



## Human MCF7 breast epithelial cells

Untreated ■ 5 uM Dimethyl Varacin for 2 h ■



**KINEXUS**

Kinetworks™ KPSS Phospho-Site Screens