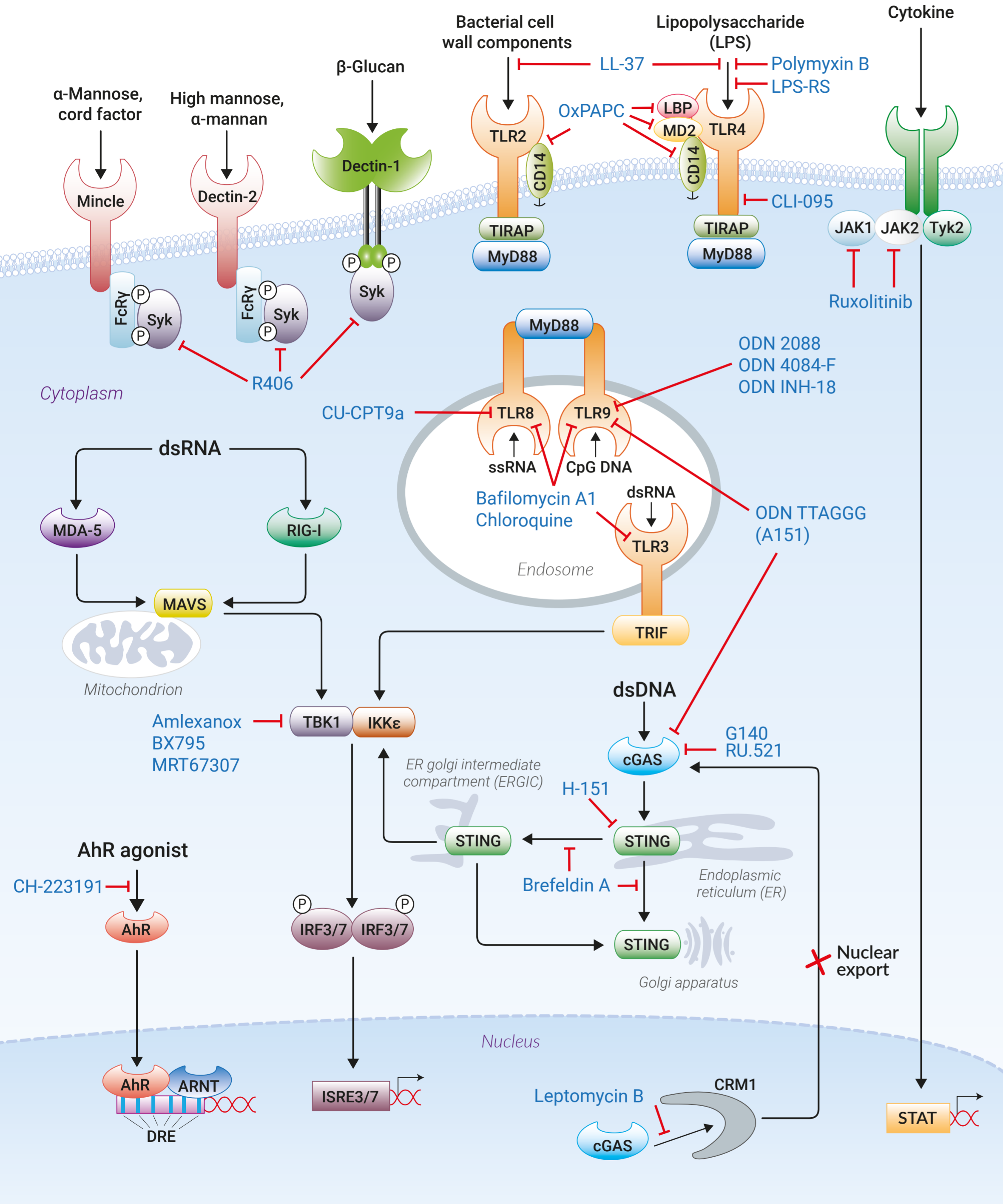
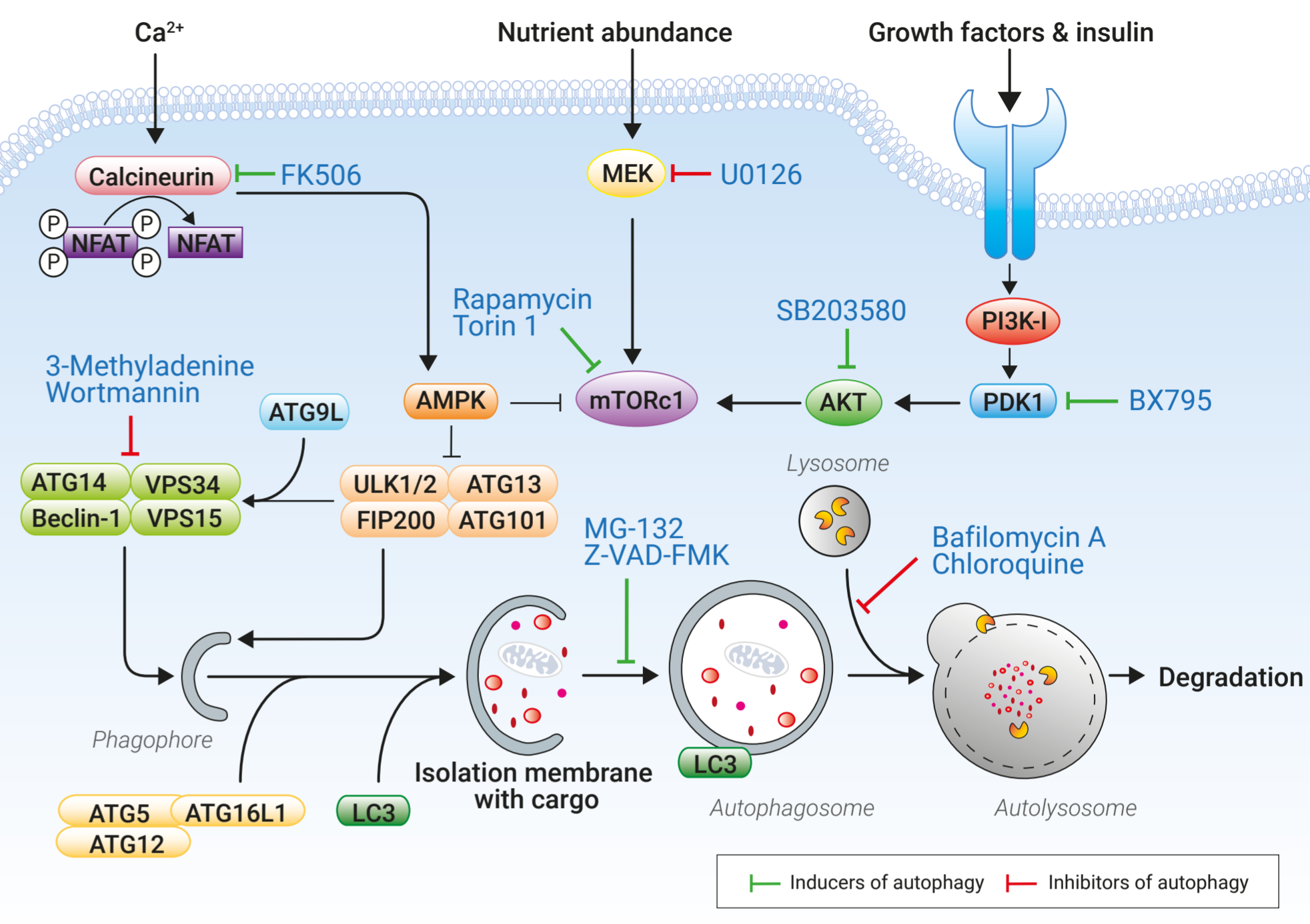


# TLR, RLR, CLR & AhR Inhibitors



# Autophagy Modulators



# Inflammasome & Regulated Cell Death Modulators

