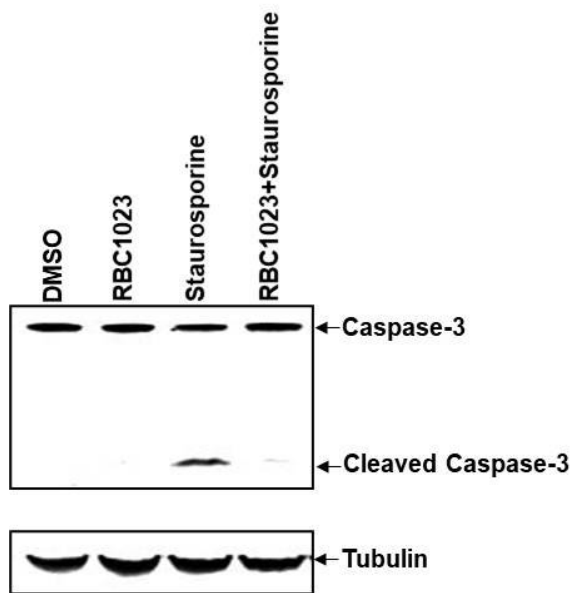


# Caspase-3 Activation Assay by Western Blot

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

## Assay Properties

<b>Cell Line</b>	NIH3T3 cell line from ATCC
<b>Assay Format</b>	Western blot
<b>Assay Protocol</b>	NIH3T3 cells were treated with DMSO, 20 $\mu$ M of RBC1023, 1 $\mu$ M of staurosporine, or 20 $\mu$ M of RBC1023 plus 1 $\mu$ 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.
<b>Readout</b>	LI-COR Odyssey Fc Imaging System
<b>Reference Compound</b>	Staurosporine
<b>Screening location</b>	Malvern, PA, USA
<b>Further information</b>	More information can be found on our website <a href="#">Apoptosis Assay Services for Drug Discovery</a> .



Let's Discover Together.

© May 2022 **Reaction Biology**. All rights reserved.

[https://www.reactionbiology.com/datasheet/caspase-3\\_activation\\_assay\\_malvern](https://www.reactionbiology.com/datasheet/caspase-3_activation_assay_malvern)

