

Instruction Manual

Catalog Number	PK-CA577-K253																																			
Description	<p>Apurinic/aprimidinic (AP) sites are one of the major types of DNA lesions formed during the course of base excision and repair of oxidized, deaminated or alkylated bases. It has been estimated that about 2×10^5 base lesions are generated per cell per day. The level of AP sites in cells can thus be a good indicator of DNA lesion and repair against chemical damage and cell aging. The DNA Damage Detection Kit utilizes the ARP (Aldehyde Reactive Probe) reagent that reacts specifically with an aldehyde group which is the open ring form of the AP sites. After treating DNA containing AP sites with ARP reagents, AP sites are tagged with biotin residues, which can be quantified using avidin-biotin assay followed by a colorimetric detection. The kit provides the necessary reagents for convenient determination of abasic sites in purified DNA sample in 96-well plate format.</p>																																			
Quantity	25 assays																																			
Kit Components	<table border="1"> <thead> <tr> <th>Component</th> <th>Quantity</th> <th>Color Code</th> </tr> </thead> <tbody> <tr> <td>ARP Solution (10 mM, in DMSO)</td> <td>0.125 ml</td> <td>Red</td> </tr> <tr> <td>TE Buffer</td> <td>30 ml</td> <td>NM</td> </tr> <tr> <td>Glycogen Solution (10 $\mu\text{g}/\mu\text{l}$)</td> <td>0.1 ml</td> <td>Blue</td> </tr> <tr> <td>0 ARP-DNA Standard (0.5 $\mu\text{g}/\text{ml}$)</td> <td>0.6 ml</td> <td>Clear</td> </tr> <tr> <td>40 ARP-DNA Standard (0.5 $\mu\text{g}/\text{ml}$)</td> <td>0.6 ml</td> <td>Yellow</td> </tr> <tr> <td>DNA Binding Solution</td> <td>10 ml</td> <td>WM</td> </tr> <tr> <td>HRP-Streptavidin</td> <td>0.10 ml</td> <td>Green</td> </tr> <tr> <td>10X Wash Buffer</td> <td>30 ml</td> <td>WM</td> </tr> <tr> <td>HRP Developer</td> <td>10 ml</td> <td>Brown/NM</td> </tr> <tr> <td>96-well Microplate (8 x 12 strips)</td> <td>1 plate</td> <td>Clear</td> </tr> </tbody> </table>	Component	Quantity	Color Code	ARP Solution (10 mM, in DMSO)	0.125 ml	Red	TE Buffer	30 ml	NM	Glycogen Solution (10 $\mu\text{g}/\mu\text{l}$)	0.1 ml	Blue	0 ARP-DNA Standard (0.5 $\mu\text{g}/\text{ml}$)	0.6 ml	Clear	40 ARP-DNA Standard (0.5 $\mu\text{g}/\text{ml}$)	0.6 ml	Yellow	DNA Binding Solution	10 ml	WM	HRP-Streptavidin	0.10 ml	Green	10X Wash Buffer	30 ml	WM	HRP Developer	10 ml	Brown/NM	96-well Microplate (8 x 12 strips)	1 plate	Clear		
Component	Quantity	Color Code																																		
ARP Solution (10 mM, in DMSO)	0.125 ml	Red																																		
TE Buffer	30 ml	NM																																		
Glycogen Solution (10 $\mu\text{g}/\mu\text{l}$)	0.1 ml	Blue																																		
0 ARP-DNA Standard (0.5 $\mu\text{g}/\text{ml}$)	0.6 ml	Clear																																		
40 ARP-DNA Standard (0.5 $\mu\text{g}/\text{ml}$)	0.6 ml	Yellow																																		
DNA Binding Solution	10 ml	WM																																		
HRP-Streptavidin	0.10 ml	Green																																		
10X Wash Buffer	30 ml	WM																																		
HRP Developer	10 ml	Brown/NM																																		
96-well Microplate (8 x 12 strips)	1 plate	Clear																																		
Applications / Assay Protocol	<p>DNA Damage Quantification Protocol:</p> <p>A. Purification of Genomic DNA: Several different methods and products are available for isolating genomic DNA. Among all the methods, the guanidine/detergent lysis method is simple, and it gives highly purified genomic DNA for the ARP-based abasic sites detection. During the purification process, avoid heating of the DNA solution. Determine the concentration and purity of the purified genomic DNA using the spectrophotometer*. Dissolve the genomic DNA in TE at concentration of 0.1 $\mu\text{g}/\mu\text{l}$. It is important for an accurate assay that the DNA concentration is adjusted exactly to 0.1 $\mu\text{g}/\mu\text{l}$. $1 \text{ OD}_{260 \text{ nm}} = 50 \mu\text{g}/\text{ml}$ for genomic DNA. The ratio of $\text{OD}_{260 \text{ nm}}/\text{OD}_{280 \text{ nm}}$ of highly purified DNA solution is 1.7 or higher. Protein contamination in the sample solution may cause a positive error. Note: Genomic DNA Isolation Kit (Cat.# PK-CA577-K281) is also available. For a positive control, cells may be treated using 10 mM H_2O_2 for 1 hour at 37°C to induce AP site formation.</p> <p>B. ARP Labeling:</p> <ol style="list-style-type: none"> Mix 5 μl of 0.1 $\mu\text{g}/\mu\text{l}$ purified sample DNA solution with 5 μl ARP Solution at the bottom of a microcentrifuge tube and incubate at 37°C for 1 hour to tag the DNA AP site. Add 88 μl TE and 2 μl Glycogen to the reaction solution, mix well. Add 0.3 ml of pure ethanol (not provided) mix well and keep at -20°C for 10 minutes. Centrifuge with microcentrifuge at the top speed for 10 minutes to precipitate the AP-site tagged DNA. Wash the pellet three times with 0.5 ml 70% ethanol. Quick spin to remove the trace amount of ethanol. Air dry the pellet for 5 minutes. The Biotin-tagged genomic DNA pellet can be used immediately or store at -20°C. The tagged DNA sample is stable for at least one year. <p>C. Determination of the number of abasic sites in DNA</p> <ol style="list-style-type: none"> Dilute the 40 ARP-DNA Standard (40 ARP sites per 10^5 bp) with 0 APR-DNA Standard to generate 200 μl each of the 0, 8, 16, 24, 32, 40 ARP-DNA solutions in microcentrifuge tubes (see table below). 																																			

ARP Number	0	8	16	24	32	40
40 ARP-DNA Standard (µl)	0	40	80	120	160	200
0 ARP-DNA Standard (µl)	200	160	120	80	40	0

- 2) Dissolve the Biotin-tagged DNA samples prepared in B with 1 ml of TE buffer (0.5 µg/ml).
- 3) Add 60 µl each of the above ARP-DNA Standards and ARP-labeled DNA samples into each well. For more accurate measurement, use three wells per sample.
- 4) Add 100 µl of the DNA Binding Solution to each well, keep the plate at room temperature overnight to allow the tagged-DNA bind on the plate surface. Keep the wells sealed.

Prepare Solutions before use:

Washing Buffer: Dilute the 10X Wash Buffer to 1X Buffer with ddH₂O (total volume 300 ml). Store this 1X Wash Buffer at room temperature.

HRP-Streptavidin Solution: Just before use, prepare 1:100 diluted working solution by diluting the 100 µl of HRP-Streptavidin with 9.9 ml 1X Wash Buffer.

- 5) Discard the DNA Binding Solution in the wells, and wash the well with 250 µl Wash Buffer 5 times.
- 6) Add 100 µl diluted HRP-Streptavidin solution to each well, and shake the plate for 1 hour at room temperature.
- 7) Discard the solution in the wells, and wash the wells with 250 µl Wash Buffer 5 times.
- 8) Add 100 µl of HRP Developer to each well, and incubate at 37°C for 1 hour.
- 9) Measure the O.D. 650 nm (within 1 hour after the incubation is ended), or add 100 µl of 1 M Sulfuric acid (or 6 M HCl) to stop the reaction. Mix well and measure O.D.450 nm. *

*Note: The value from O.D.450 nm will roughly double of that from O.D.650 nm.

10) Prepare the calibration curve using the data obtained with standard ARP-DNA solutions. Apply your sample DNA readings to the Standard Curve. The ARP numbers are the abasic sites per 10⁵ bp in the genomic DNA samples. Compare the numbers of AP sites in treated samples vs control samples to determine the level of DNA Damage.

Intended Use	For in vitro research use only. Not for diagnostic or therapeutic procedures.
Storage & Stability	Store kit at -20°C.

PromoCell GmbH

Sickingenstr. 63/65
69126 Heidelberg
Germany

Email: info@promokine.info
www.promokine.info

North America

Phone: 1 – 866 – 251 – 2860 (toll free)
Fax: 1 – 866 – 827 – 9219 (toll free)

Deutschland

Telefon: 0800 – 776 66 23 (gebührenfrei)
Fax: 0800 – 100 83 06 (gebührenfrei)

France

Téléphone: 0800 90 93 32 (ligne verte)
Téléfax: 0800 90 27 36 (ligne verte)

United Kingdom

Phone: 0800 – 96 03 33 (toll free)
Fax: 0800 – 169 85 54 (toll free)

Other Countries

Phone: +49 6221 – 649 34 0
Fax: +49 6221 – 649 34 40