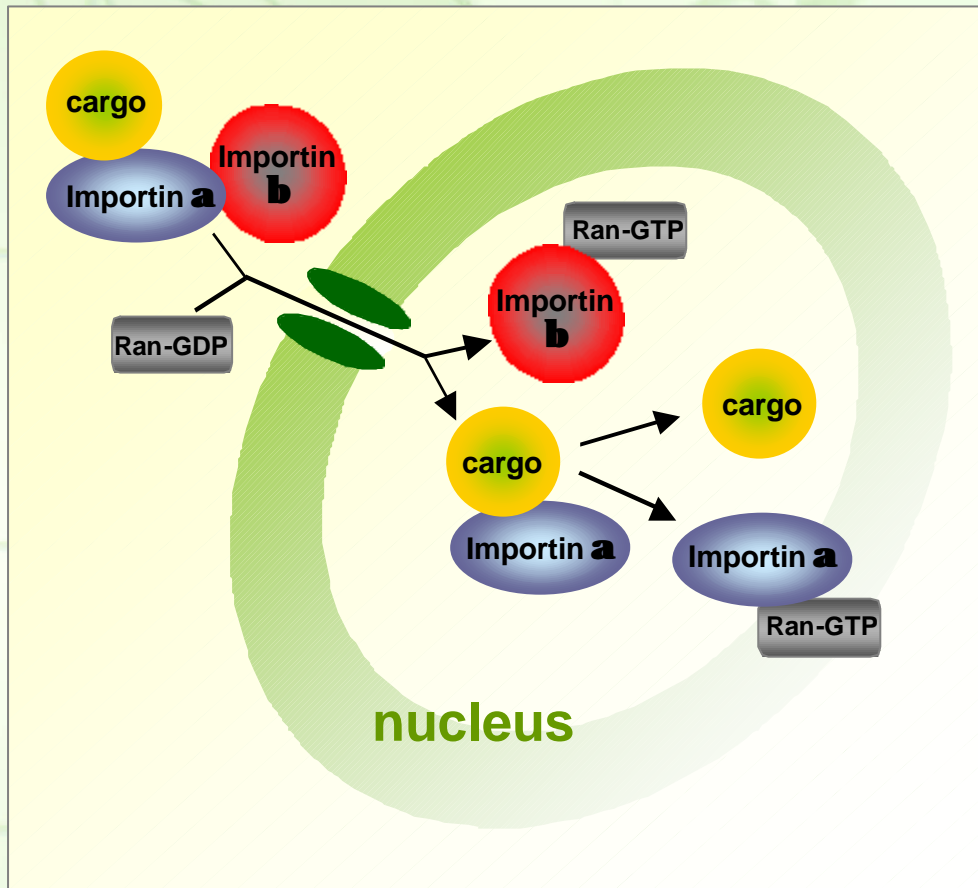


Importin **a** and **b** - Nuclear import factors



Simplified illustration of the transport mechanism of importin **a** and **b** into the nucleus

Importin **a** and **b**

- ▶ essential for nuclear protein import
- ▶ belong to a separate class of Ran-GTP binding proteins.
- ▶ Importin **a** binds directly to the nuclear localization signal (NLS) of substrates
- ▶ Importin **b** binds the **a**-NLS complex or other adaptors to mediate the uptake through the nuclear pore.

Jena Bioscience GmbH offers active Importin **a** and **b** suitable for

- ▶ in vitro nuclear import assays
- ▶ nucleocytoplasmic transport pathways regulation under different cellular conditions
- ▶ characterization of the auto-inhibitory sequence of importin **a**
- ▶ interaction studies with importin **b** or the importin **a/b** heterodimer

References:

- Furuta M. et al. (2004) Heat-shock induced nuclear retention and recycling inhibition of importin alpha. *Genes Cells*. 9:429
- Freedman N.D. et al. (2004) Importin 7 and importin alpha/importin beta are nuclear import receptors for the glucocorticoid receptor. *Mol. Biol. Cell*. 15:2276
- Schwindling S.L. et al. (2004) Mutation of a CK2 phosphorylation site in cdc25C impairs importin alpha/beta binding and results in cytoplasmic retention. *Oncogene*. 23:4155