

| PATHWAY ASSAYS |                      |                           |                      |
|----------------|----------------------|---------------------------|----------------------|
| Target         | Pathway              | Assay Measures            | Select for Profiling |
| Cdc25a         | DNA Damage           | Cdc25a Degradation        |                      |
| IκB            | NFκB                 | IκB Degradation           |                      |
| p53            | DNA Damage           | p53 Nuclear Translocation |                      |
| PI3            | PI3                  | FOXO Translocation        |                      |
| AKT            | PI3                  | FOXO Translocation        |                      |
| Wnt            | Wnt-Frizzled Pathway | β-Catenin Translocation   |                      |
| GSK3β          | Wnt-Frizzled Pathway | β-Catenin Translocation   |                      |