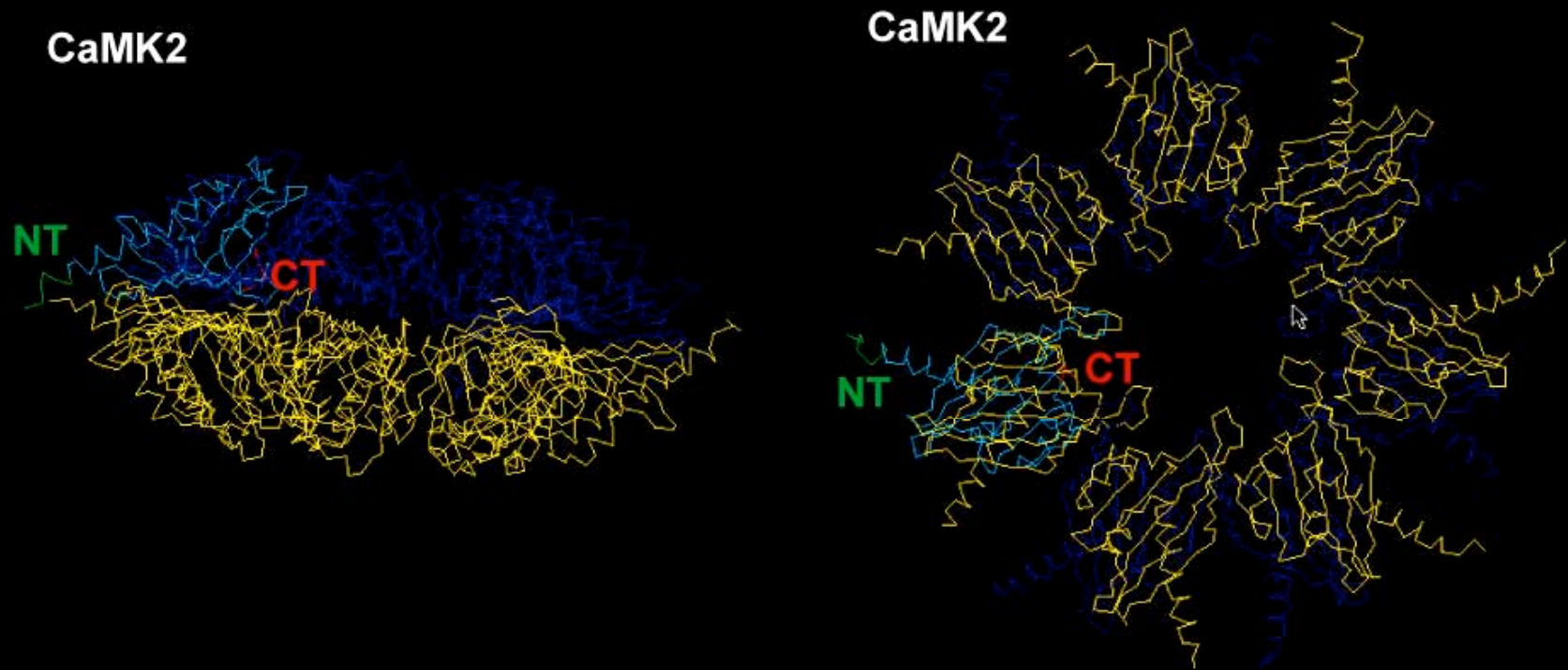


Stick Model Views of Calmodulin-Dependent Kinase 2 Holoenzyme Structure



KINEXUS



www.kinexus.ca