



Neurite Tracing Application Module

for MetaMorph NX Microscopy Automation and Image Analysis Software

New in MetaMorph® NX Software 2.5, the Neurite Tracing Application Module simplifies the meticulous tasks associated with neuronal research. For simple analyses, quickly segment, then measure overlapping neurons and their outgrowths acquired with a one-wavelength image. For more advanced, multifaceted analyses, the module can just as quickly also identify spines, perform a Sholl Analysis to obtain precise data about neuron outgrowth patterns, and identify nuclei in images acquired with a nuclear stain.

Benefits

- Effortlessly conduct simple and advanced neurite analyses with one versatile tool
- Instantly delineate intricate details of elaborate, overlapping neuronal networks with color-coded segmentation
- Accurately analyze images of any size acquired using transmitted light or a fluorescent probe