

Caspase-3 Activation Assay by Western Blot

[Apoptosis assays](#) are available for single and multi-time point analysis as well as real-time live-cell imaging.

Assay Properties	
Cell Line	NIH3T3 cell line from ATCC
Assay Format	Western blot
Assay Protocol	NIH3T3 cells were treated with DMSO, 20 μ M of RBC1023, 1 μ M of staurosporine, or 20 μ M of RBC1023 plus 1 μ M of staurosporine for 6 hours. Caspase-3 activation and processing was measured by Western blotting with anti-caspase-3 antibody. Tubulin was used as a protein loading control.
Readout	LI-COR Odyssey Fc Imaging System
Reference Compound	Staurosporine
Screening location	Malvern, PA, USA
Further information	More information can be found on our website Apoptosis Assay Services for Drug Discovery .

