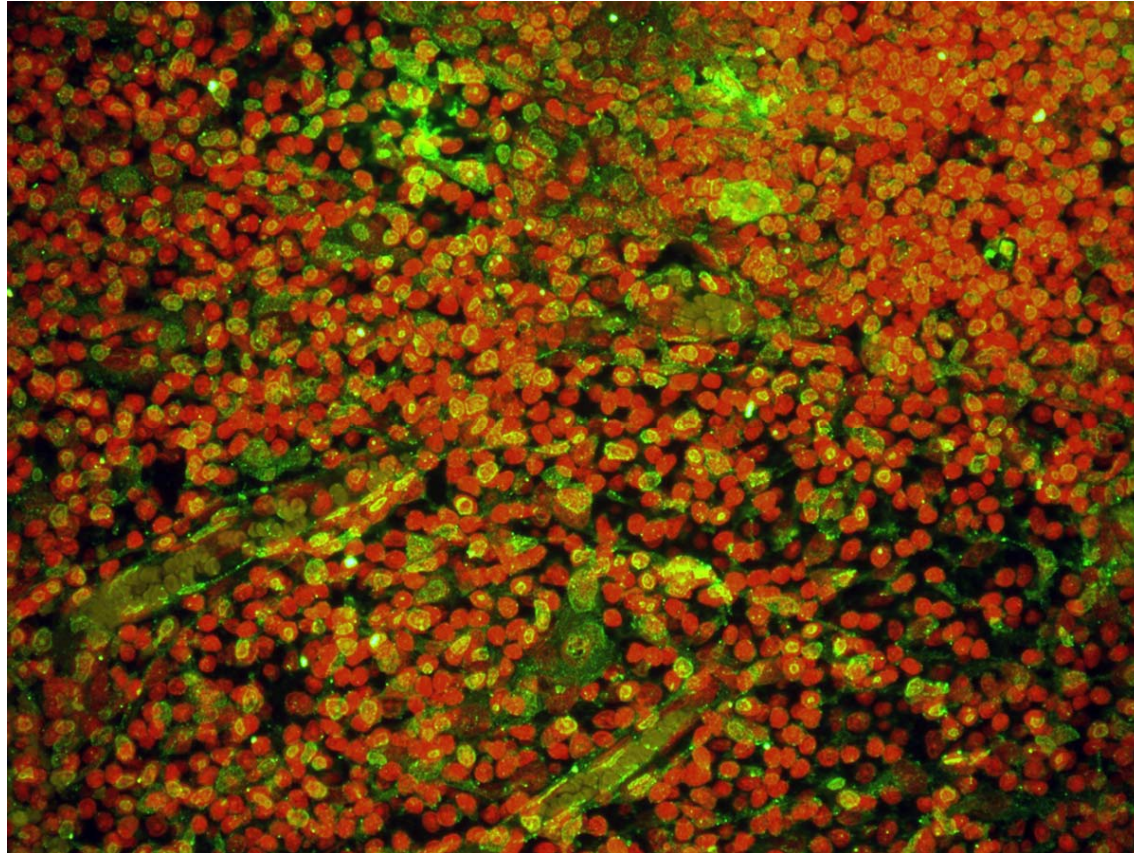


8-oxo-dG Monoclonal Antibody Staining of a Rat Thymus Section



Paraffin embedded rat thymus sections were stained with 8-oxo-dG monoclonal antibody (clone 2E2) (Cat# 4354-MC-50) at a 1:250 dilution and detected by Alexa-488, which appears fluorescent green. 8-oxo-dG is present in the cytoplasm and nucleus of most but not all cells. Sections were counterstained with 7-Aminoactinomycin D (7AAD), which labels nuclear DNA. Images were captured at 40X magnification.