



# Assay kits for the ImageXpress Pico system

Easy-to-use, reliable EarlyTox™ Cell Viability Assay Kits for ImageXpress® Pico Automated Cell Imaging System

These assay kits are a family of fluorescence-based reagents for the assessment of cell viability, cell proliferation, and apoptosis events using mammalian cells. The kits employ a no-wash, homogenous assay protocol that enables characterization of a full concentration-response profile of test compounds.

Assay kit	Application	Biological marker	Key component	Part number (*Explorer/Bulk)
EarlyTox Live/Dead Assay kit	Live/Dead cell assay	Plasma membrane integrity	Calcein AM, Ethidium Homodimer-III	R8340/R8341
EarlyTox Live Cell Assay kit	Live cell/cell proliferation assay	Plasma membrane integrity	Calcein AM	R8342/R8343
EarlyTox Caspase-3/7 NucView 488 Assay kit	Caspase 3/7 apoptosis assay	Caspase 3/7 activity	NucView 488 Caspase-3 Substrate in PBS; Masking reagent	R8350/R8351
EarlyTox Caspase-3/7-D NucView 488 Assay kit	Caspase 3/7 apoptosis assay	Caspase 3/7 activity	NucView 488 Caspase-3 Substrate in DMSO; Masking reagent	R8348

\*Explorer kit is sufficient for two 96-well plates and bulk kit is enough for ten 96-well plates.

## Contact Us

Phone: [+1.800.635.5577](tel:+18006355577)  
 Web: [www.moleculardevices.com](http://www.moleculardevices.com)  
 Email: [info@moldev.com](mailto:info@moldev.com)

Check our website for a current listing of worldwide distributors.