

Caspase Activation Assay (Apoptosis)

Cat #: CAS_DEVD

Cells: Mouse Fibroblast Cells, 3T3

Description:

3T3 Cells were treated with Staurosporine or Methotrexate for overnight in the presence of 1% Fetal Bovine Serum. Lysate of the 3T3 cells were analyzed for the activity of cleavage of Ac-DEVD-AMC substrate.

Reference Compound IC50:

Assay	Caspase Activation (Ac-DEVD-AMC) (nM)
Staurosporine	2.1
Methotrexate	179

Reference Compound EC50 for
Caspase Activation in 3T3 Cells

