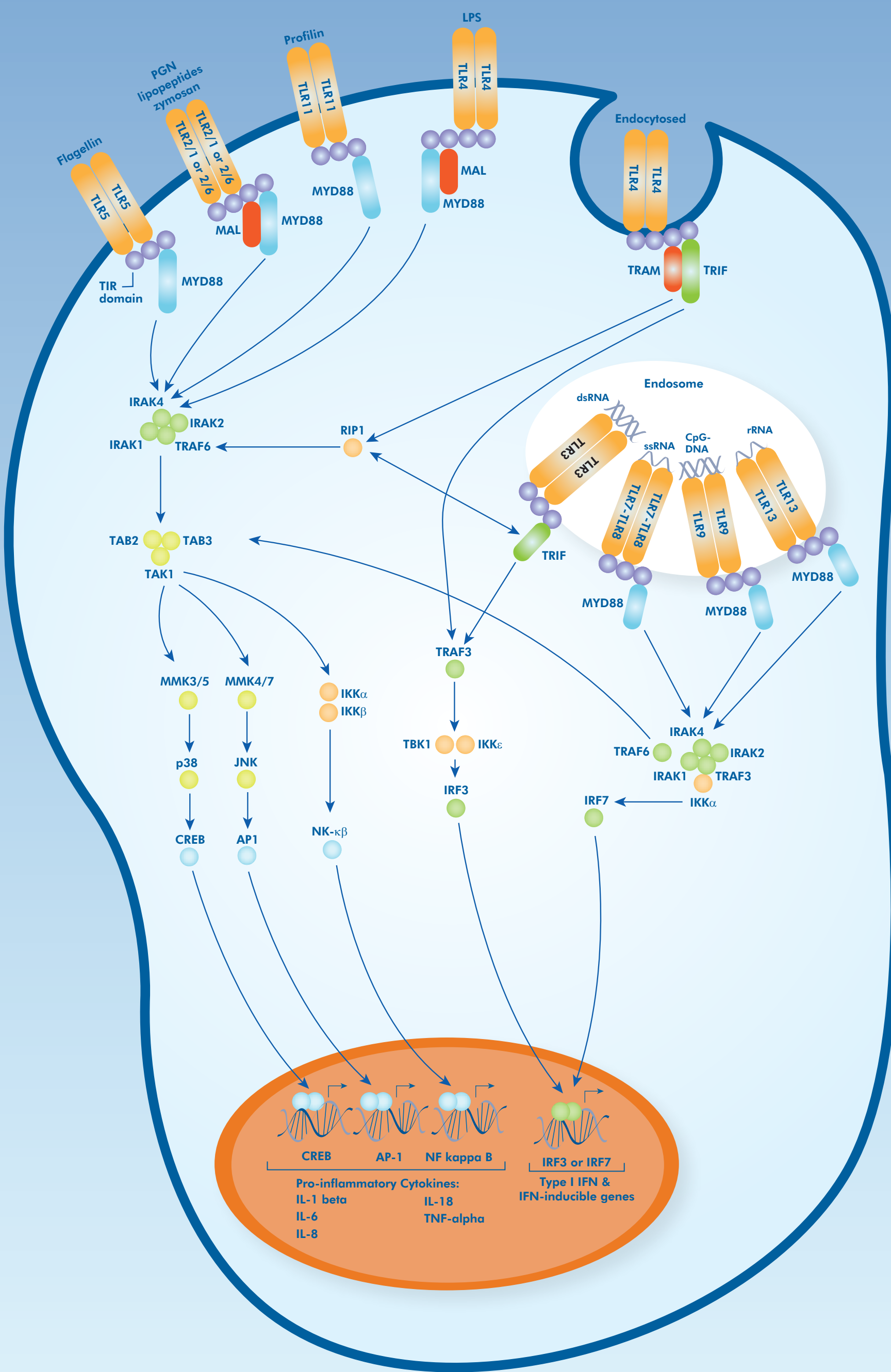
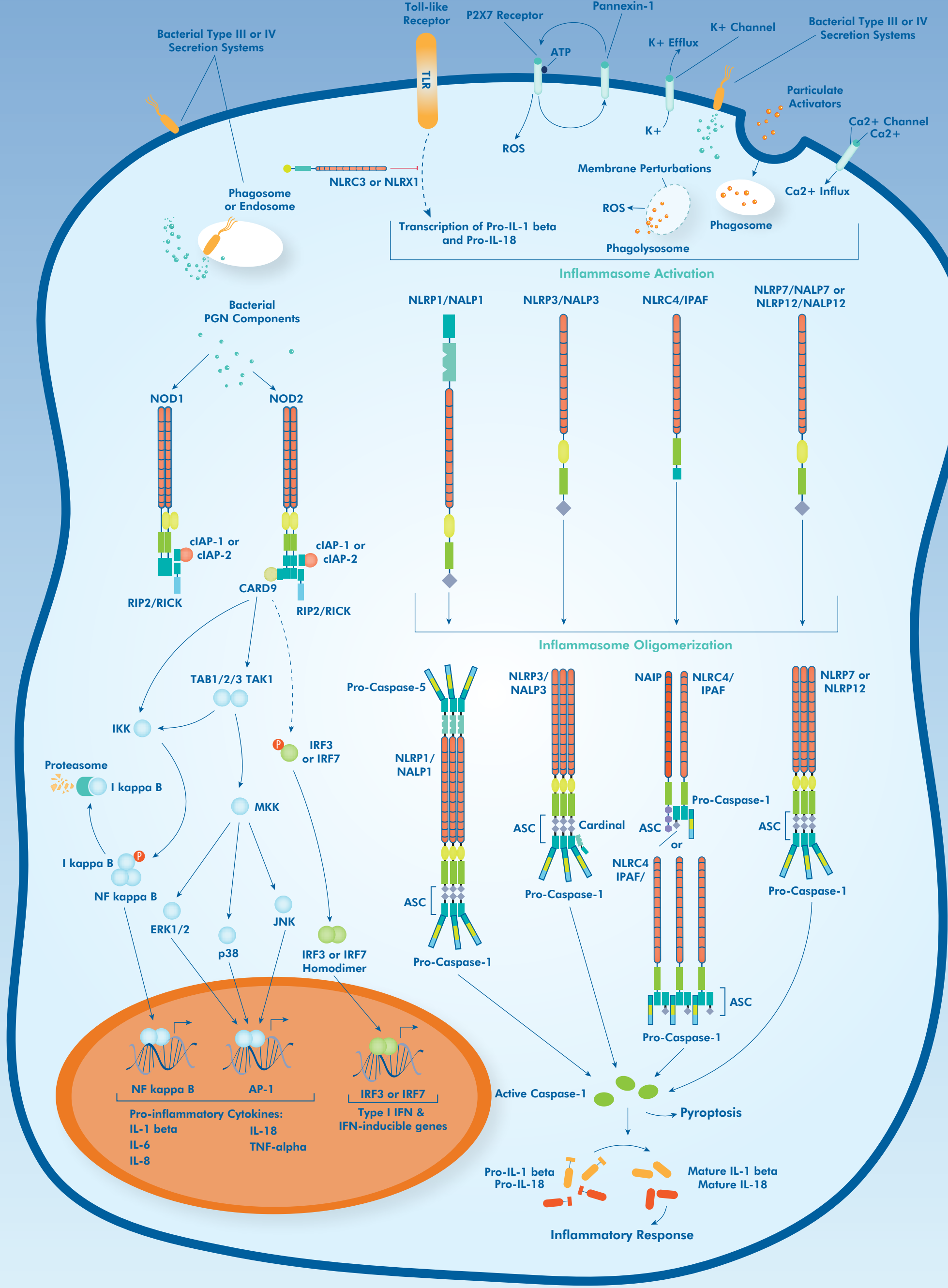


TOLL-LIKE RECEPTOR (TLRs)



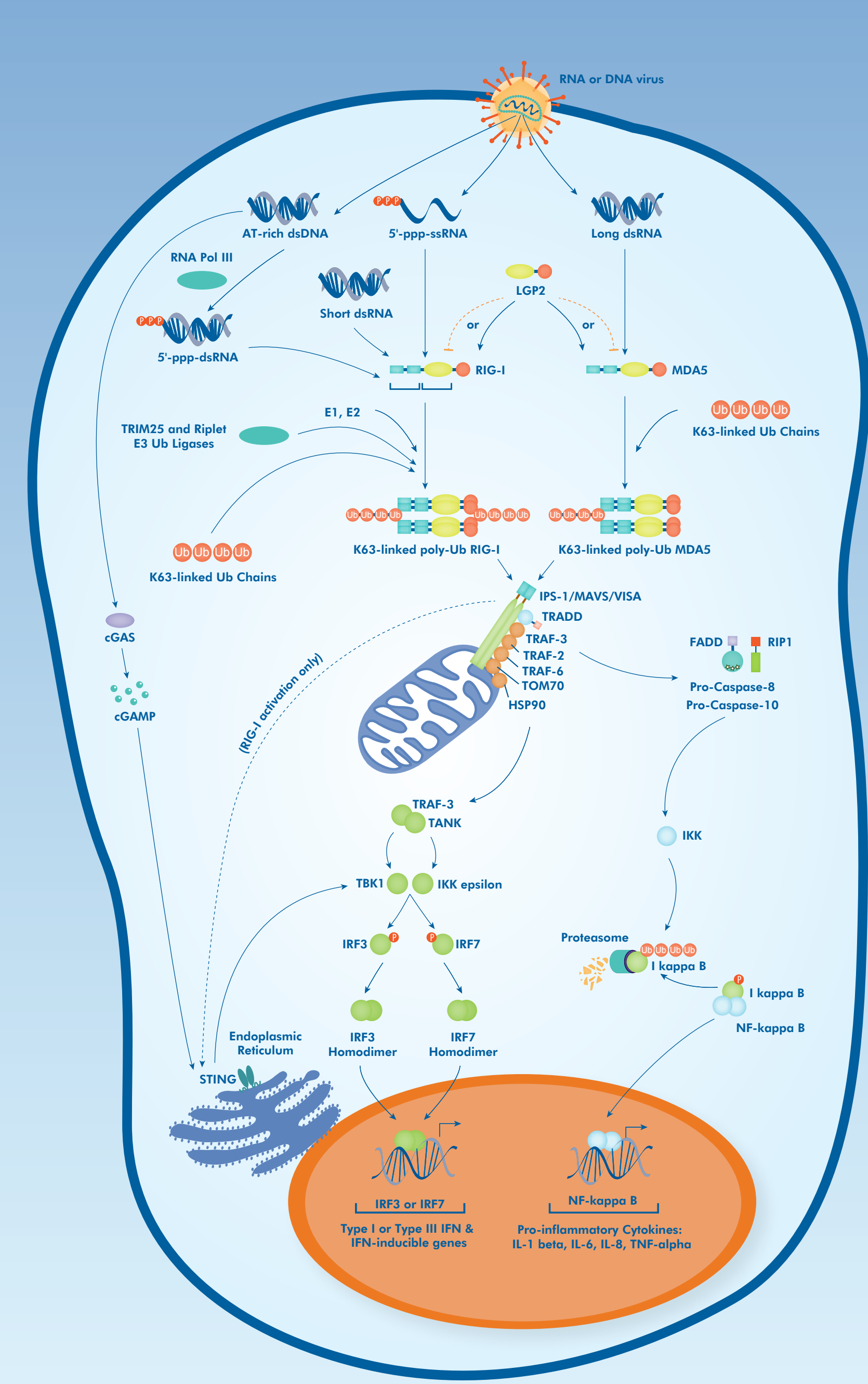
Domain Key
 Toll-Interleukin-1 Receptor Domain (TIR)
 MyD88-adaptor-like protein (MAL)
 TIR domain-containing adaptor-inducing IFN-β (TRIF)
 TRIF-related adaptor molecule (TRAM)

NOD-LIKE RECEPTORS (NLRs)



Domain Key
 Caspase Recruitment Domain (CARD)
 Unknown N-terminal Domain
 Pyrin Domain (PYD)
 Nucleotide-binding/Oligomerization Domain (NACHT or NOD)
 NACHT-Associated Domain (NAD)
 Function to Find Domain (FIIND)
 Leucine-Rich Repeat Domain (LRR)

RIG-I-LIKE RECEPTORS (RLRs)



Domain Key
 Caspase Recruitment Domain (CARD)
 RNA Helicase Domain
 Repressor Domain
 Death Domain (DD)
 Death Effector Domain (DED)
 Kinase Domain