				●	
<b>4418-APC-020</b>	Anti-γ-H2AX	poly. (AP)	20 µl	H,M	n/a		●		●	●
<b>4418-APC-100</b>	Anti-γ-H2AX	poly. (AP)	100 µl	H,M	n/a		●		●	●
<b>4422-AMC-100</b>	Anti-XPF	2B3	100 µl	H	IgG <sub>2a</sub>		●			●
<b>4424-AMC-100</b>	Anti-ERCC1	2C6	100 µl	H	IgG <sub>2a</sub>		●			●

## Apoptosis Detection Antibodies

Catalog #	Description	Clone	Size	Specificity	Subtype	IHC	WB	IP
<b>2280-MC-100</b>	Anti-Bax	YTH-5B7	100 µg	M	IgG <sub>1</sub>	●*	●	●
<b>2281-MC-100</b>	Anti-Bax	YTH-6A7	100 µg	H, M, R	IgG <sub>1</sub>	●*	●	●
<b>2282-MC-100</b>	Anti-Bax	YTH-2D2	100 µg	H	IgG <sub>1</sub>	●*	●	●
<b>2291-MC-100</b>	Anti-Bcl-2	YTH-8C8	100 µg	H	IgG <sub>1</sub>		●	●
<b>2300-MC-100</b>	Anti-Bcl-X <sub>L</sub>	YTH-2H12	100 µg	H,M,R	IgG <sub>2a</sub>	●	●	●
<b>2305-PC-020</b>	Anti-Cleaved Caspase 3	poly. (AP)	20 µl	H,M	n/a	●	●	●
<b>2305-PC-100</b>	Anti-Cleaved Caspase 3	poly. (AP)	100 µl	H,M	n/a	●	●	●

## Control

Catalog #	Description	Clone	Size	Specificity	Subtype	IHC	WB	IP
<b>2275-PC-020</b>	Anti-G3PDH	polyclonal	20 µl	H,M	n/a		●	
<b>2275-PC-100</b>	Anti-G3PDH	polyclonal	100 µl	H,M	n/a		●	

IHC-Immunohistochemistry  
 WB-Western Blot  
 IP-Immunoprecipitation  
 E-ELISA  
 ICC-Immunocytochemistry

Mm-Mammalian  
 H-Human  
 M-Mouse  
 U-Species Independent

AP-Affinity Purified HA-Hamster  
 Y-Yeast  
 D-*D. melanogaster*  
 X-*X. laevis*

S-Monkey  
 FB-Free Base  
 R-Rat

\* In a cell system that overexpressed BAX protein

## Trevigen, Inc.

8405 Helgerman Court, Gaithersburg, MD 20877 USA

Phone: 1-800-TREVIGEN (1-800-873-8443) • 301-216-2800 • Fax: 301-560-4973 • Email: [info@trevigen.com](mailto:info@trevigen.com) • <http://www.trevigen.com>