



Kinetworks™ KPSS 1.3 Phospho-Site Screen Analysis
of Maturing *Xenopus laevis* Oocytes at GVBD

