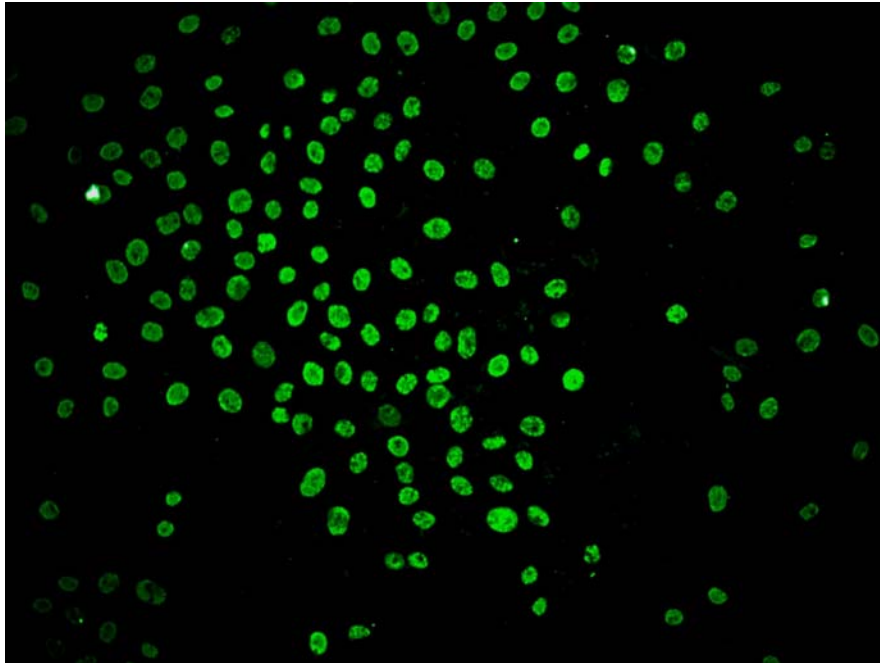
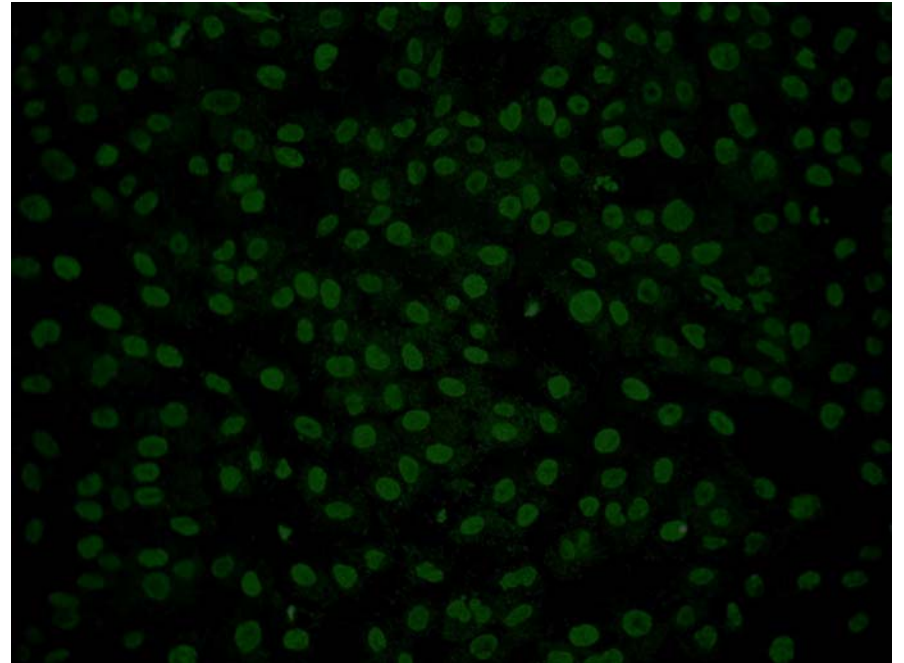


Data Images for 8-oxo-dG Antibody



A



B

H₂O₂ treated (A) and untreated (B) cells stained with 8-oxo-dG antibody according to the protocol using AlexaFluor 488 anti-mouse antibody.