

ab63801 – Anti-Rad51 Antibody

Anti-Rad51 antibody (ab63801) is a rabbit polyclonal antibody that is used to detect Rad51 in Western Blot, ICC/IF. Suitable for Human, *Xenopus laevis* samples.

For overview and additional information please visit: <http://www.abcam.com/ab63801>

Storage and Stability

The antibody may be stored at -20°C. Avoid freeze-thaw cycles. The antibody may be stored at +4°C for up to 1 – 2 weeks.

Materials Supplied

Item	Quantity	Storage Condition
Anti-Rad51 antibody	100 µL	-20°C

Protocol for the use of anti-Rad51 (human) antibody, rabbit serum, for immunostaining to observe foci-formation on chromosome after incubation of DNA-damaged cells as follows. (provided by Prof. K Suzuki)

1. Incubate cells on glass slips for at least 24 hours.
2. Wash cells with PBS once.
3. Fix cells with cold methanol for 10 min on ice.
4. Wash cells with PBS.
5. Incubate cells with 0.5% Triton X-100 for 5 min on ice.
6. Wash cells with PBS extensively.
7. Incubate cells with anti-Rad51 antibody diluted (1:100) in TBS-DT (TBS: 20 mM Tris-HCl, pH7.6, 137 mM NaCl, containing 0.1% Tween-20 and 5% Skim milk).
8. Wash cells with PBS.
9. Incubate cells with Alexa594-labeled anti-rabbit antibody diluted (1:1000) in TBS-DT

In addition to the above;

1. Different fixation methods have to be employed depending on types of cells.
2. Different blocking reagents have to be tried in case of high background.
3. In the absence of DNA-damage, Rad51 is localized diffusely in whole nucleus but Rad51 accumulates as foci on chromosome with double strand DNA breaks or stalled replication forks

For technical support contact information, visit: www.abcam.com/contactus

Copyright © 2025 Abcam, All Rights Reserved. All information / detail is correct at time of going to print.
Version 1c | 2025-July-08