



Figure 2. Expression of EA restricted to the cytoplasm. Confocal images showing immunolabeling of the EA protein using PathHunter Anti-EA Antibody (PN, 93-0246) in the PathHunter CHO-K1 EA-Cyto Cell Line demonstrating the specific localization of the EA protein within the cytoplasm and the homogeneous expression of EA in the clonally derived cell population.