

Visualization and High-throughput Quantification of Akt Activity in Live-Cell Neuroinflammatory Models

J. Trigg^{1*}, G. Lovell¹, A. Ward¹, J. Rauch², G. Filonov², S. L. Alcantara¹, and T. Dale¹

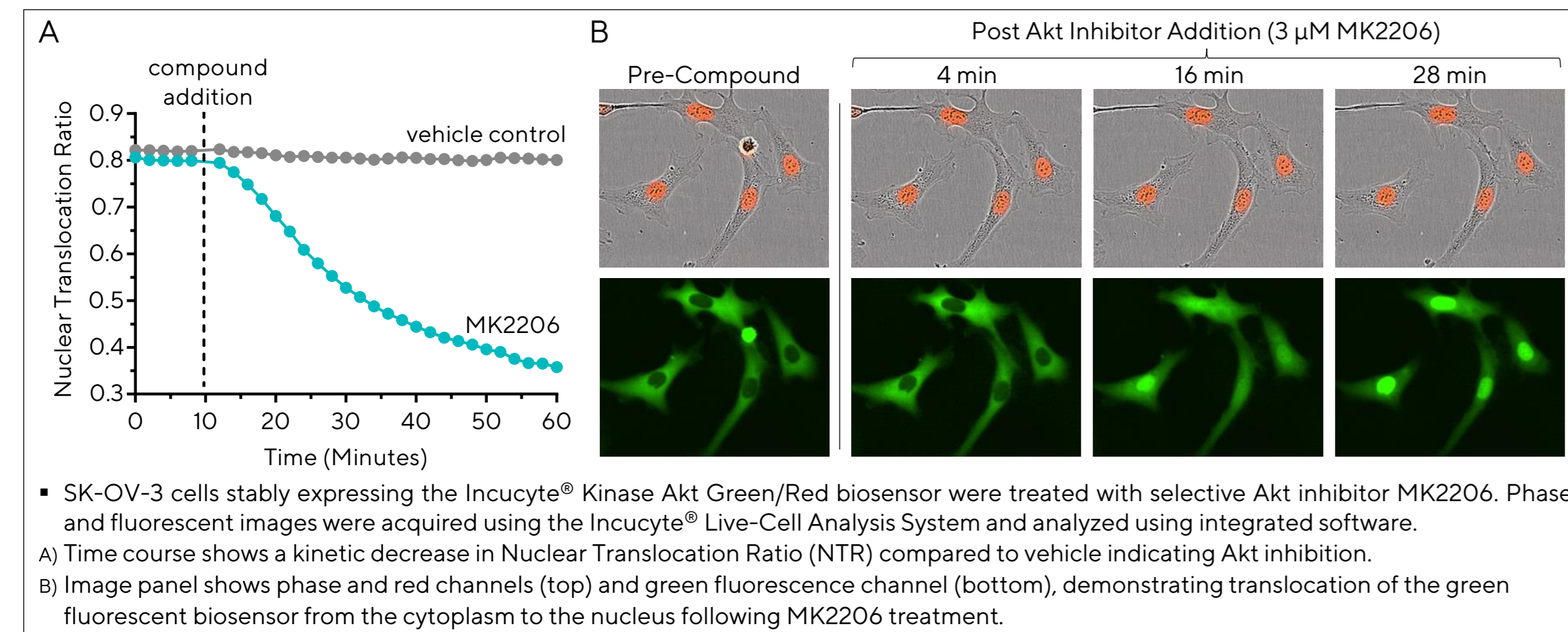
Sartorius Corporation; ¹Royston UK, ²Ann Arbor MI
*Corresponding author: Jasmine.Trigg@sartorius.com

Introduction

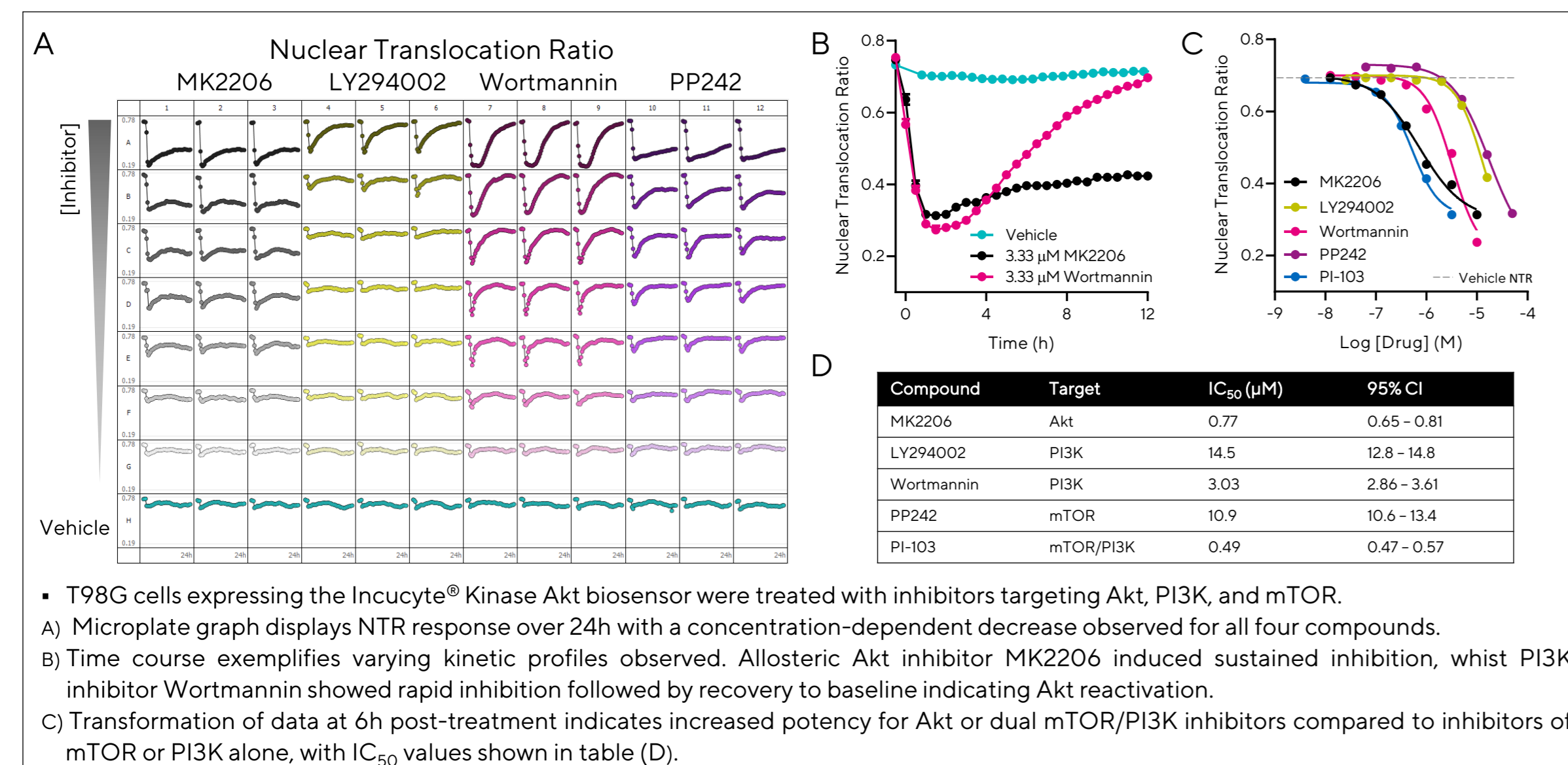
- Chronic neuroinflammatory states are associated with the development of several neurodegenerative diseases. The PI3K/Akt signaling pathway has been implicated in these disease processes and holds therapeutic promise as a target to modulate neuroinflammatory responses.
- Here we demonstrate a robust *in vitro* assay to assess dynamic changes in Akt activity in real-time.
- To monitor Akt we used astrocytic and immune cells stably expressing the Incucyte[®] Kinase Akt Green/Red biosensor, a genetically-encoded fluorescent kinase translocation reporter

- whose subcellular localization is dependent on phosphorylation by Akt.
- We investigated the effects of inhibitors targeting the PI3K/Akt pathway and observed differential time- and concentration-dependent responses. We also examined the effects of immunocompetent cell activation on Akt using LPS stimulation and inflammatory cytokines.
- These data exemplify the Incucyte[®] Kinase Akt Assay as a powerful live-cell approach for assessing Akt activity in neuroinflammatory models and its amenability to screening of therapeutic candidates.

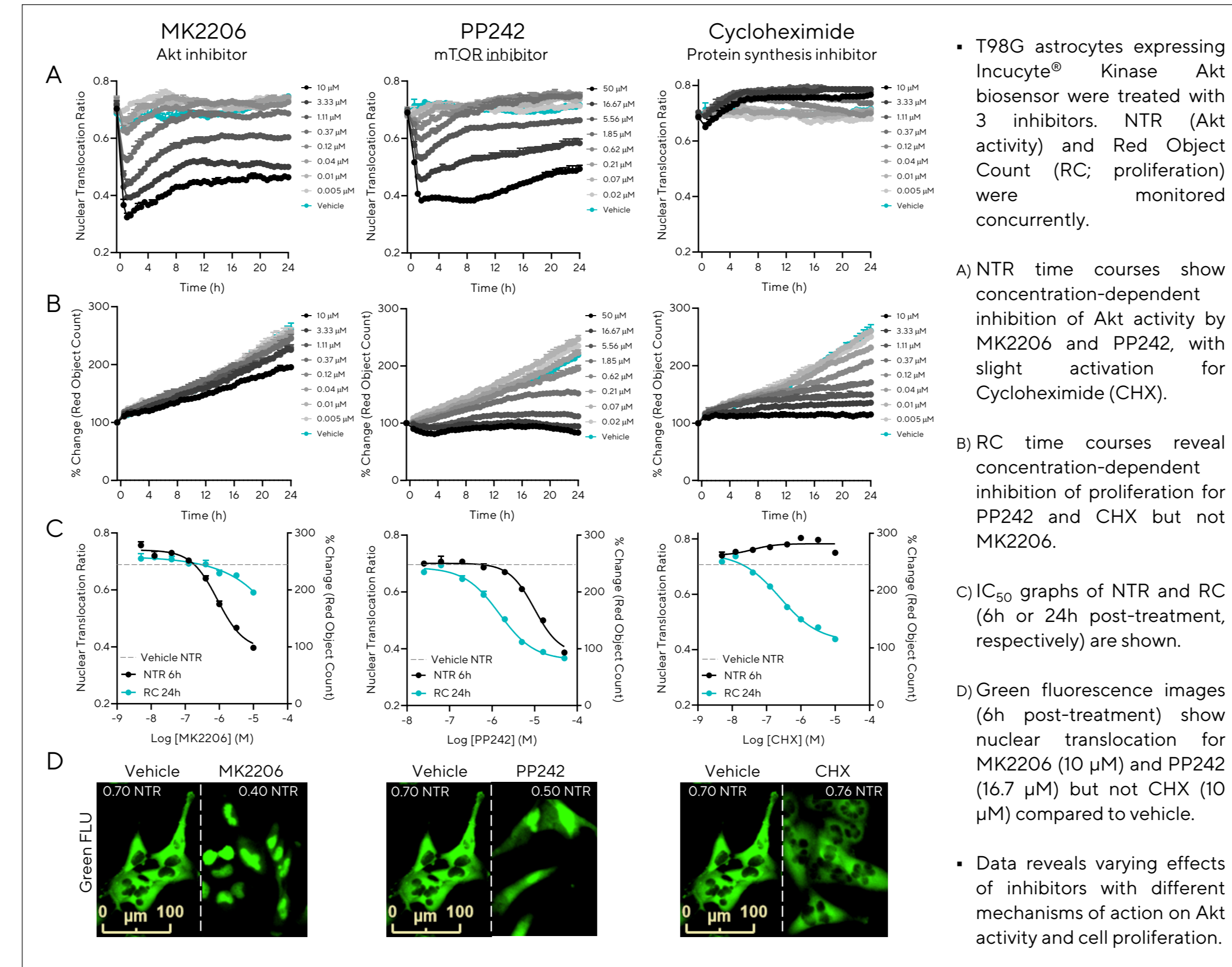
Live-cell imaging of Akt kinase activity



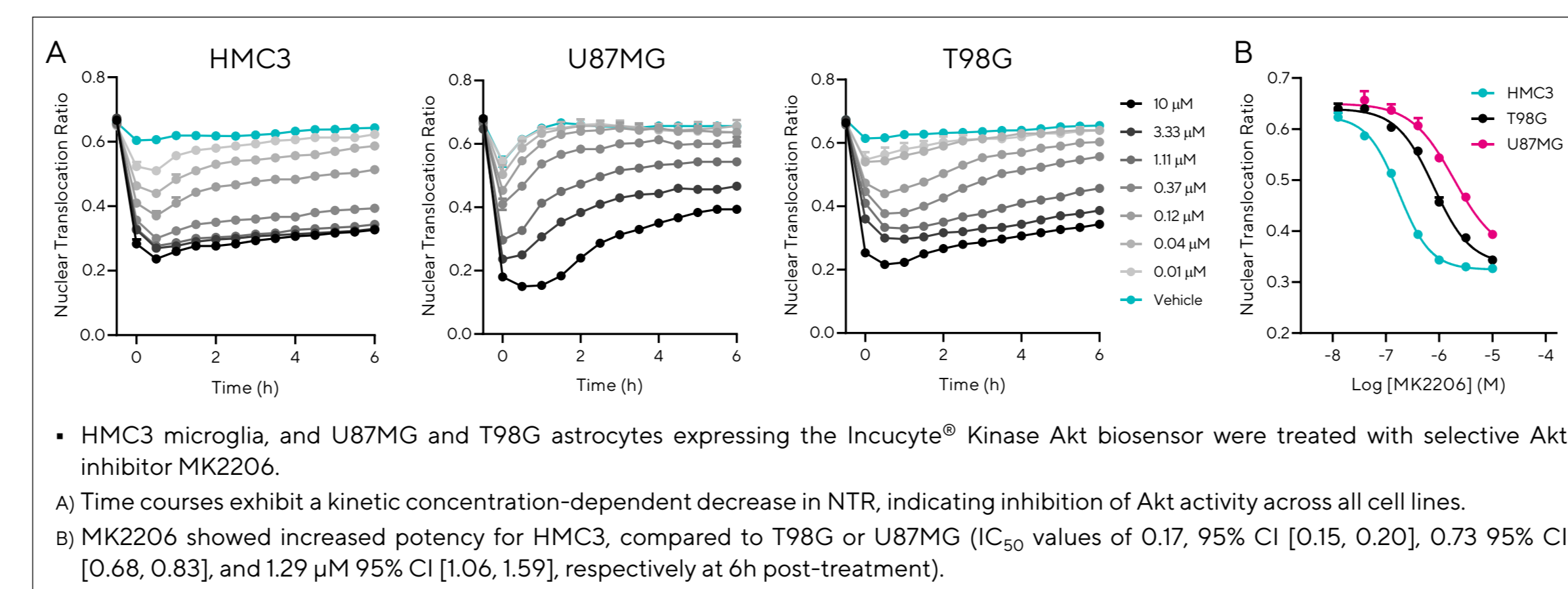
Direct and indirect inhibitory compound effects on Akt activity



Concurrent quantification of Akt activity and cell proliferation



Comparison of Akt inhibitor MK2206 across neural cell types



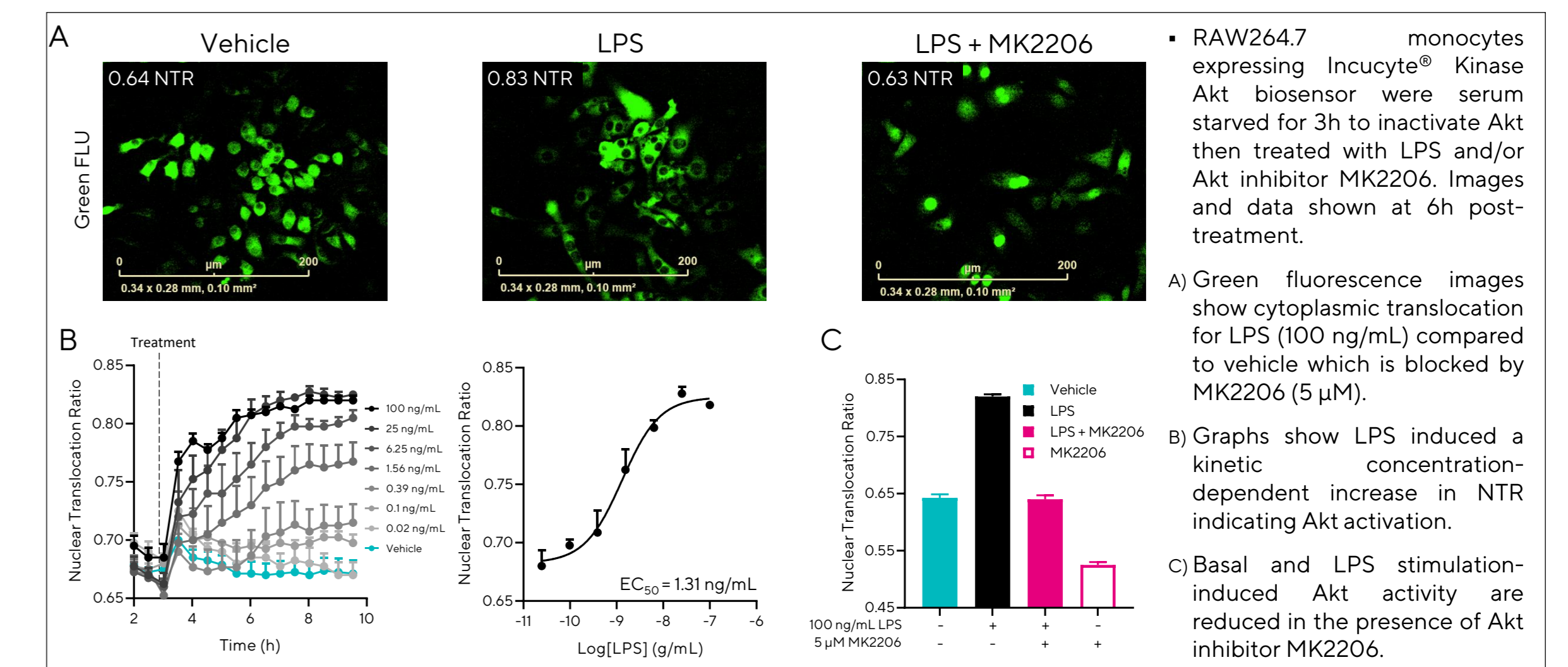
Incucyte[®] Live-Cell Imaging and Analysis Solutions

Incucyte[®] Live-Cell Analysis System
A fully automated HD phase-contrast and multi-color fluorescence system that resides within a standard cell incubator for optimal cell viability. Designed to scan plates and flasks repeatedly over time.

Incucyte[®] Software
Fast, flexible and powerful control hub for continuous live-cell analysis comprising image acquisition, processing and data visualization.

Incucyte[®] Reagents and Consumables
A suite of non-perturbing cell labelling and reporter reagents. Includes nuclear-targeted fluorescent proteins for cell counting plus no-wash cell health reagents for apoptosis, cytotoxicity, and many more.

Activation of Akt in LPS-stimulated immune cells



Inflammatory effect of cytokines on Akt activity

