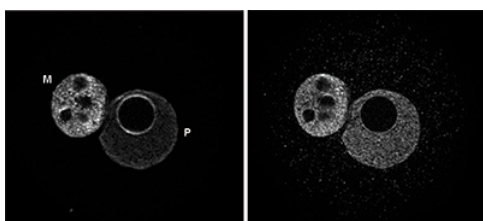


**Figure 4. FISH using the Diagenode monoclonal antibody directed against 5-mC**

To detect methylated chromosomal regions, FISH was performed on metaphase chromosomes from HeLa cells using the Diagenode monoclonal antibody against 5-mC (Cat. No. C15200003). The cells were blocked in metaphase by treatment with colcemid (0.1 µg/ml) for 1 - 2 hours, fixed overnight at -20°C with ethanol/glacial acetic acid and treated with 2N HCl for 30' at room temperature. Subsequently, the cells were blocked with PBS containing 1% BSA and 0.1% Triton X-100 and stained with the 5-mC antibody (left) diluted 1:1,000 in blocking solution, followed by an anti-mouse antibody conjugated to Alexa594. The middle panel shows staining of the chromosomes with DAPI. A merge of the two stainings is shown on the right.



**Figure 5. FISH using the Diagenode monoclonal antibody directed against 5-mC**

To detect methylated chromosomal regions, FISH was performed on metaphase chromosomes from HeLa cells using the Diagenode monoclonal antibody against 5-mC (Cat. No. C15200003). The cells were blocked in metaphase by treatment with colcemid (0.1 µg/ml) for 1 - 2 hours, fixed overnight at -20°C with ethanol/glacial acetic acid and treated with 2N HCl for 30' at room temperature. Subsequently, the cells were blocked with PBS containing 1% BSA and 0.1% Triton X-100 and stained with the 5-mC antibody (left) diluted 1:1,000 in blocking solution, followed by an anti-mouse antibody conjugated to Alexa594. The middle panel shows staining of the chromosomes with DAPI. A merge of the two stainings is shown on the right.