

SUPPLEMENTAL METHODS AND FIGURES 1 -3

SUPPLEMENTAL METHODS

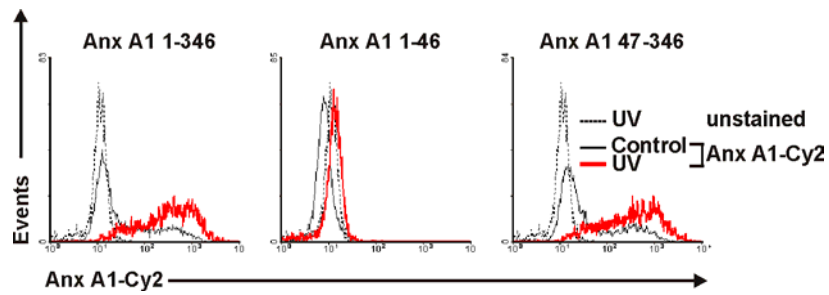
Preparation of primary human PHA-stimulated lymphoblasts, monocytes and neutrophils

Human PBMCs and neutrophils were prepared from heparinized blood by double-Ficoll gradient purification (Histopaque 1.119 g/ml, Sigma and Ficoll-Paque 1.077 g/ml, GE Healthcare). Primary human monocytes were positively selected from PBMCs with anti-CD14 magnetic beads (Miltenyi, Bergisch Gladbach, Germany) according to the manufacturer's recommendation. PHA-stimulated lymphoblasts were prepared by cultivating monocyte-depleted human PBMCs for 2 days in RPMI-1640 supplemented with 10% heat-inactivated fetal calf serum, 100 units of penicillin/ml, 0.1 mg streptomycin/ml, 10 mM HEPES and 0.5 µg/ml PHA (Biochrom, Berlin, Germany) and 5 further days in PHA-free medium.

Flow cytometric triple staining for PS exposure, membrane permeabilization and anx A1 externalization

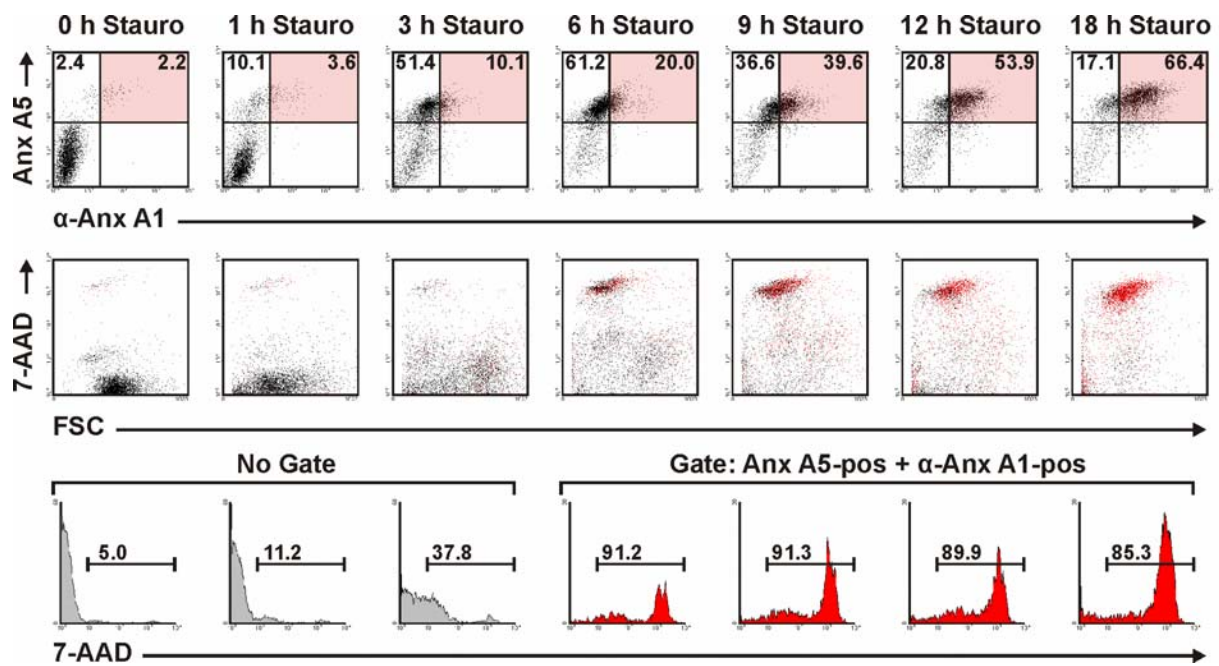
Staining with the GFP-certified apoptosis/necrosis detection kit (Enzo Life Sciences, Loerrach, Germany) was performed according to the manufacturer's instructions. Anti-anx A1 staining was combined with the procedure by pre-incubating the cells with anti-anx A1 for 30 min and subsequently adding the secondary anti-mouse-IgG-Cy2 antibody to the anx A5-Cy3/7-AAD staining mixture.

SUPPLEMENTAL FIGURE LEGENDS



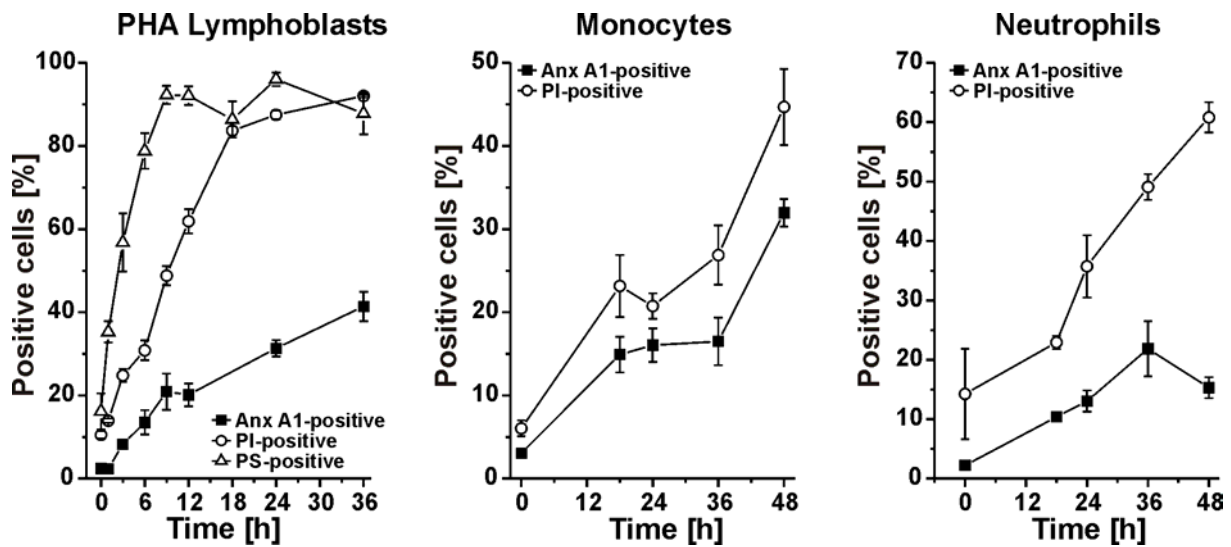
Supplemental Figure 1: Apoptotic cells bind Cy2-labeled anx A1 and an anx A1 core fragment.

THP-1 cells were left untreated or UV-irradiated with 10 mJ/cm². After 12 h of incubation cells were stained with purified Cy2-labeled anx A1 mutants (aa 1-346, aa 1-46 and aa 47-346) and analyzed by flow cytometry. Representative histograms of triplicates are shown (analysis gate was set on PI-negative cells).



Supplemental Figure 2: Time course of phosphatidylserine exposition, membrane permeabilization and anx A1 externalization during apoptosis and secondary necrosis.

Jurkat cells were stimulated with 2.5 μ M staurosporine (Stauro) for the indicated times in order to undergo apoptosis followed by secondary necrosis. Subsequently, cells were stained with anti-anx A1/anti-mouse-IgG-Cy2, anx A5-Cy3 and 7-AAD and analyzed by flow cytometry. *Upper panel:* Dotplots showing anti-anx A1 vs. anx A5 staining. *Middle panel:* Dotplots showing FSC vs. 7-AAD. Cells positive for anti-Anx A1 and anx A5 staining (upper right quadrant in upper panel) are displayed in red. *Lower panel:* Histograms of 7-AAD fluorescence. The first 3 histograms display the whole population analyzed, whereas the last 4 histograms are gated on the population positive for anti-Anx A1 and anx A5 (upper right quadrant in upper panel). Representative plots of triplicates are shown.



Supplemental Figure 3: Exposure of phosphatidylserine and loss of plasma membrane integrity precede anx A1 externalization in primary human PHA-stimulated lymphoblasts, monocytes and neutrophils.

Primary human PHA-stimulated lymphoblasts, monocytes and neutrophils were stimulated with 2.5 μ M staurosporine to undergo apoptosis and secondary necrosis. Cell surface exposure of phosphatidylserine (PS) was detected by anx A5-FITC staining. Anx A1 externalization and PI exclusion were measured as in Fig. 2A. The percentages of PS-positive, anx A1-positive and PI-positive cells (mean values \pm standard deviations of triplicates) are shown.