

General Information

H2AX is phosphorylated rapidly and widely upon DNA double-strand breaks. Among H2AX histone variants, H2AX with phosphorylated Ser-139 is called γ H2AX. As a DNA damage marker, γ H2AX is monitored to investigate genotoxicity and carcinogenicity of chemicals. Recent studies suggest that γ H2AX is a marker for cellular senescence. The DNA Damage Detection Kit includes all necessary components, an anti- γ H2AX antibody, secondary antibody (fluorophore labeled), and blocking solution for staining which allows ready-to-use immunodetection of γ H2AX.

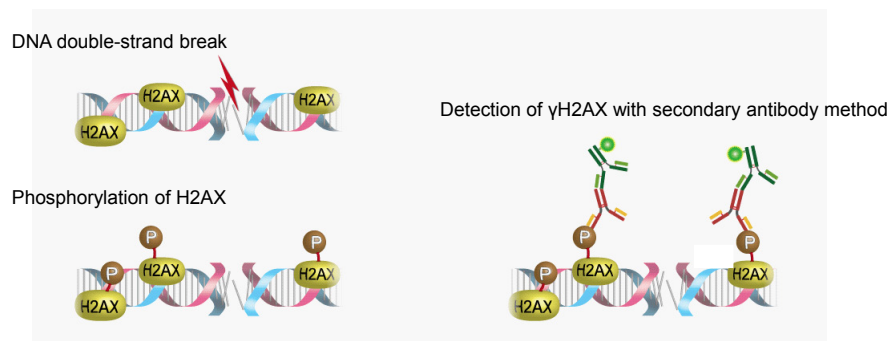


Figure 1. Detection of γ H2AX

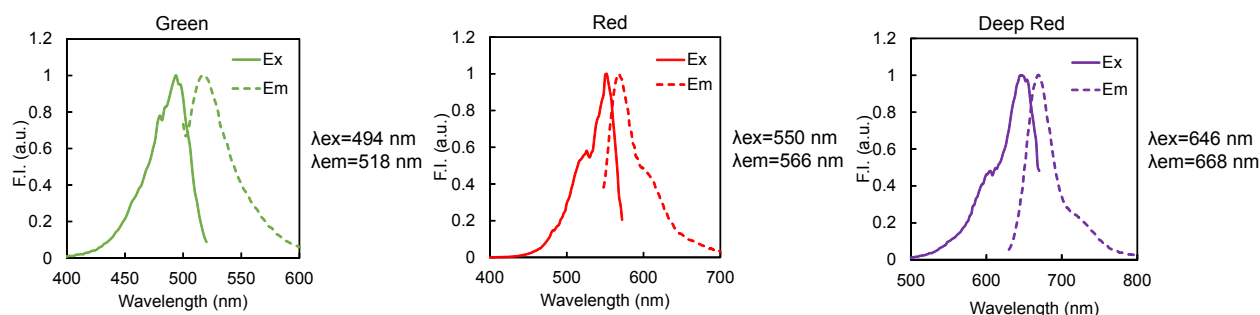


Figure 2. Excitation and emission spectra of fluorophores

Kit Contents

- Anti γ H2AX antibody (Host ; Mouse) x 1
- ※ The anti- γ H2AX antibody used in the kit is manufactured by MAB Institute, Inc.
- Secondary antibody - Green, Red or Deep Red (Host ; Goat) x 1
- Blocking Solution x 1

Storage Condition

Store at 0–5 °C

Required Equipment and Materials

- ddH₂O (double-deionized water)
- Phosphate Buffered Saline (PBS)
- Micropipette
- Paraformaldehyde (PFA)
- Triton-X
- 250 mM HEPES (pH7.4)

Preparation of Solutions

Preparation of the Anti γ H2AX antibody stock solution

Add 100 μ L ddH₂O to the Anti γ H2AX antibody and dissolve by pipetting to prepare the Anti γ H2AX antibody stock solution.

*Store the Anti γ H2AX antibody stock solution at 0–5 °C .The prepared antibody solution is stable at 4°C for 3 weeks.

Preparation of the Anti γ H2AX antibody staining solution

Dilute the Anti γ H2AX antibody stock solution by 50 times with the Blocking Solution.

*Prepare working solution fresh each day.

Preparation of the Secondary antibody staining solution

Dilute the Secondary antibody - Green, Red or Deep Red 50 times with the Blocking Solution.

*Prepare the working solution fresh each day.

Below the table shows the number of staining possible with a kit.

Staining volume	Number of staining possible
100 μ L	50
200 μ L	25
2 mL	2

1. Seed cells on a dish, chamber slide or microplate for fluorescent imaging and culture overnight at 37°C in a 37°C incubator equilibrated with 95% air, 5% CO₂.
2. Discard the supernatant and wash the cells with PBS.
3. Remove the PBS and add 250 mmol/L HEPES (pH 7.4) containing 4% PFA and 0.1% Triton X-100 for fixation. Incubate the cells at room temperature for 5 minutes .
4. Discard the supernatant and wash the cells with PBS twice.
5. Remove the PBS and add PBS containing 1% Triton X-100. Incubate the cells at room temperature for 20 minutes.
6. Discard the supernatant and wash the cells with PBS twice.
7. Add the Blocking Solution and incubate the cells at room temperature for 20 minutes .
8. Discard the supernatant and wash the cells with PBS twice.
9. Remove the PBS and add the Anti γ H2AX antibody staining solution. Incubate the cells at room temperature for 60 minutes .
10. Discard the supernatant and wash the cells with PBS twice.
11. Remove the PBS and add the Secondary antibody staining solution. Incubate the cells at room temperature for 60 minutes .
12. Discard the supernatant and wash the cells with PBS twice.
13. Observe the cells under a fluorescence microscope.

Usage Example

Detection of γ H2AX in HeLa cells treated with doxorubicin

1. HeLa cells were seeded on a μ -slide 8 well (ibidi) and cultured overnight at 37°C in a 37°C incubator equilibrated with 95% air, 5% CO₂.
2. After discarding the supernatant, medium containing 0.5 μ mol/L doxorubicin was added. The cells were incubated overnight at 37°C in the 37°C incubator equilibrated with 95% air, 5% CO₂.
3. After discarding the supernatant, the cells were washed with PBS.
4. The PBS was removed and 200 μ L 250 mmol/L HEPES (pH 7.4) containing 4% PFA and 0.1% Triton X-100 was added for fixation. The cells were incubated at room temperature for 5 minutes.
5. After discarding the supernatant, the cells were washed with PBS twice.
6. The PBS was removed and 200 μ L PBS containing 1% Triton X-100 was added. The cells were incubated at room temperature for 20 minutes.
7. After discarding the supernatant, the cells were washed with PBS twice.
8. The PBS was removed and 200 μ L Blocking Solution was added. The cells were incubated at room temperature for 20 minutes.
9. After discarding the supernatant, the cells were washed with PBS twice.
10. The PBS was removed. The Anti γ H2AX antibody staining solution was added. The cells were incubated at room temperature for 60 minutes.
11. After discarding the supernatant, the cells were washed with PBS twice.
12. The PBS was removed. The Secondary antibody staining solution was added. The cells were incubated at room temperature for 60 minutes.
13. After discarding the supernatant, the cells were washed with PBS twice.
14. The cells were observed under a fluorescence microscope.

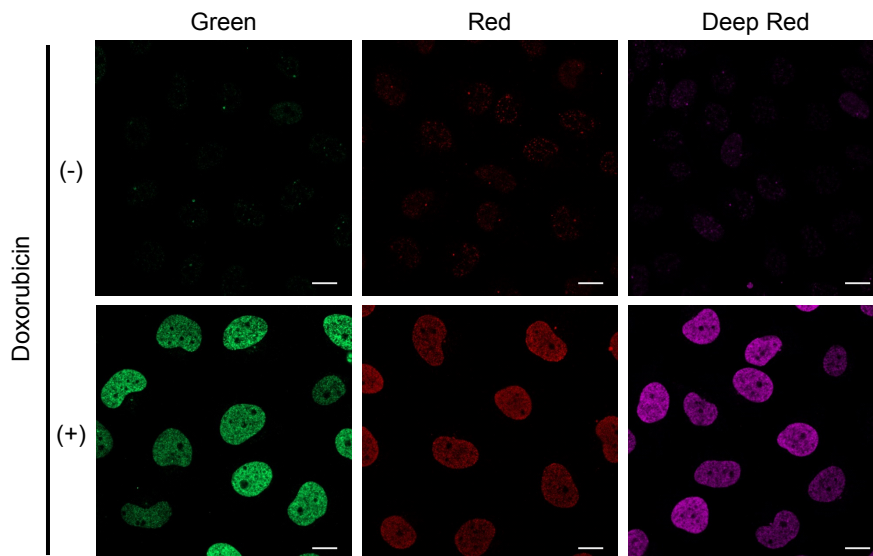


Figure 3. Detection of γ H2AX in HeLa cells treated with doxorubicin

Green Ex/Em = 488 nm/ 500–550 nm

Red Ex/Em = 561 nm/ 570–620 nm

Deep Red Ex/Em = 640 nm/ 650–700 nm

Scale bars 20 μ m

If you need more information, please contact Dojindo technical service.

Dojindo Laboratories

2025-5 Tabaru, Mashiki-machi, Kamimashiki-gun, Kumamoto
861-2202, Japan Phone: +81-96-286-1515 Fax: +81-96-286-1525
E-mail: info@dojindo.co.jp Web: www.dojindo.co.jp

Dojindo Molecular Technologies, Inc.

Tel: +1-301-987-2667 Web: <http://www.dojindo.com/>

Dojindo EU GmbH

Tel: +49-89-3540-4805 Web: <http://www.dojindo.eu.com/>

Dojindo China Co., Ltd

Tel: +86-21-6427-2302 Web: <http://www.dojindo.cn/>

G265: DNA Damage Detection - γ H2AX - Green

G266: DNA Damage Detection - γ H2AX - Red

G267: DNA Damage Detection - γ H2AX - Deep Red