

# Cell Counting Kit-8 (WST-8\*) Assay for Cell Viability

