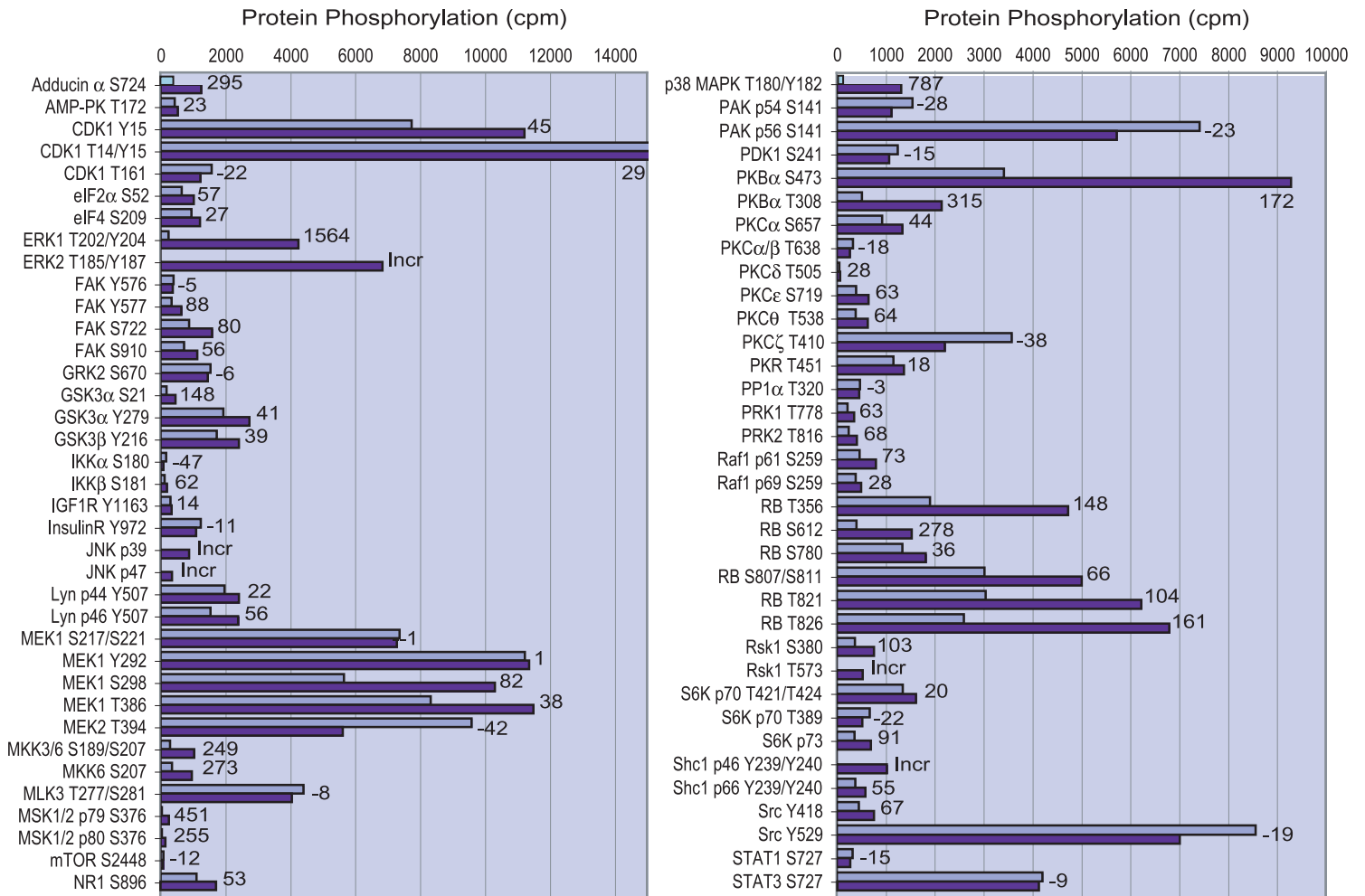


Dimethyl Varacin Effects on Protein Phosphorylation in MCF7 Cells



The combined results from the Kinetworks™ KPSS 1.3, 2.0, 3.0, and 4.0 screen analyses of the control (light bars) and those from the drug-treated (dark bars) human MCF7 breast epithelial cells were averaged and then graphed. The percentage difference between the 5 μ M dimethylvaracin for 2 h-treated and control cells is indicated at the end of the bars. Only those target phosphoproteins that were detected with these screens are shown.