



Figure 5. Immunofluorescence using the Diagenode antibody directed against Cas9

HeLa cells expressing Cas9 under the control of the tight TRE promoter were fixed in methanol at -20°C , permeabilized with acetone at -20°C and blocked with PBS containing 2% BSA. The cells were stained with the Cas9 antibody (cat. No. C15310258) diluted 1:1000, followed by incubation with a goat anti-rabbit secondary antibody coupled to AF594. Nuclei were counter-stained with Hoechst 33342. Figure 5 shows the result in the presence (left) or absence (right) of doxycycline.