

UNIVERSITY OF COPENHAGEN

Faculty of Health and Medical Sciences



Structure, function and pharmacology of G protein-coupled receptors:  
State-of-the-art and future challenges

Hans Bräuner-Osborne

Molecular Pharmacology Group  
Department of Drug Design and Pharmacology  
[www.molpharm.net](http://www.molpharm.net)  
[hbo@sund.ku.dk](mailto:hbo@sund.ku.dk)



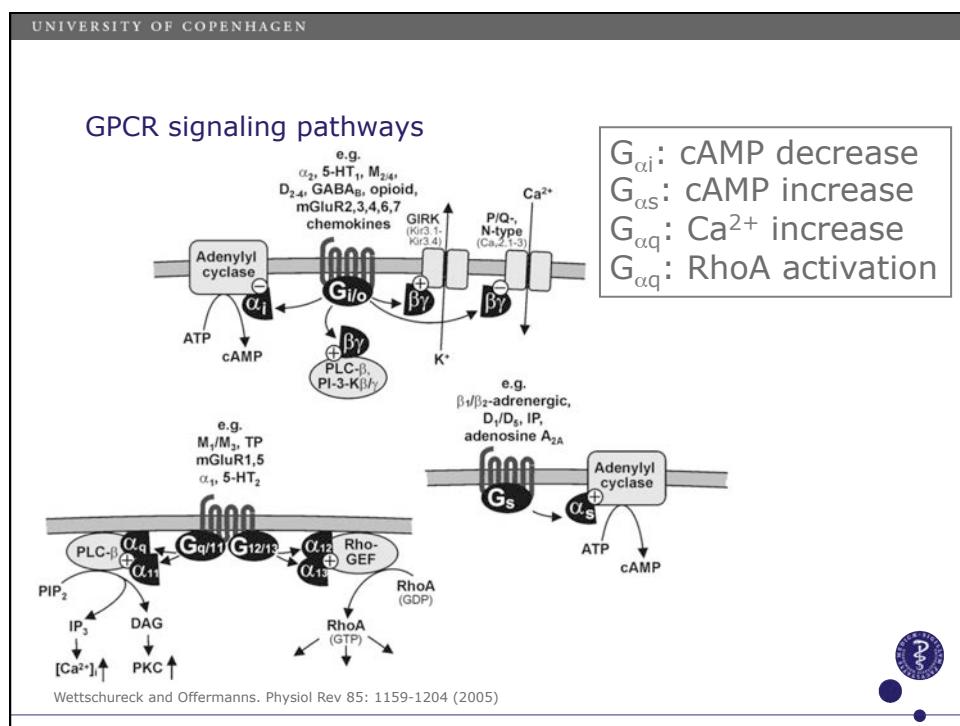
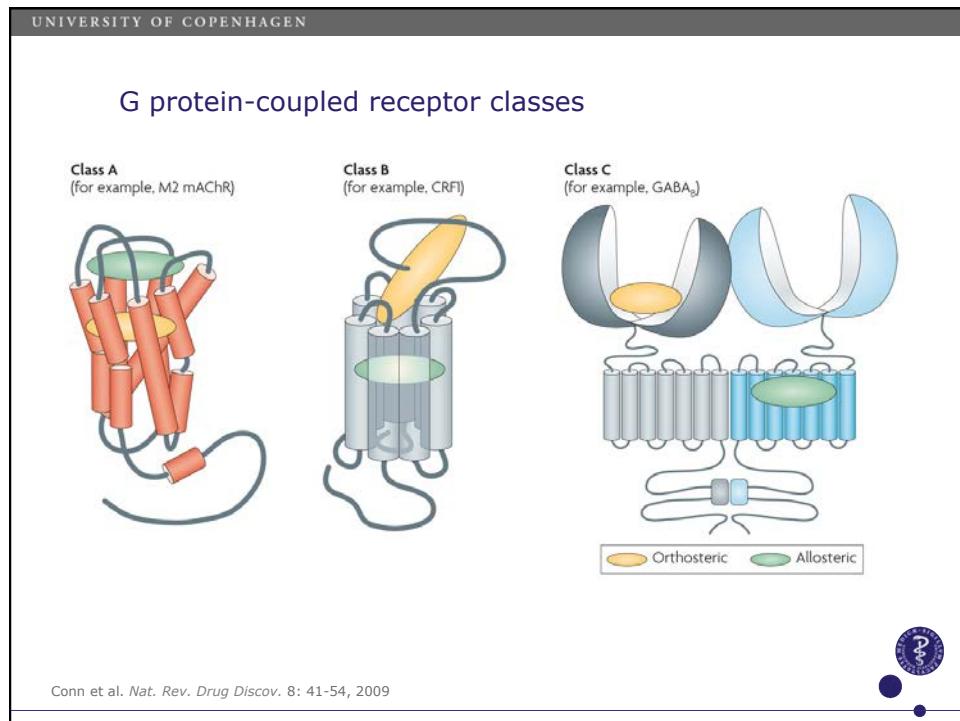
UNIVERSITY OF COPENHAGEN

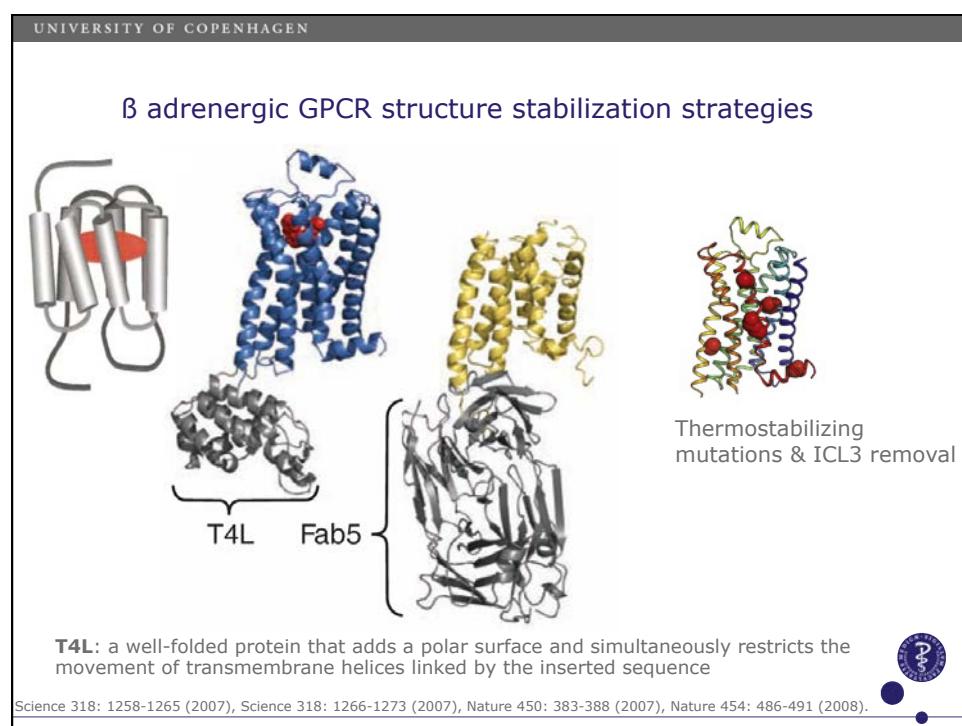
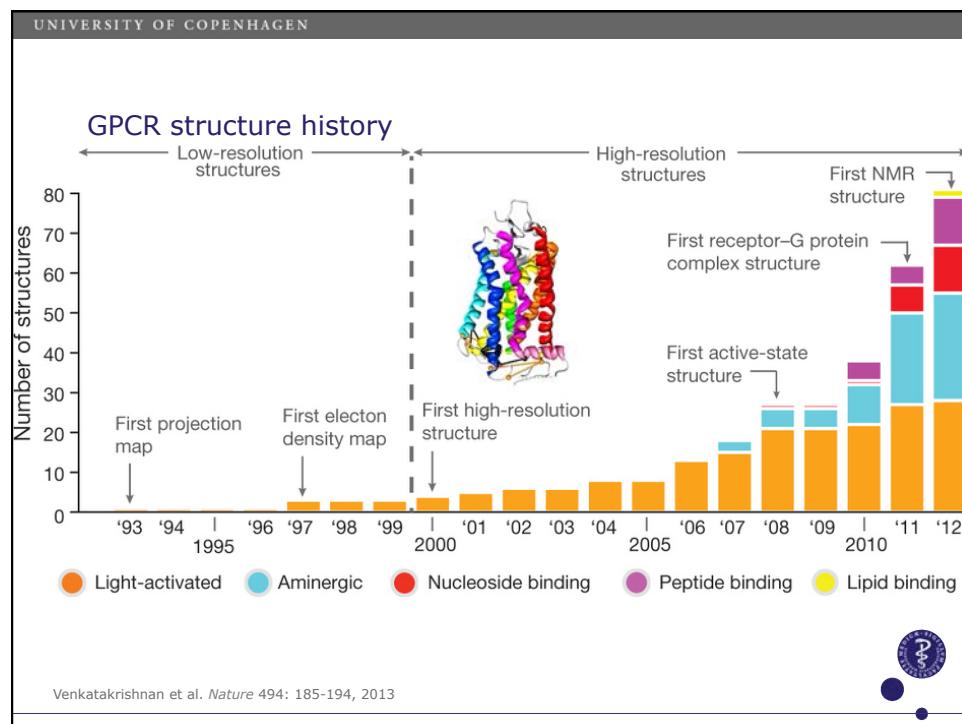
## Overview

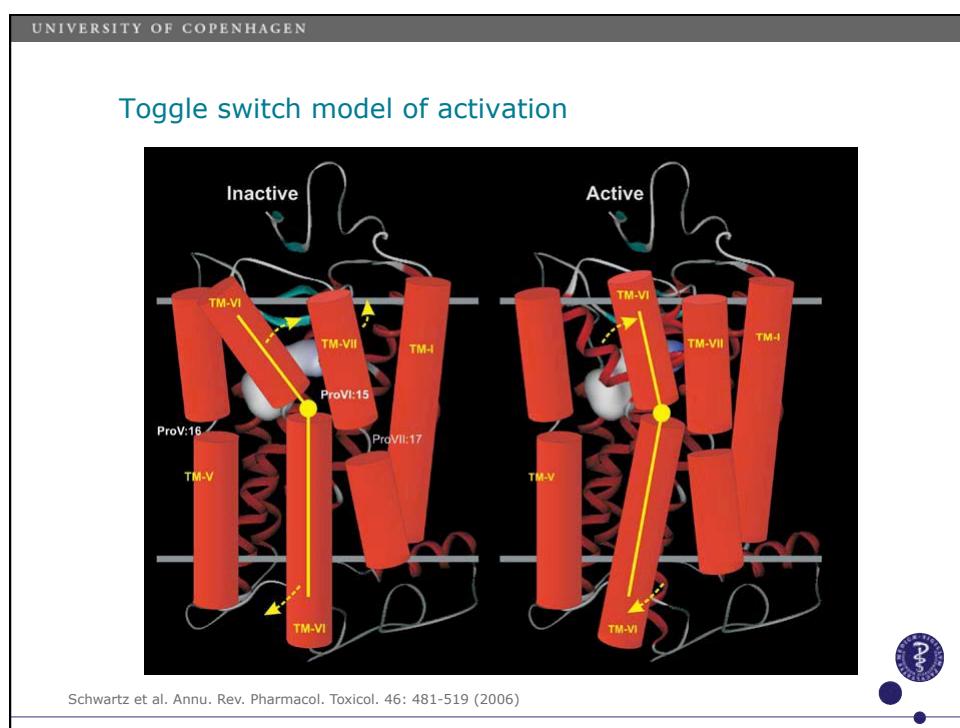
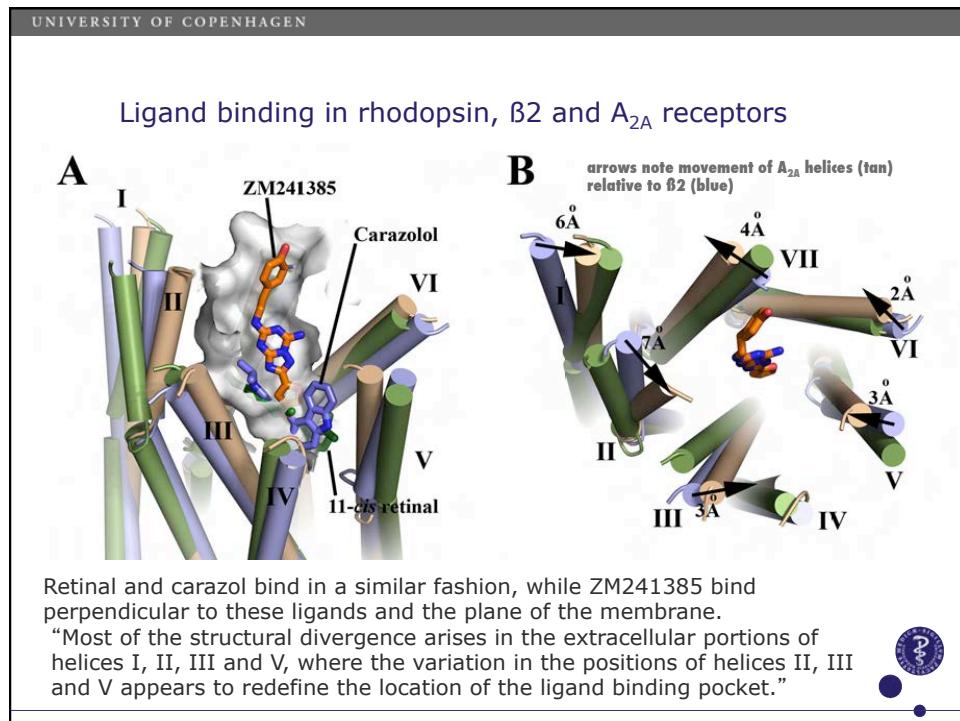
- GPCR structures
- Biased signaling
- Real-time receptor activation measurements
- Orphan receptors





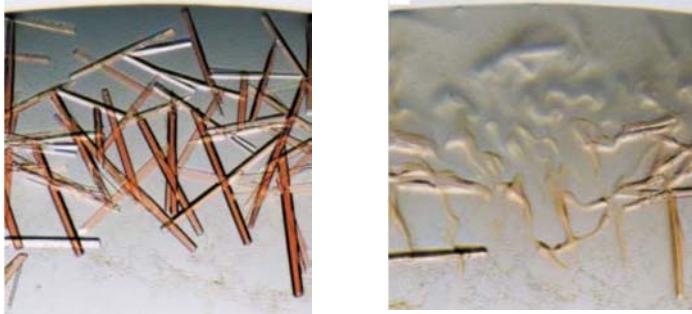






UNIVERSITY OF COPENHAGEN

GPCR structure - active conformation?



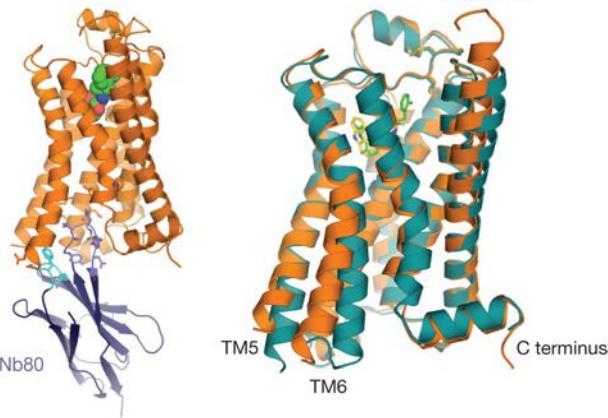
Crystals dissolve upon 1 min radiation with visible light.  
Indicates conformational/structural change.

Okada and Palczewski. Curr. Opin. Struct. Biol. 11: 420-426 (2001)



UNIVERSITY OF COPENHAGEN

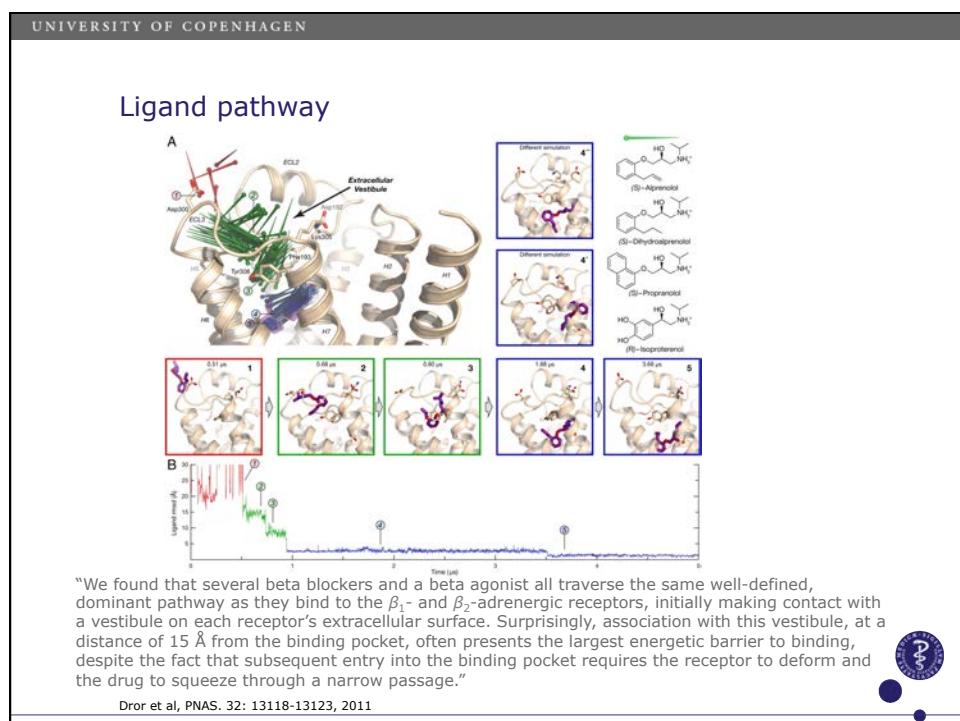
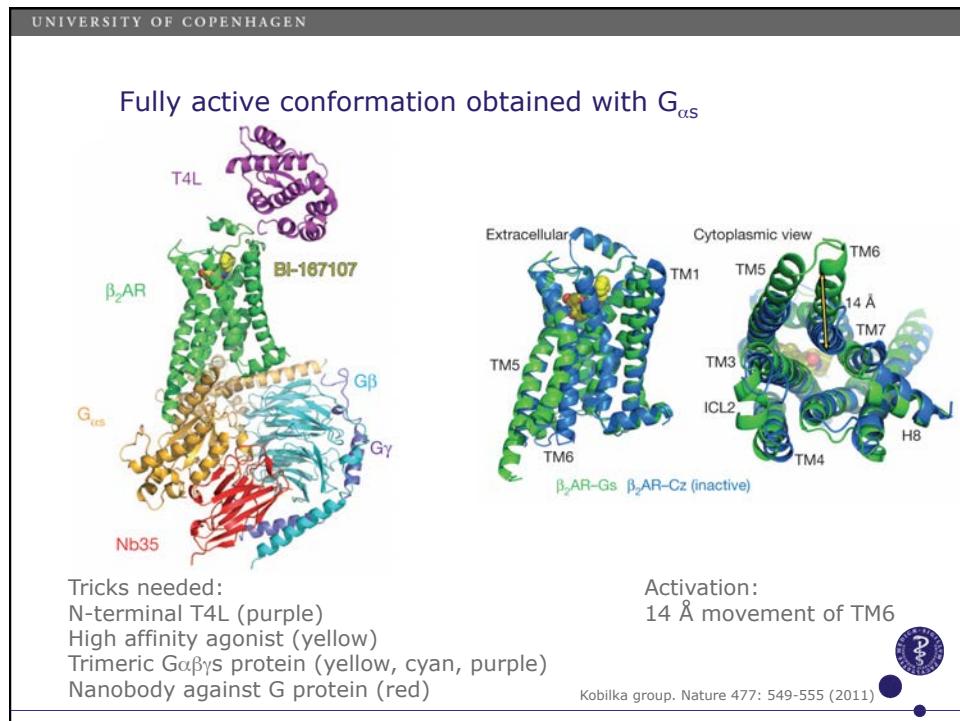
"Active" conformation by agonist and nanobody

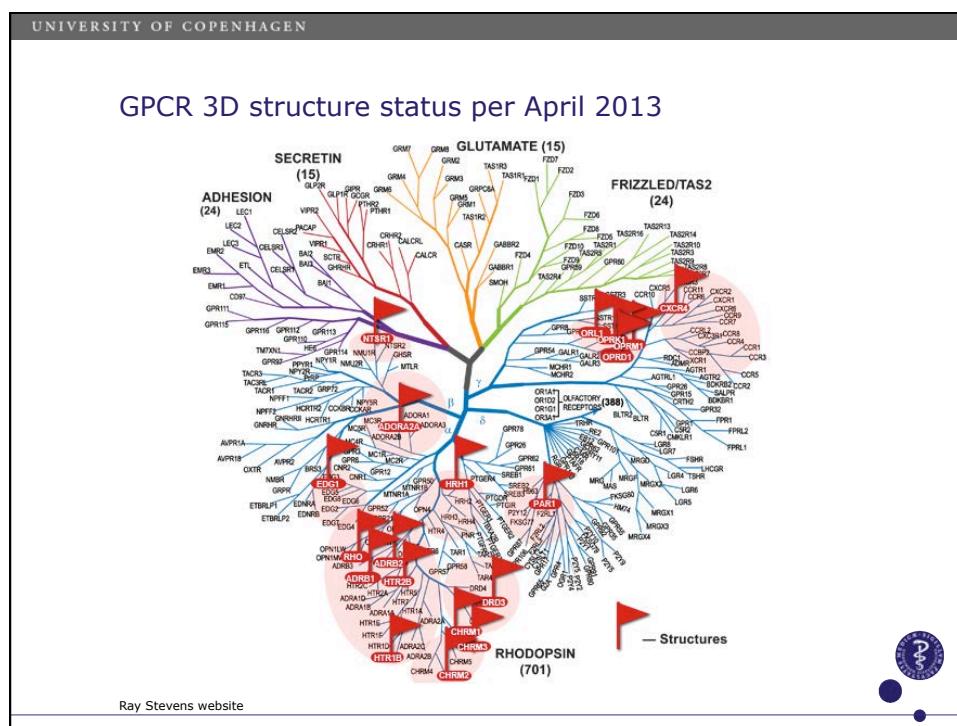
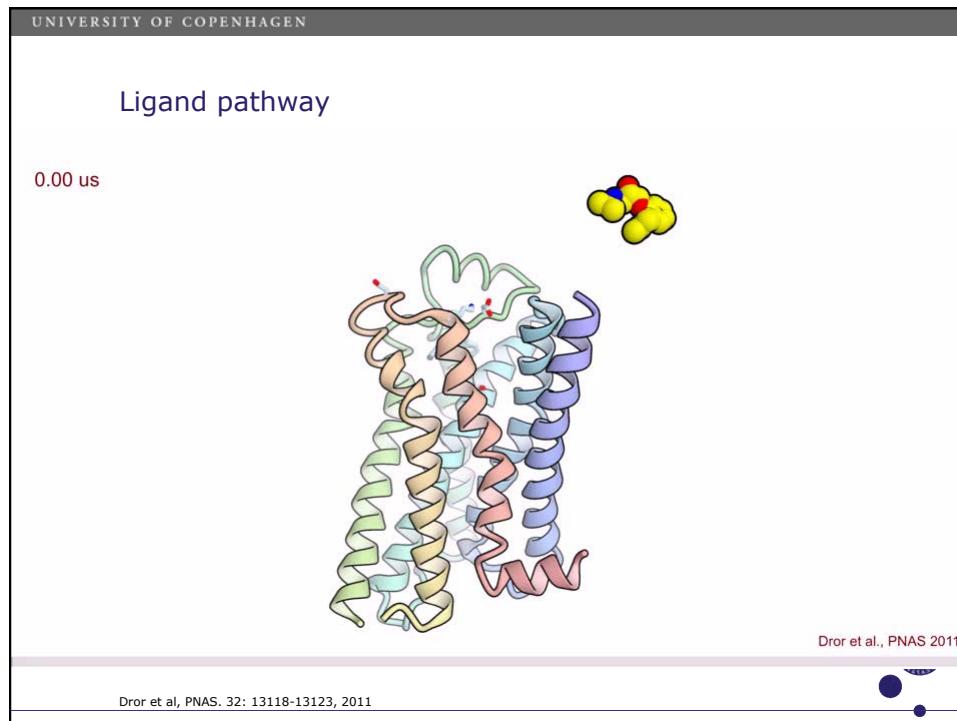


Orange: B2-T4L-Nanobody with full agonist (in green)  
Cyan: B2-T4L with inverse agonist (in yellow)  
Not possible to make agonist structure in the absence of nanobody.  
TM5 top moves 2 Å inward, TM3-5-6-7 hydrogen bond network rearranged.  
TM6 bottom moves 11 Å outward, TM5 & TM7 bottoms rearranged

Kobilka group. Nature 469: 175-180 (2011)



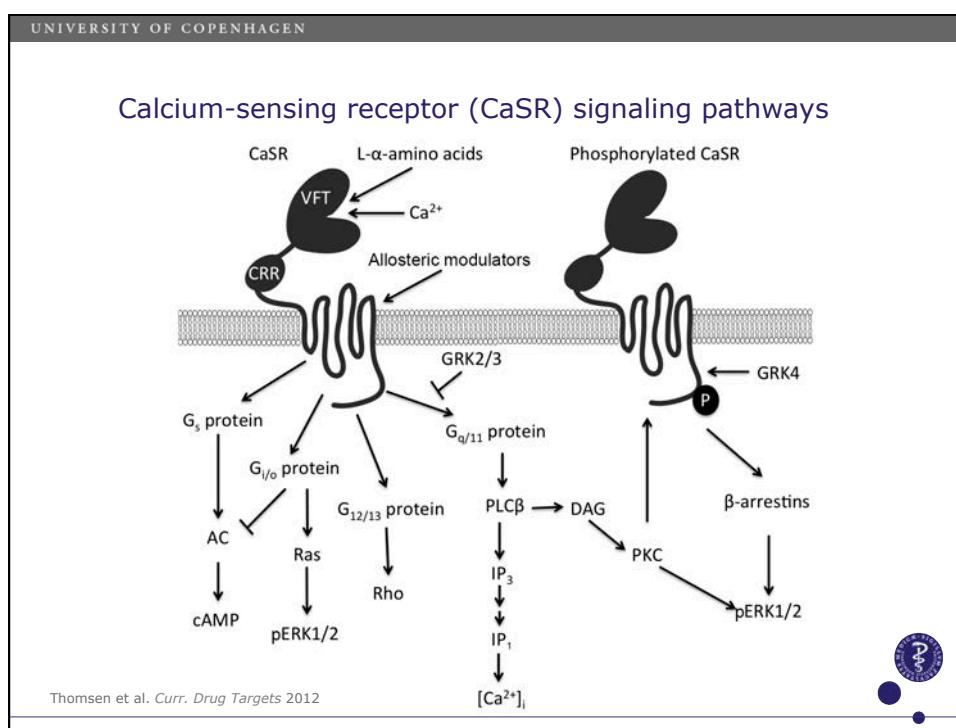


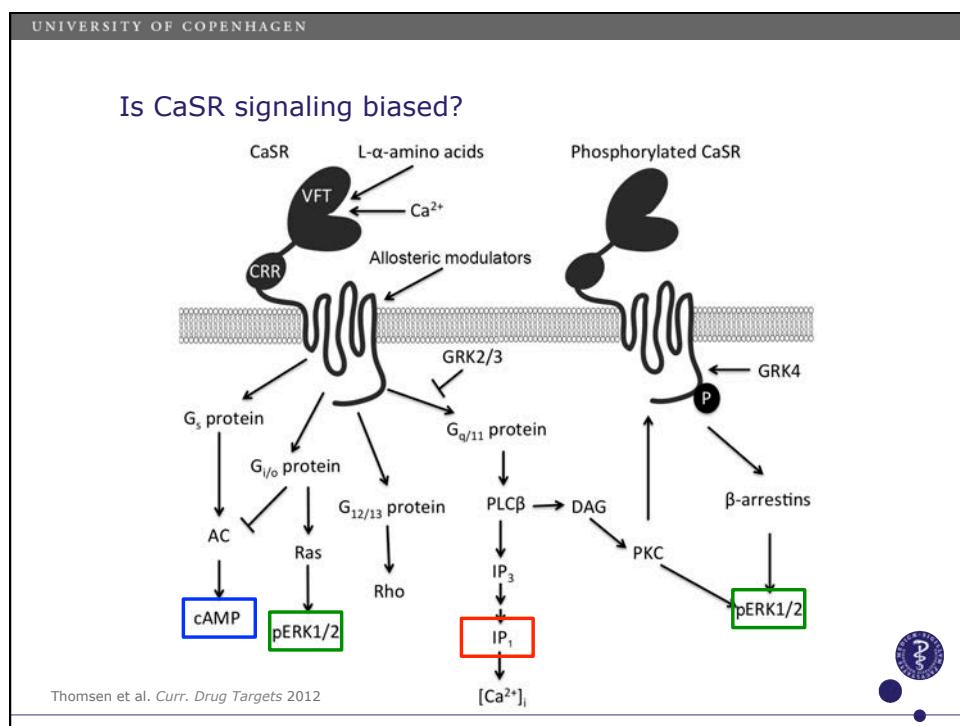
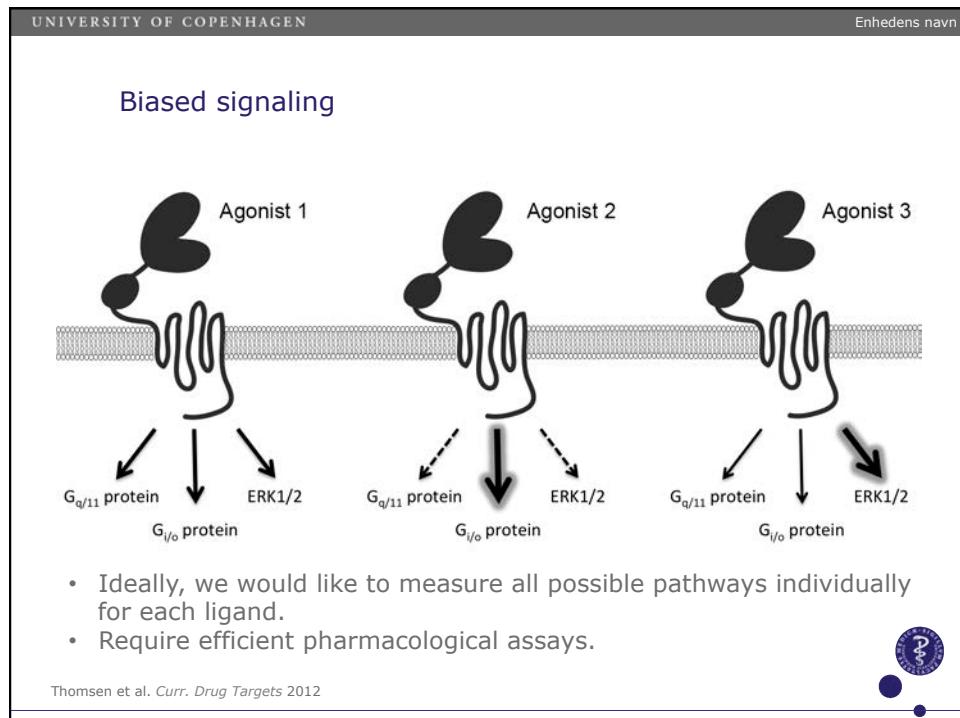


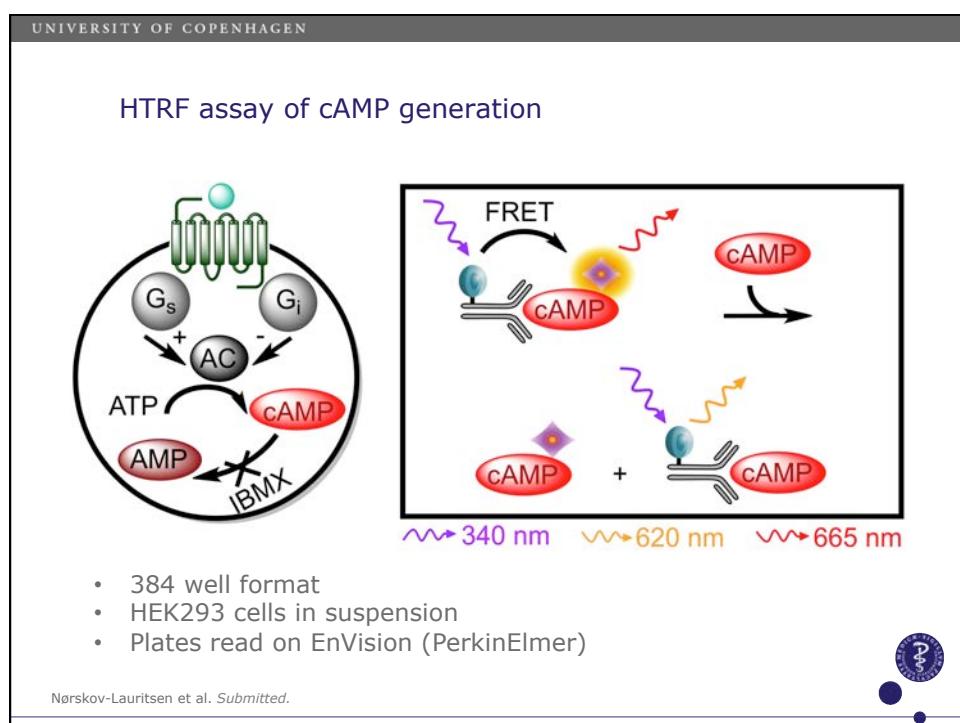
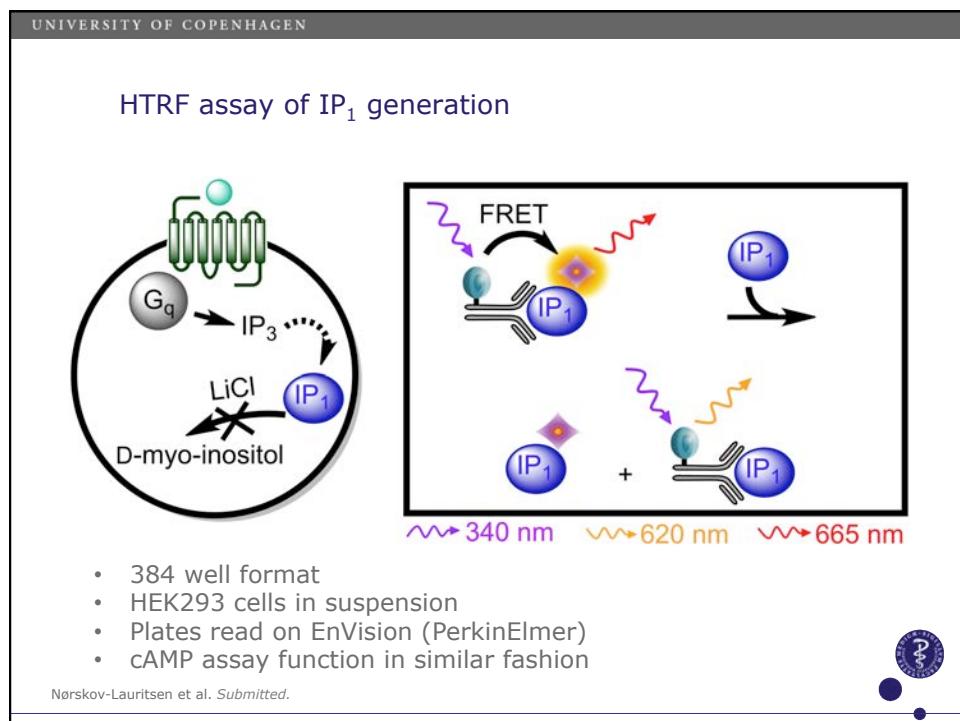
UNIVERSITY OF COPENHAGEN

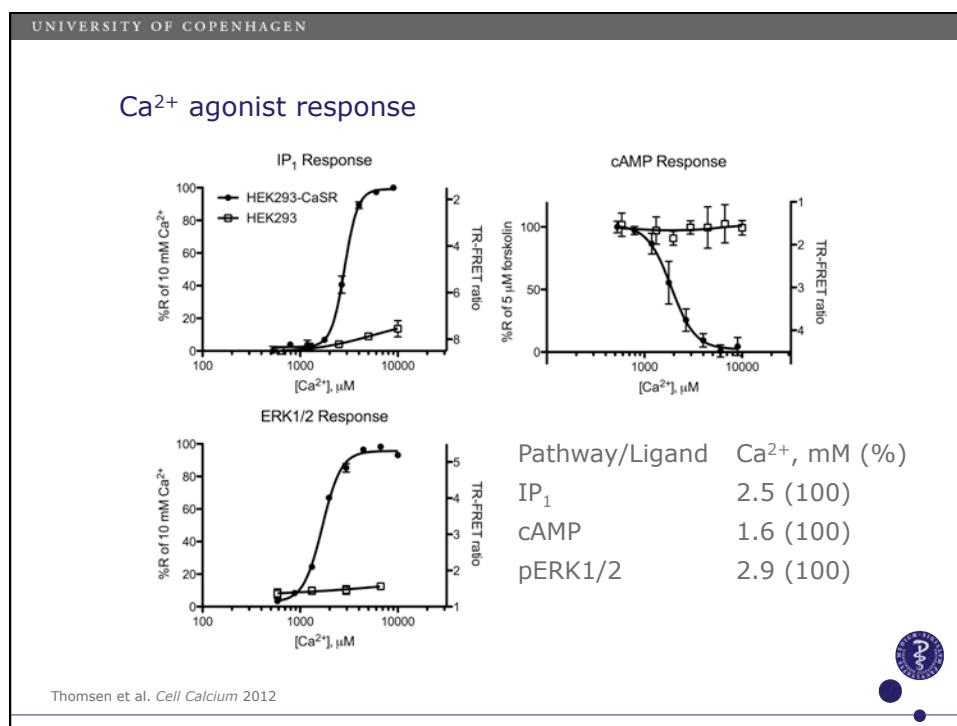
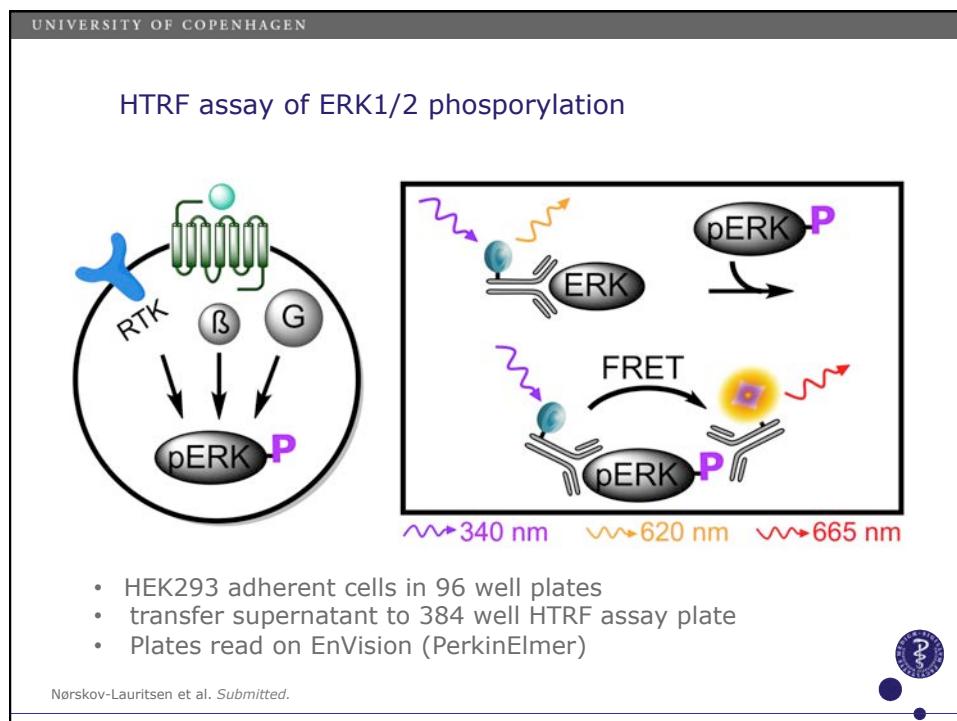
## Overview

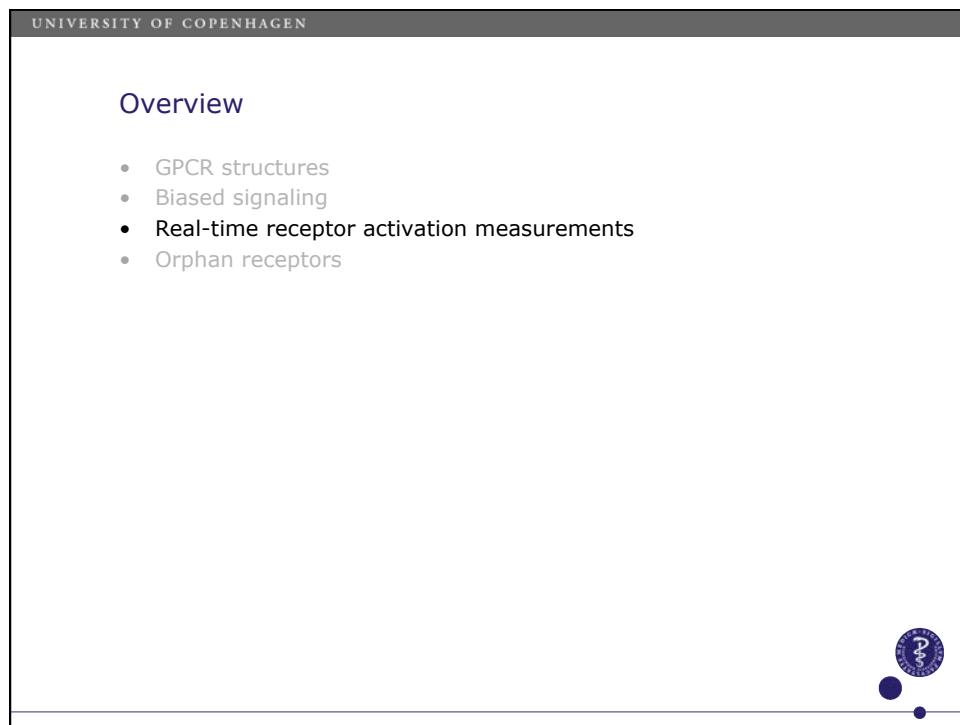
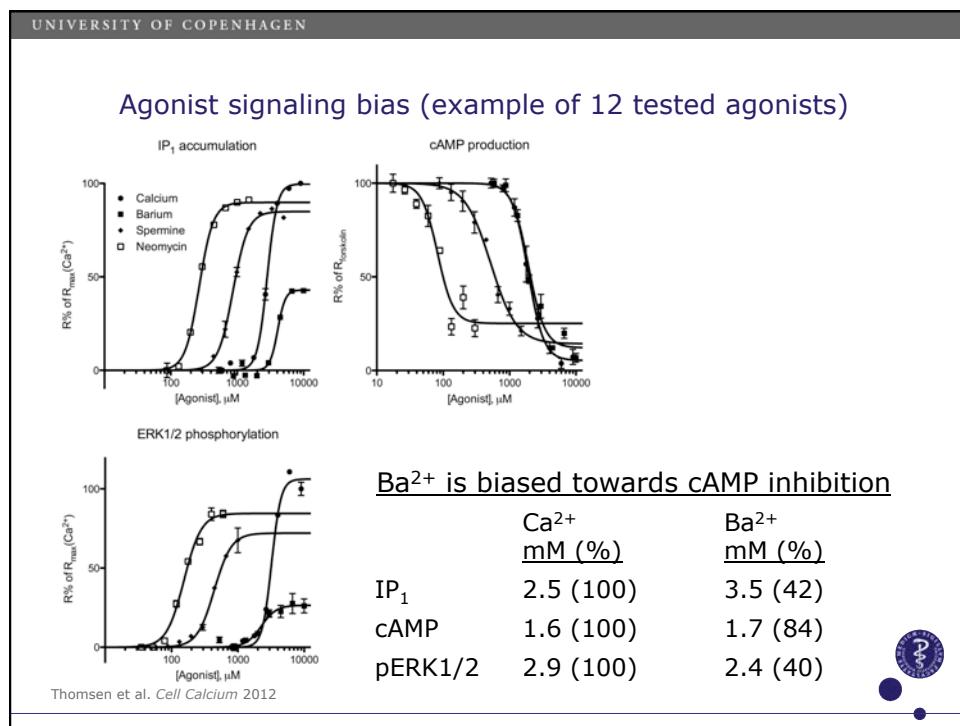
- GPCR structures
- Biased signaling
- Online receptor activation measurements
- Orphan receptors

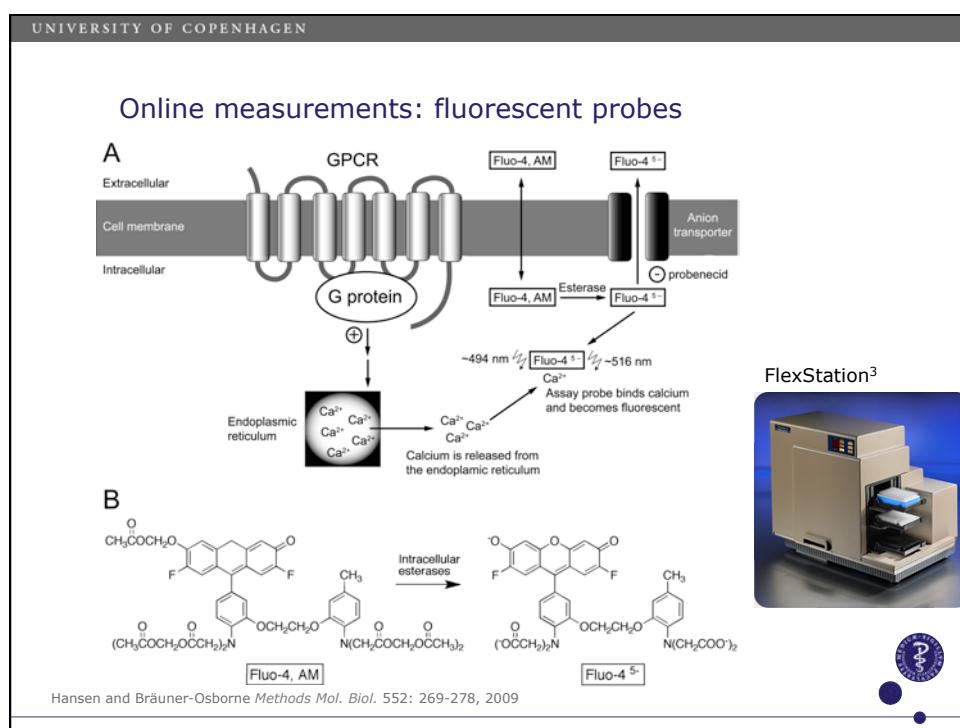
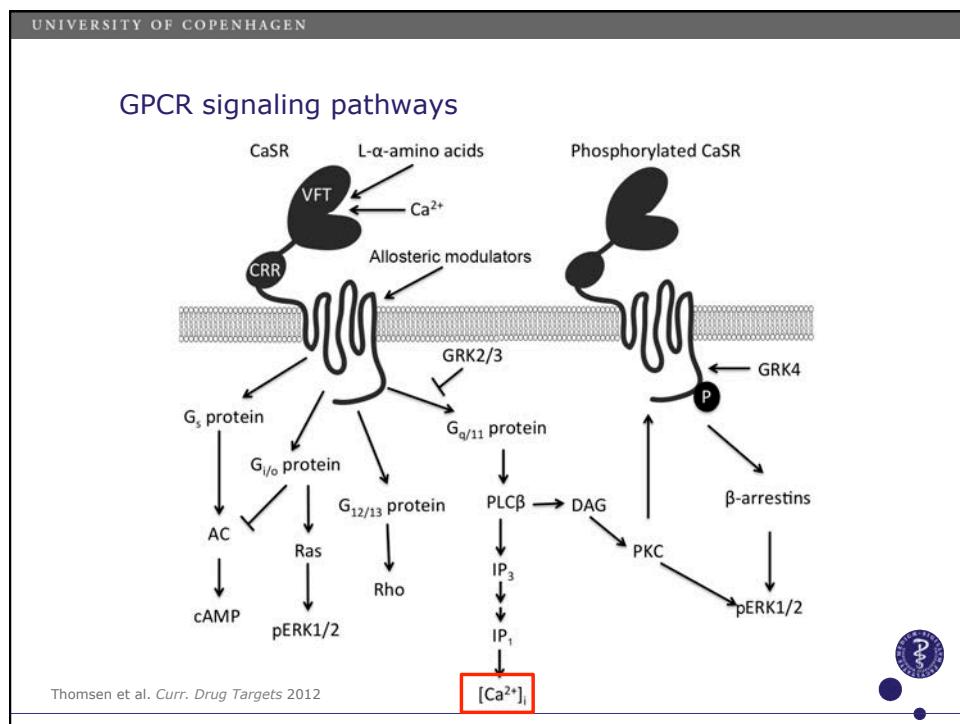



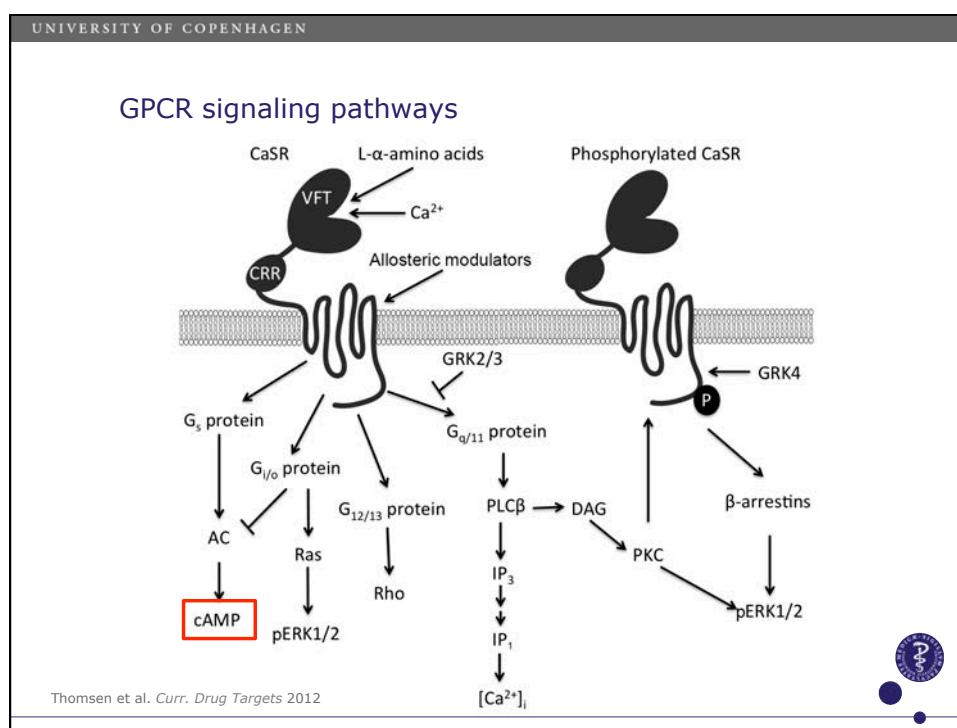
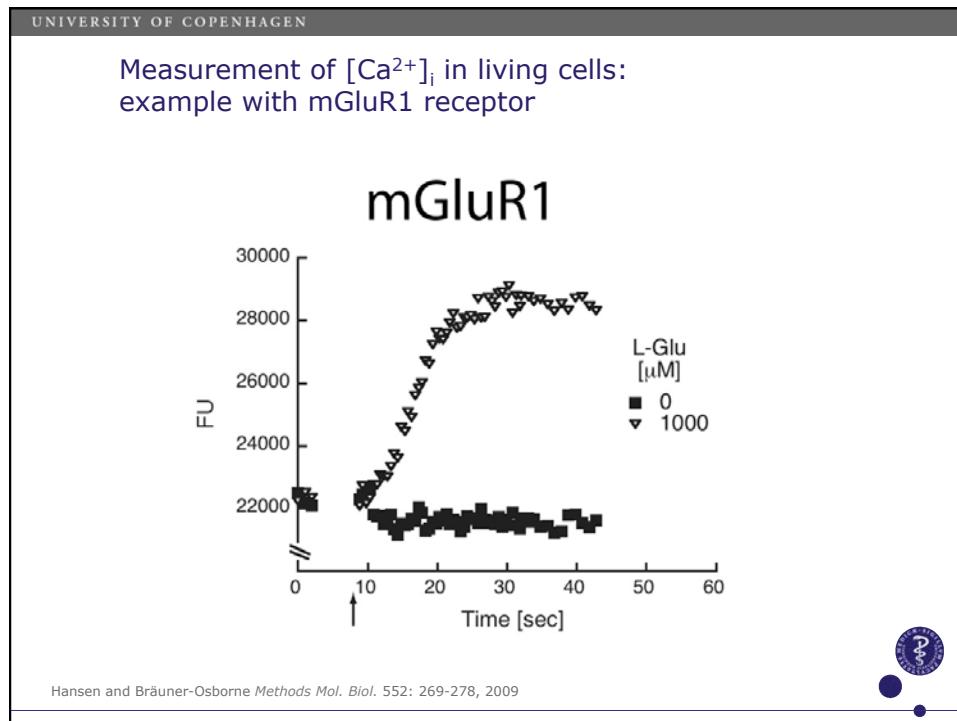


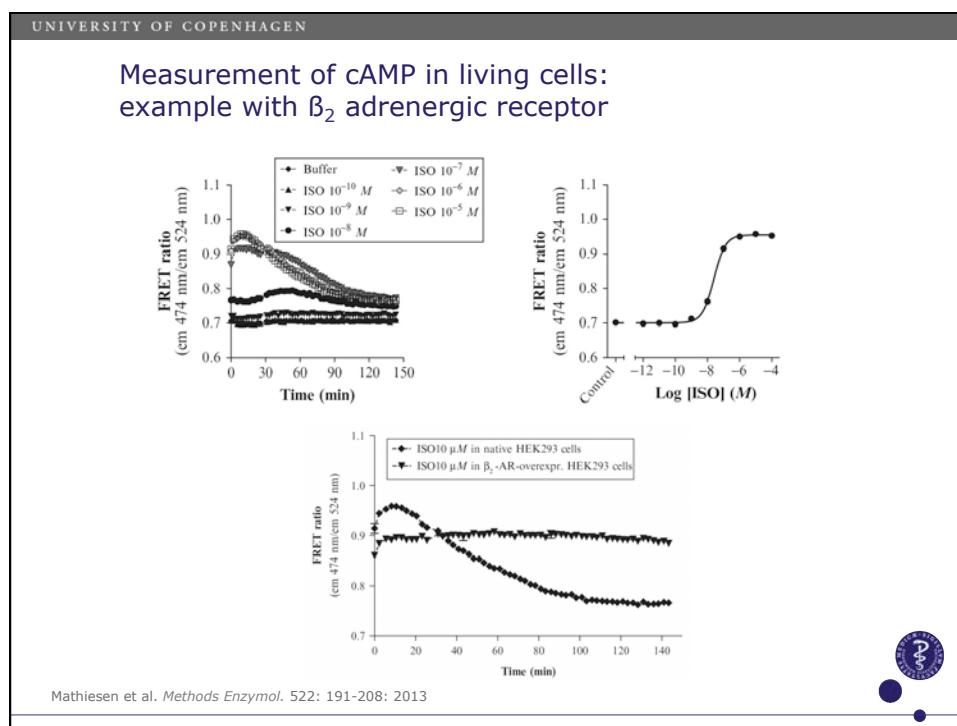
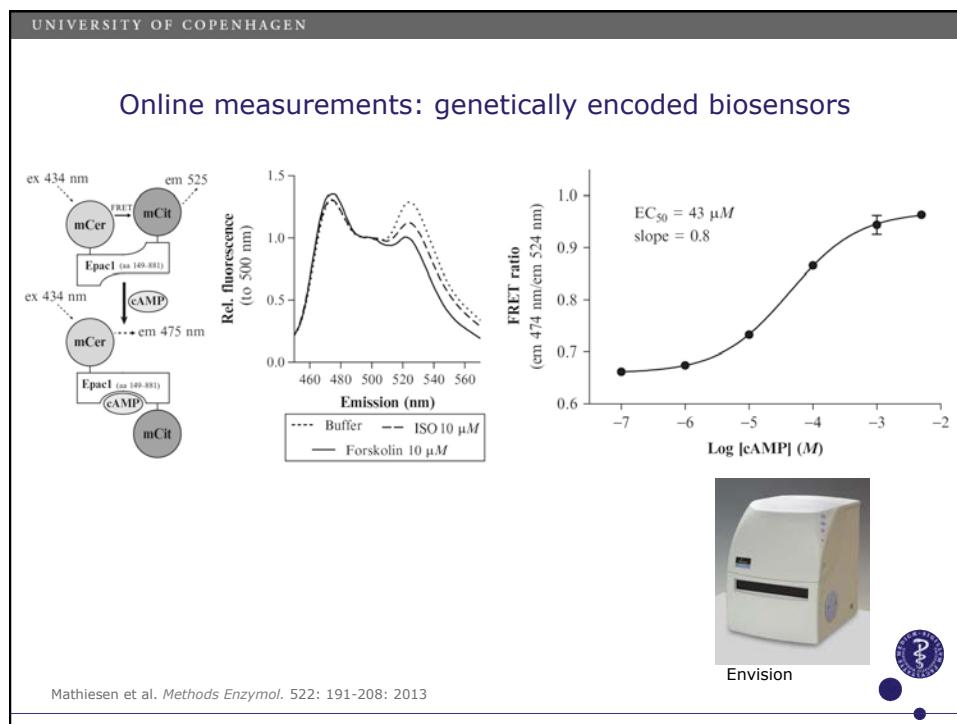


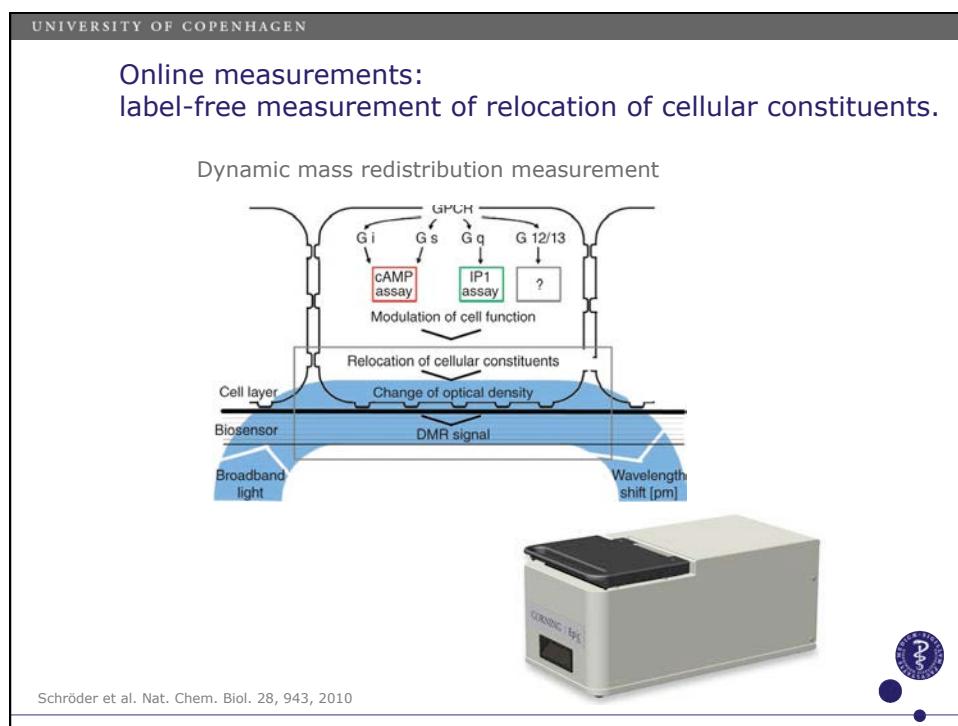
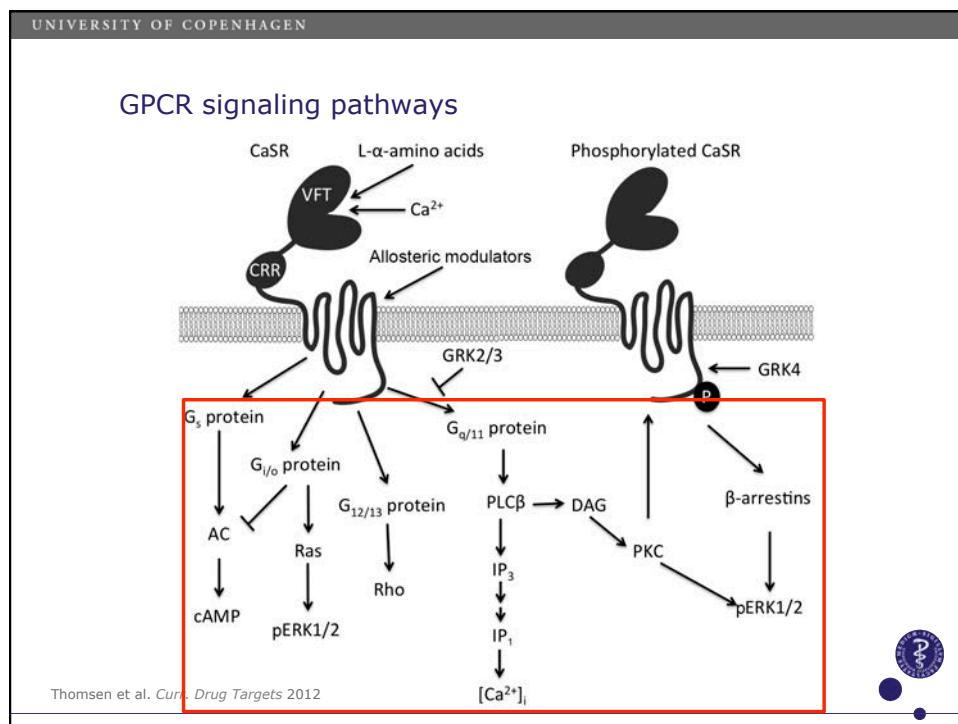


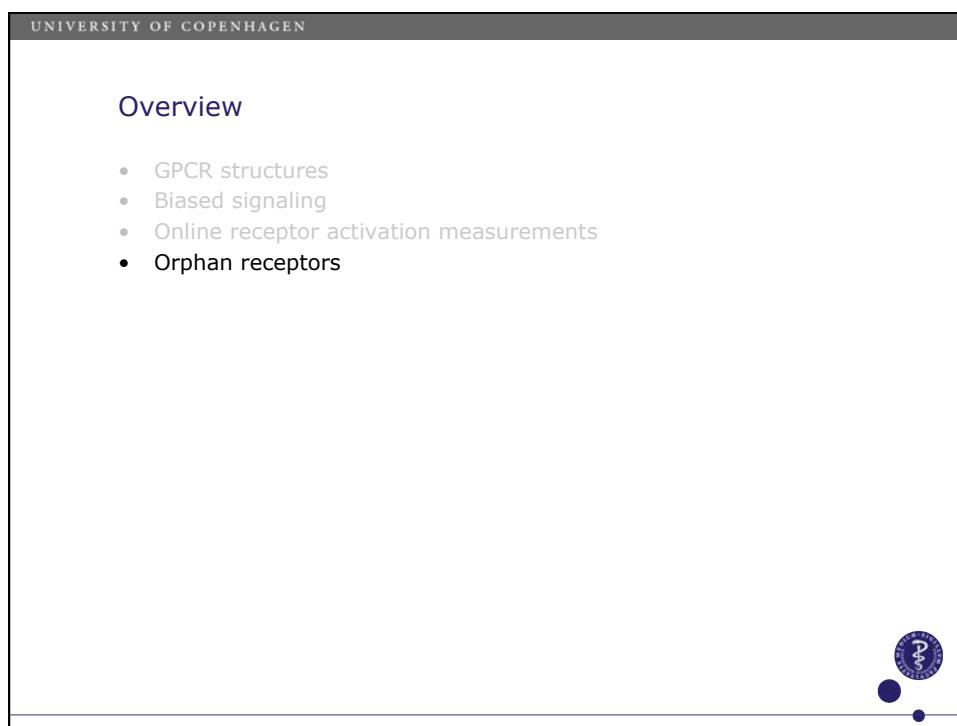
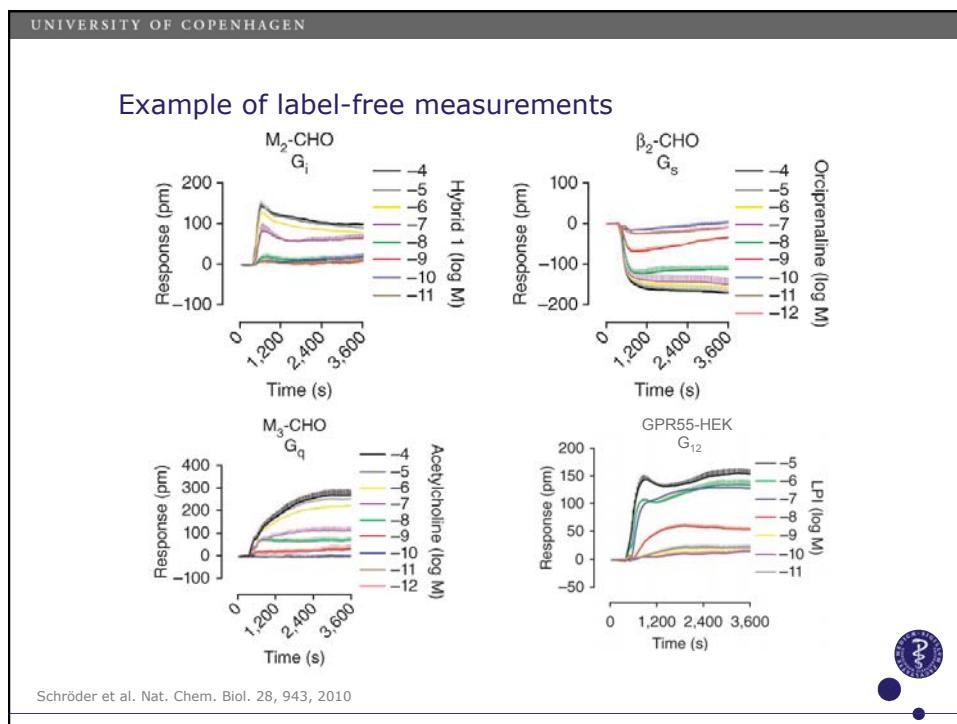


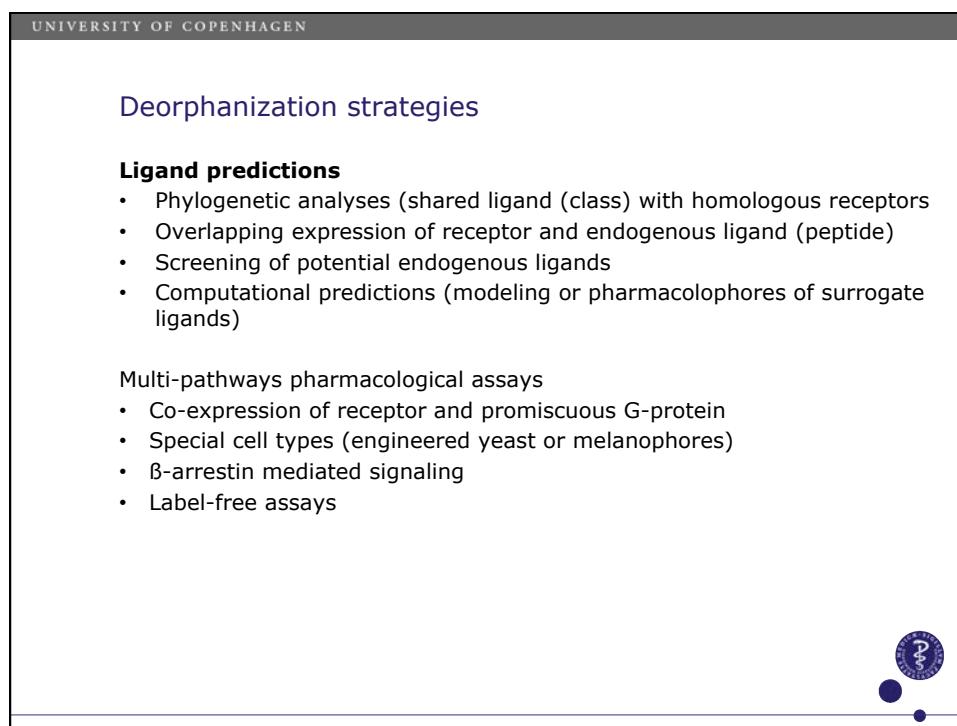
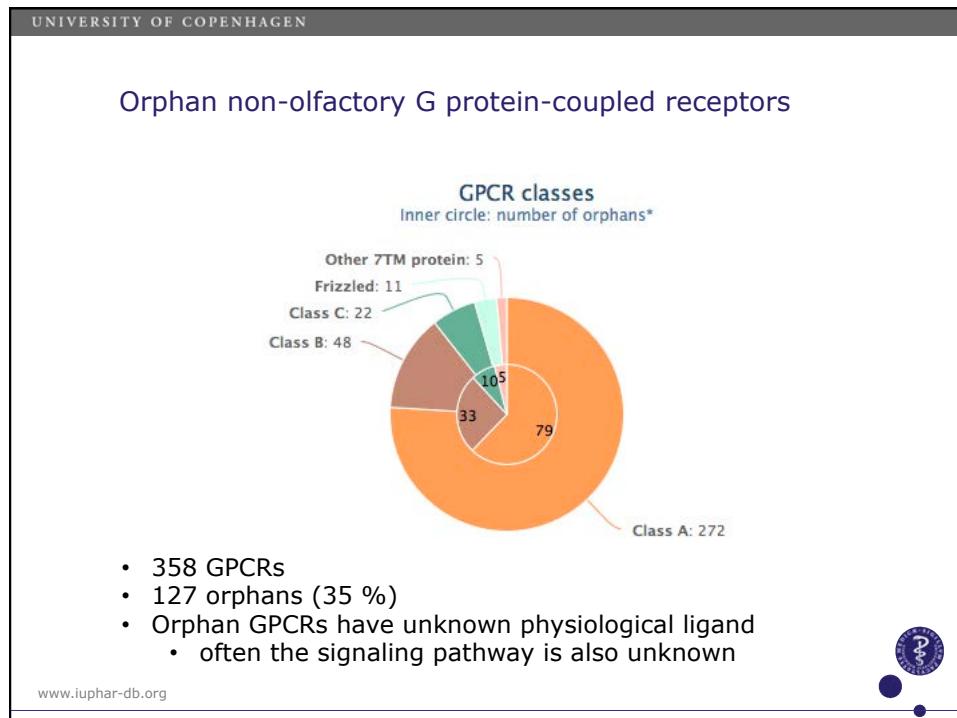


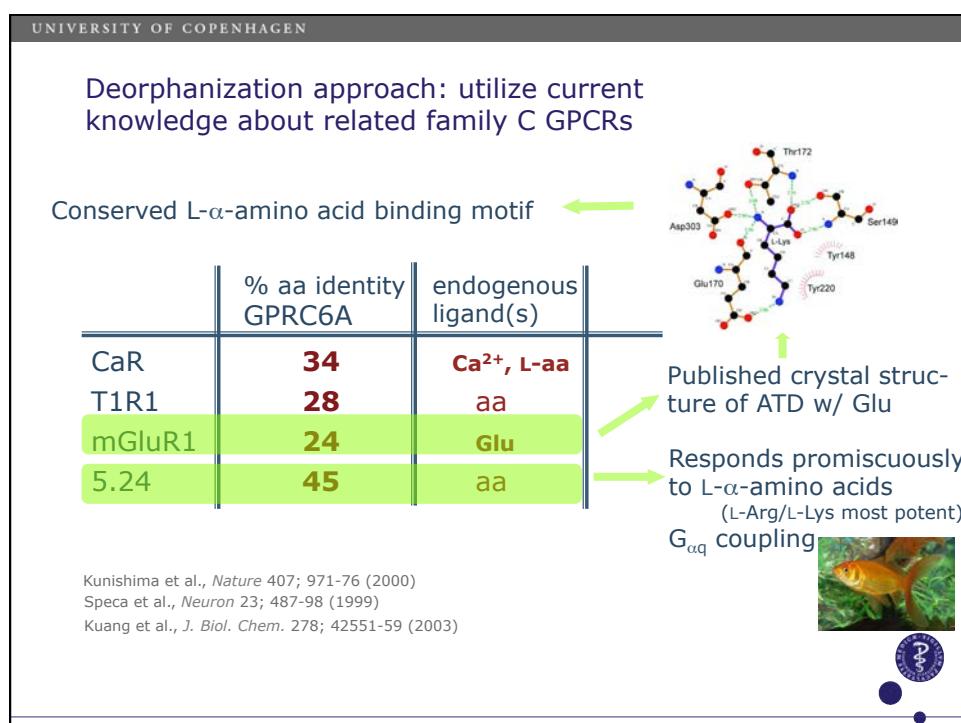
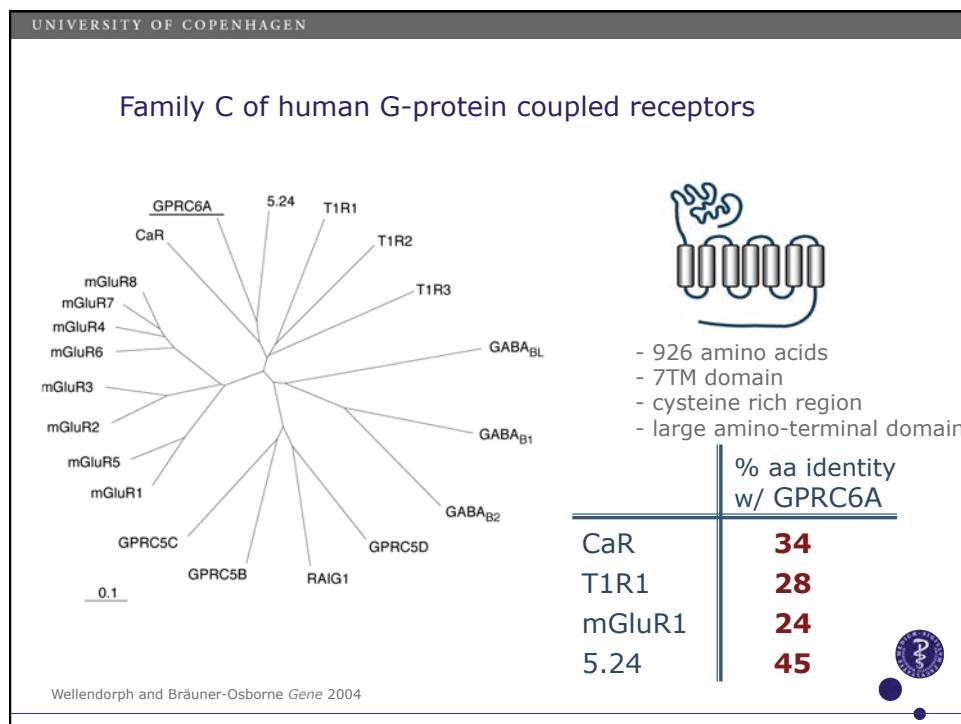












UNIVERSITY OF COPENHAGEN

## Development of pharmacological assay

**Problem:**

- mammalian cell culture media contains high levels of putative agonists

**Solution:**

Expression in *Xenopus* oocytes

