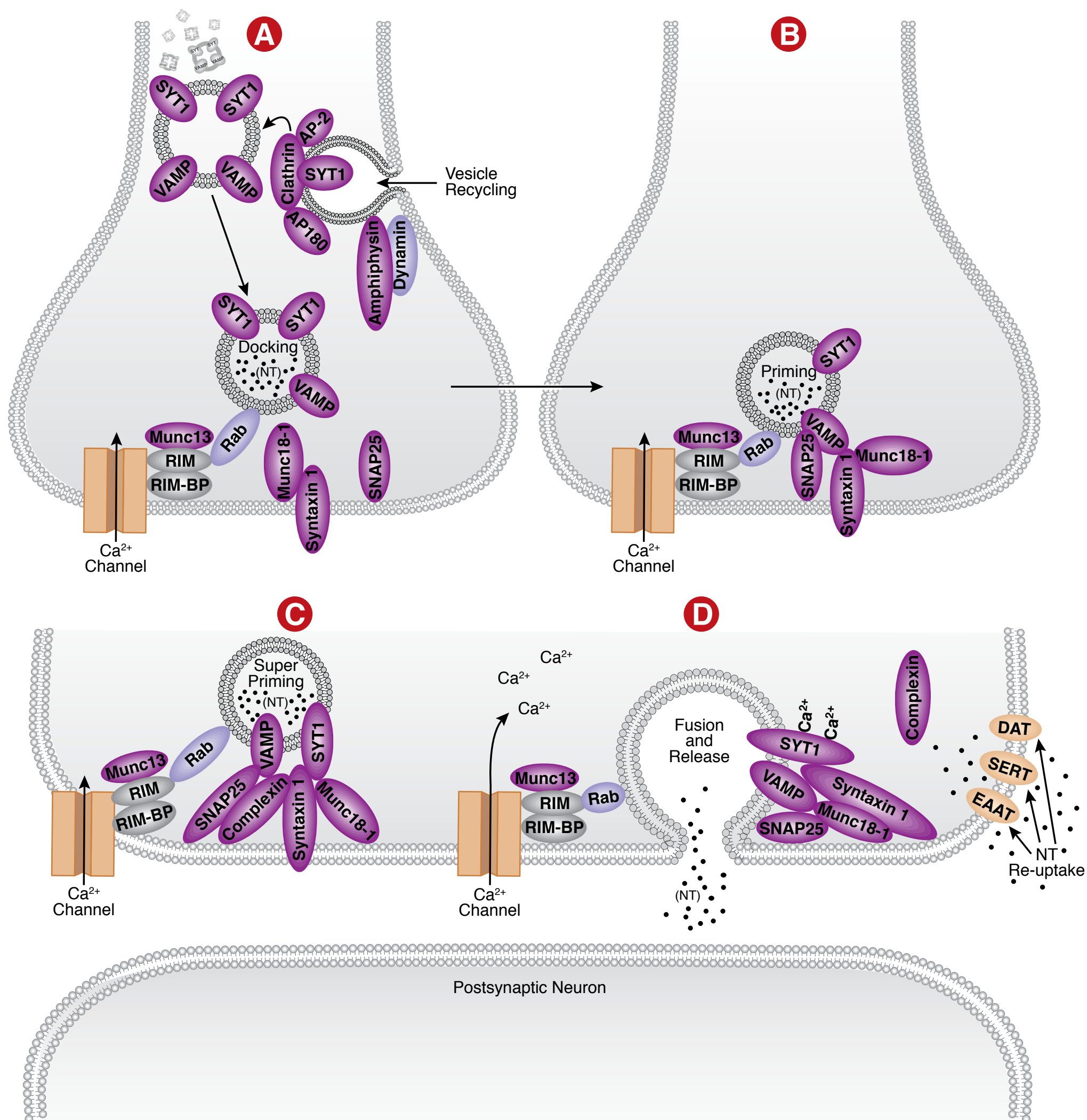


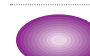


Vesicle Trafficking in Presynaptic Neurons: Synchronous Release



Pathway Diagram Key

-  Ion Channel/Transport
-  GTPase
-  Vesicle Formation/Trafficking

Select Reviews

This pathway diagram has been assembled by key opinion leaders in the field of Vesicle Trafficking in Presynaptic Neurons in collaboration with Cell Signaling Technology (CST) scientists. CST Pathways are reviewed and updated frequently to provide succinct and current overviews of the subject area. We would like to thank Prof. Taulant Bacaj, Stanford School of Medicine, Stanford, CA for contributing to this diagram.

Blakely, R.D. and Edwards, R.H. (2012) *Cold Spring Harb. Perspect. Biol.* 4, a005595. | Jahn, R. and Fasshauer, D. (2012) *Nature* 490, 201–207. | Saheki, Y. and De Camilli, P. (2012) *Cold Spring Harb. Perspect. Biol.* 4, a005645. | Südhof, T.C. (2013) *Neuron* 80, 675–690. | Südhof, T.C. (2013) *Nat. Med.* 19, 1227–1231. | Südhof, T.C. (2012) *Neuron* 75, 11–25.