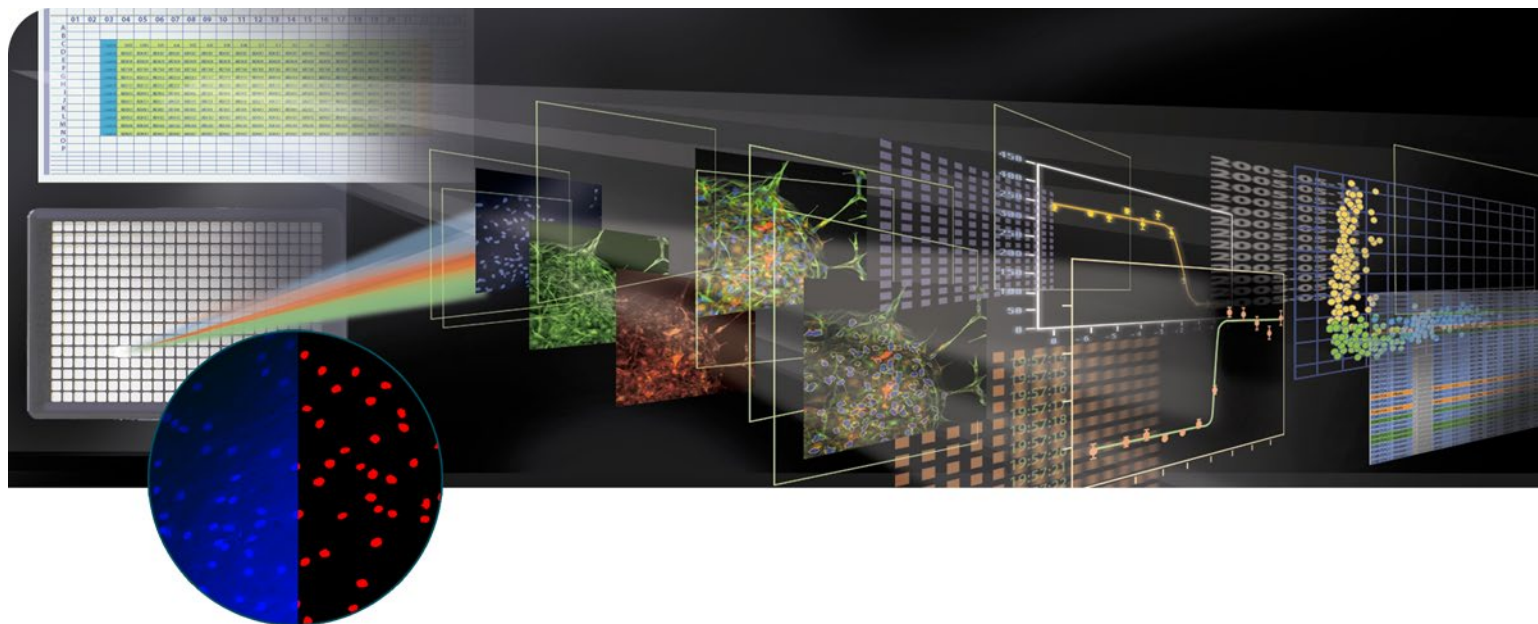


MetaXpress Software Count Nuclei Application Module

ANALYSIS SOFTWARE DROP-IN FOR METAXPRESS SOFTWARE



→ COUNT NUCLEI OF MOST CELL TYPES

→ ACCURATE SEGMENTATION OF TOUCHING CELLS

→ RUNS ACROSS ENTIRE DATA SETS AUTOMATICALLY

→ ADAPTIVE BACKGROUND CORRECTION FOR IMPROVED SEGMENTATION

The Count Nuclei Application Module for MetaXpress® Software from Molecular Devices is designed to automate accurate counting of nuclei of most types of cells.

The module accurately counts nuclei even when the background fluorescence in one imaged area is greater than the nuclei fluorescence in another area, providing superior background compensation compared to simple thresholding.

The module utilizes Adaptive Background Correction (ABC) which adapts the detection algorithm to the local intensity ranges between and within cells to provide the most robust segmentation available in an image-based screening system.

ABC enables probe detection even with highly variable background fluorescence within a single image.

This simple yet powerful module has a user-friendly interface to minimize set-up efforts, while providing the flexibility to customize the settings to the type of cells used in a specific experiment. Analysis set-up can be done in two easy steps, and the settings can be applied to an entire plate with one click.

