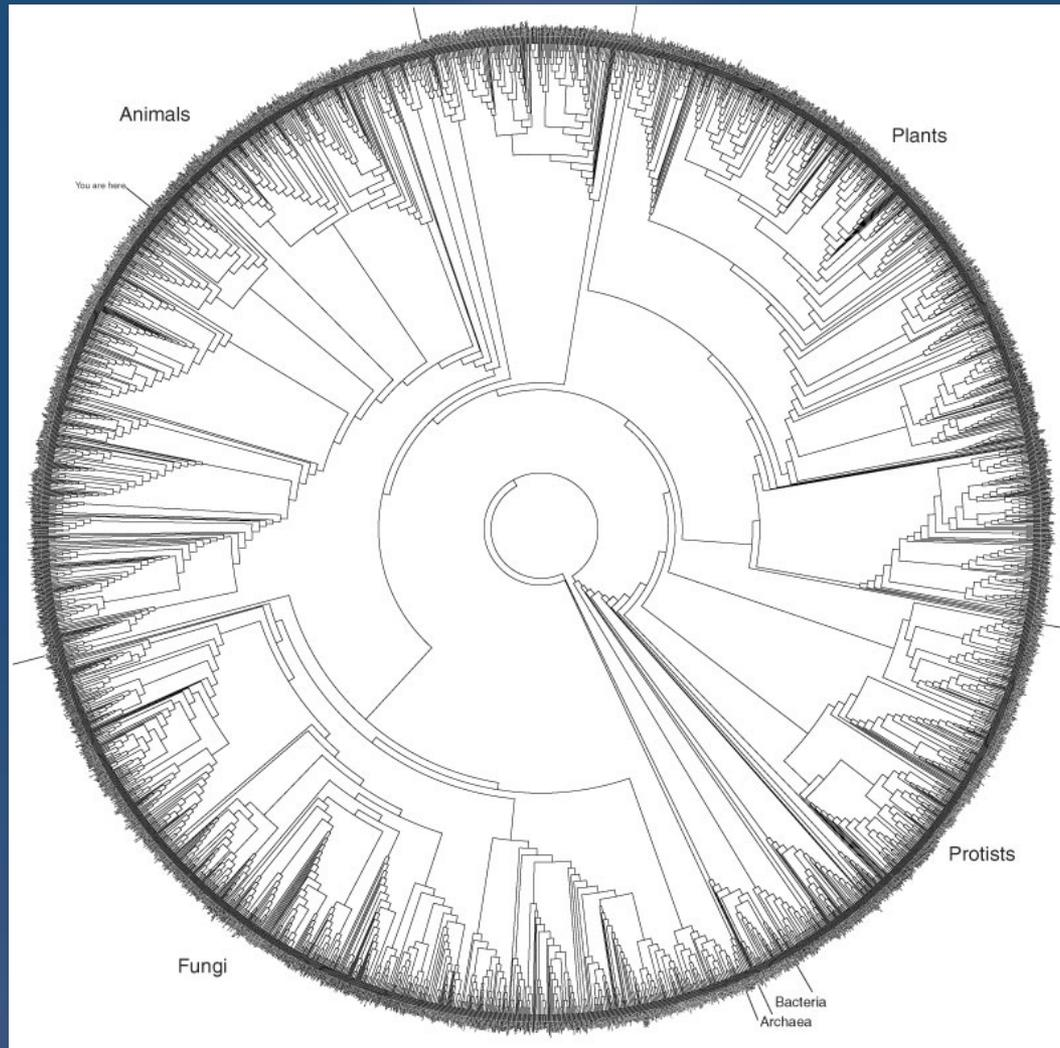




# The use of mtDNA analysis to identify animal species

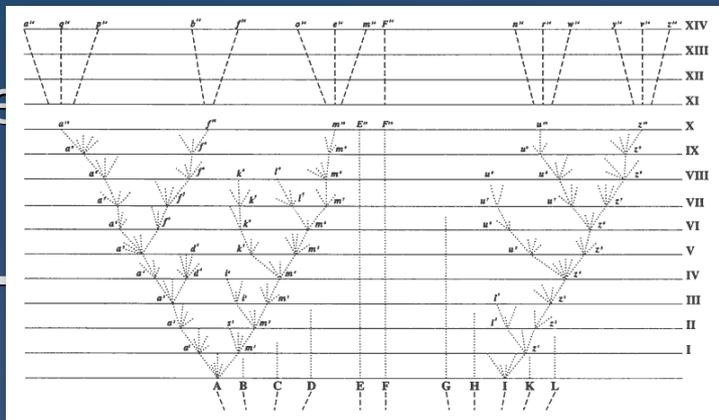


# Introduction to molecular phylogeny



All life forms on earth share a common origin, and their ancestries can be traced back to one or a few organisms that lived approx. 4 billion years ago.

→ all animals, plants and bacteria are related by descent to



organisms are descended from more recent  
C. Darwin – *Origin of Species* (1859)  
Illustration of 'Descent with modification'

common ancestors than are distantly related ones

# Phylogenetic hypothesis



- **Phylogenetic hypothesis**

all living beings share a common ancestor, from which they diverged by accumulating DNA substitutions.

→ two homologous DNA sequences will be more similar

if their common ancestor is recent than if the common ancestor

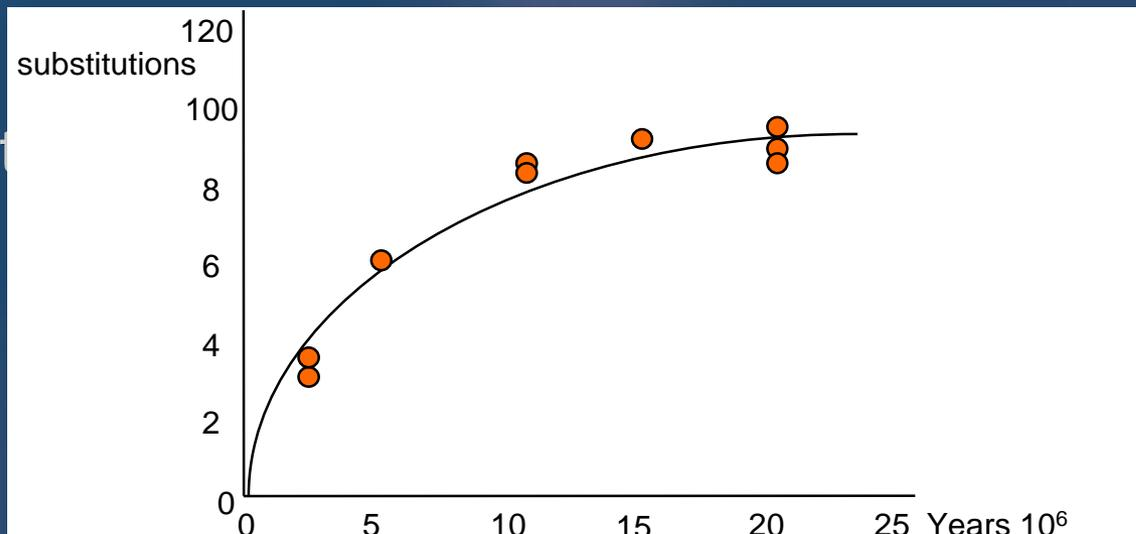
# Phylogenetic hypothesis



## Distance measures for nucleotide sequences

- *Observed differences*: count the nr. of nucleotide sites

at which t



- Sequences separated by 20 Myr are no more different than sequences that diverged 15 Myr
- Relationship not linear, but deflected downwards due to multiple hits (most of the sites changing have already changed before)

# Animal mtDNA



Non-recombining

Fast-evolving

Uniparental inheritance

High copy number

Ubiquitous

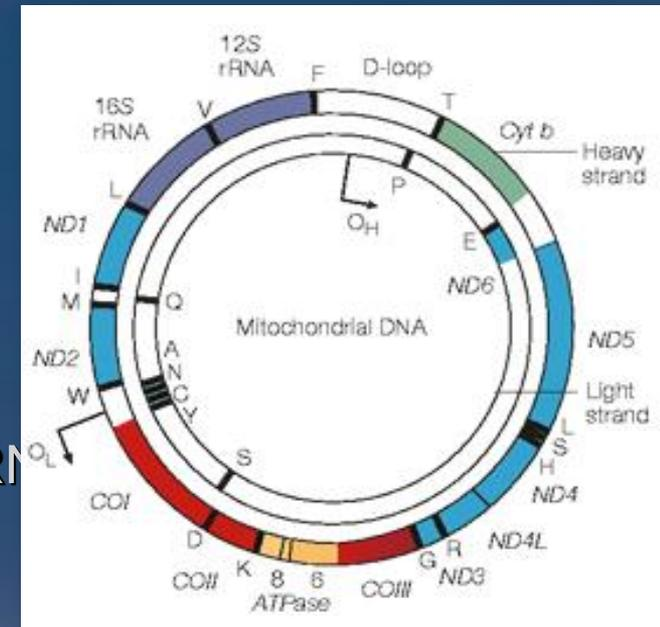
Simple structure

37 genes coding for 13 proteins, 22 tRNAs, 2 rRNAs

Gene order  $\pm$  conserved

Small size [human: 16.5kb]

Large database available



A milestone in evolutionary genetics, mtDNA-RFLP analysis inaugurated and dominated the field of phylogeography in the 1980s.

Still widely employed, direct sequencing of individual genes, regular or chip-based sequencing of whole mtDNA genomes (*mitogenomics*).

# Animal mtDNA



## Why using the mitochondrial genome for species ID?

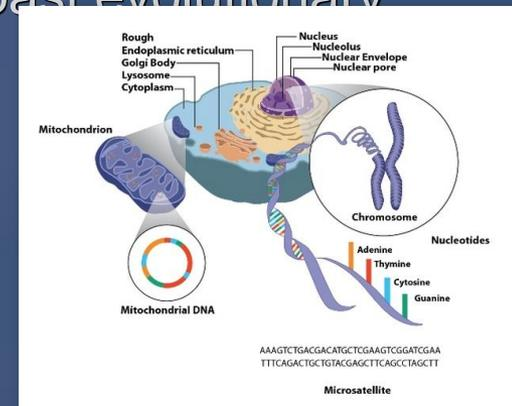
Rapid evolution: mtDNA is *polymorphic* at the intraspecific level, but less than at the interspecific level.

One quarter the effective population size (haploid and uniparental) of nuclear loci means populations *diverge quickly* with respect to nDNA.

*Non-recombining, maternal* inheritance so haplotypes can be ordered phylogenetically into a *gene genealogy* interpretable as the matriarchal component of an *organismal pedigree* (good record of past evolutionary events).

Easy to amplify (many mitochondria per cell).

Large database available.



# Animal mtDNA



*Proc. Natl. Acad. Sci. USA*  
Vol. 86, pp. 6196–6200, August 1989  
Evolution

## Dynamics of mitochondrial DNA evolution in animals: Amplification and sequencing with conserved primers

(cytochrome *b*/12S ribosomal DNA/control region/evolutionary genetics/molecular phylogenies)

T. D. KOCHER\*, W. K. THOMAS\*, A. MEYER\*<sup>†‡</sup>, S. V. EDWARDS\*<sup>†‡</sup>, S. PÄÄBO\*, F. X. VILLABLANCA<sup>†‡</sup>,  
AND A. C. WILSON\*

Departments of \*Biochemistry and <sup>†</sup>Zoology, and <sup>‡</sup>Museum of Vertebrate Zoology, University of California, Berkeley, CA 94720

Universal mtDNA animal primers:

Cytochrome *b*, 12S rRNA and control region

Phylogenetics

Systematics

Population genetics

Forensics

(...)



Species ID

# Animal mtDNA



## Evolution of the Cytochrome *b* Gene of Mammals

David M. Irwin, Thomas D. Kocher,\* and Allan C. Wilson

Division of Biochemistry and Molecular Biology, University of California, Berkeley, CA 94720, USA

J Mol Evol (1991) 32:128–144

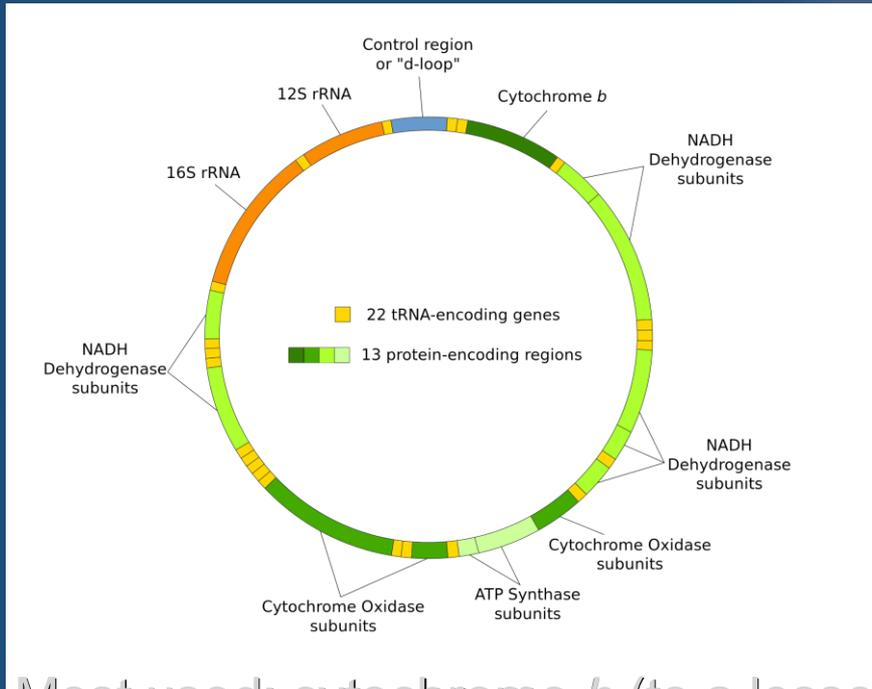
mtDNA genetic code

Substitution rates depend on:

codon position: 1st, 2nd and 3rd positions ratio 10:1:35

structure/function model: residues in transmembrane region evolve faster

# Animal mtDNA

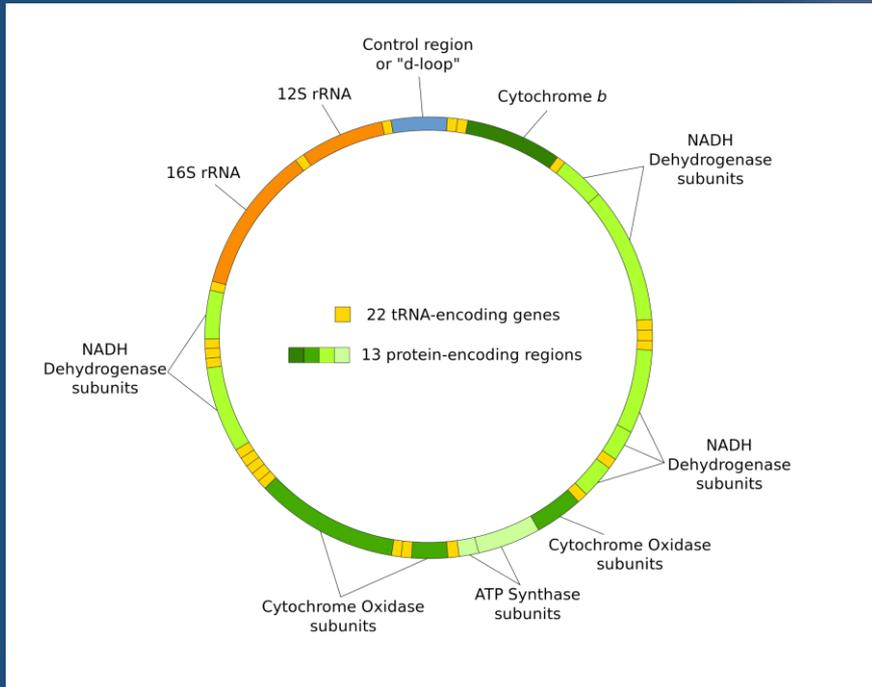


Most used: cytochrome *b* (to a lesser amount 12S and 16S rRNA, COI\*)

higher degree of variability among than within species,  
availability of universal primer pairs amplifying virtually any vertebrate  
spp,  
substantial amount of DNA sequences deposited in DNA databases

\*DNA barcoding

# Animal mtDNA



Most used: cytochrome *b* (to a lesser amount 12S and 16S rRNA, COI\*)

RESTRICTION: limited variability as a coding region, therefore:

inability to differentiate closely related species;  
identical within all vertebrate lineages, impeding the discrimination by direct DNA sequencing of different spp. in mixed samples when using a single universal primer pair

# The use of mtDNA analysis to identify animal species



Alignment of 200 bp of the mtDNA control region of three canid species (*Canis lupus*:  
*C. latrans*: coyote; *Vulpes vulpes*: red fox)

```
Canis lupus      TCCAGGTA AACCTTCTTCCCTC-CCCTATGTACGTCGTGCATTAATGGTTTGCCCATGCATAT-AAGCATGTACATAATATTATATTCTTACATAGGA
Canis latrans   ....AA-...T.....-----..C.....C.....-.....-.....CCT.....
Vulpes vulpes   C.GC.C.----.---.AAAA..TG.....C..C.A.....T.....TAC.....AA.....A..

Canis lupus      CATATCAACTC-AATCTCACAAATTCATTGATCTATCAACAGTAA-TCAAATGCATATCACTTAGTCCAATAAGGGCTTAATCACCATGCCTCGAGAAACC
Canis latrans   ....CTC.ACTT..C..T...G....C.....-.....TGG.T.....C.....
Vulpes vulpes   ....CT.TG.TT.....T.....AATC.CTATC.GG....-..T-ATG.C....CG.....G.....A..T.....
```

# The use of mtDNA analysis to identify animal species



Alignment of 200 bp of the mtDNA *cyt b* gene in 11 specimens of shrews belonging to the genus *Sorex*

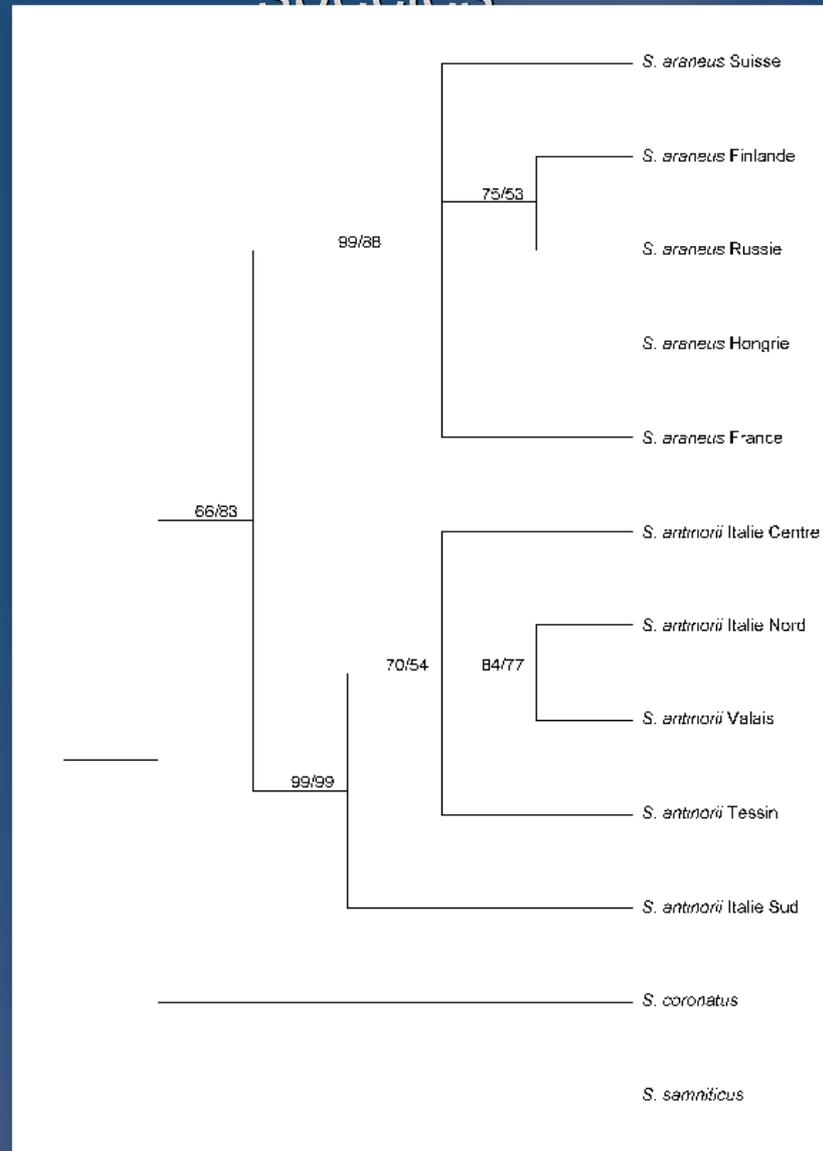
```

ara Suisse      TTTCATCCTAACCCCTTACATCCCTAGTATTATTCTCCCCAGACTTATTAGGAGACCCAGACAACTATATACCTGCAAATCCCCTCAATACACCACCCCA
ara Finlande   .....C.....
ara Hongrie     .....C.....G....
ara Russie     .....T.....C.....
ara France     .....C.....
anti Italie Centre .....G.....C.C.....C.....
anti Valais    .....G.....C.C.....C.....
anti Italie Nord .....G.....C.C.....C.....
anti Tessin    .....G.....C.C.....C.....
anti Italie Sud .....G.....G.....C.C.....C.....
coronatus     .C.....G.....G.....C.C.....
    
```

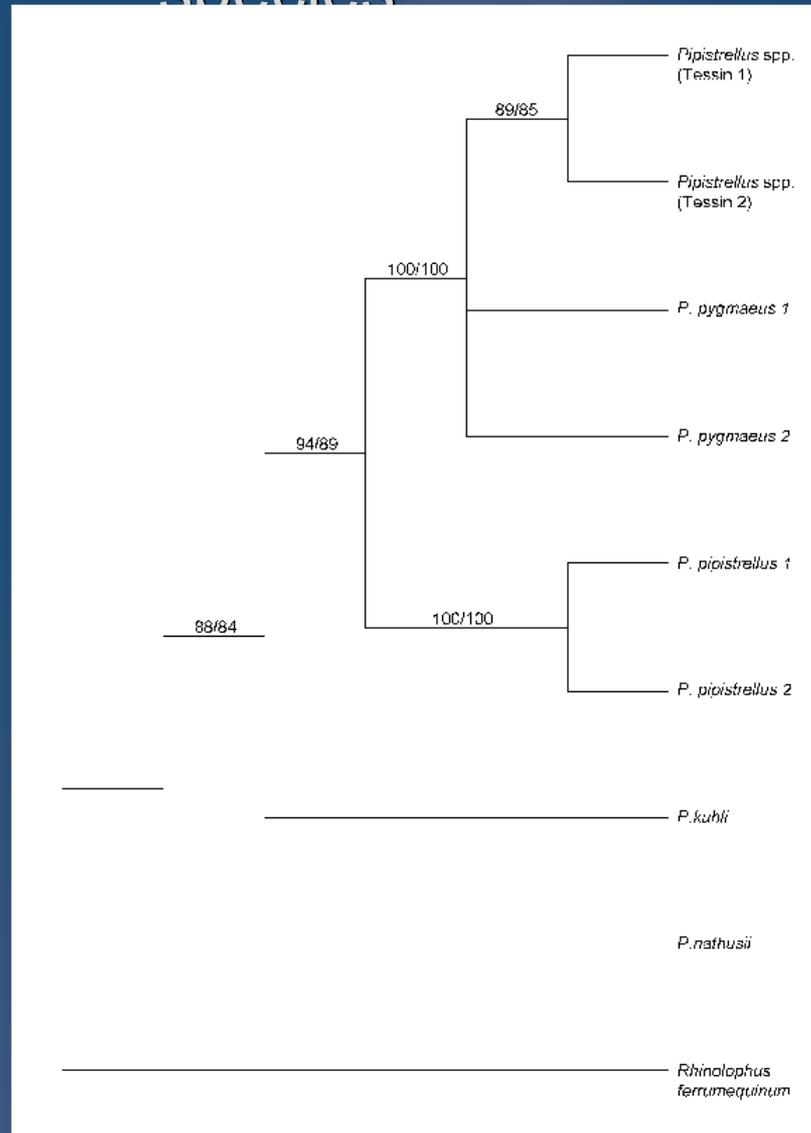
```

ara Suisse      TATTAAACCAGAATGGTATTTTCTATTGCTTACGCCATTCTACGATCTATCCCTAACAACTAGGGGGCGTCCCTAGCACTAGCCCTATCAATTTTAATT
ara Finlande   .....A.....T.....C.....
ara Hongrie     .....A.....T.....
ara Russie     .....A.....T.....
ara France     .....A.....T.....
anti Italie Centre .....A.C.....A.....T.....C.....
anti Valais    .....A.....A.....T.....C.....
anti Italie Nord .....A.....A.....T.....C.....
anti Tessin    .....A.....A.....T.....C.....
anti Italie Sud .....A.C.....A.....T.....C.....
coronatus     .....A.....C.....T.....T.....T.G.....T.....
    
```

# The use of mtDNA analysis to identify animal species



# The use of mtDNA analysis to identify animal species



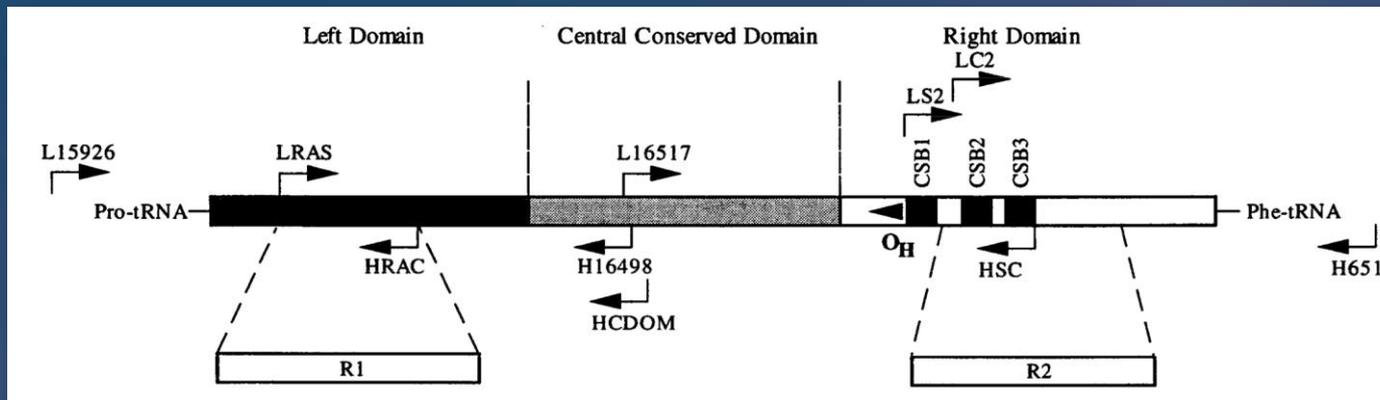
# Animal mtDNA



A single non-coding region: control region

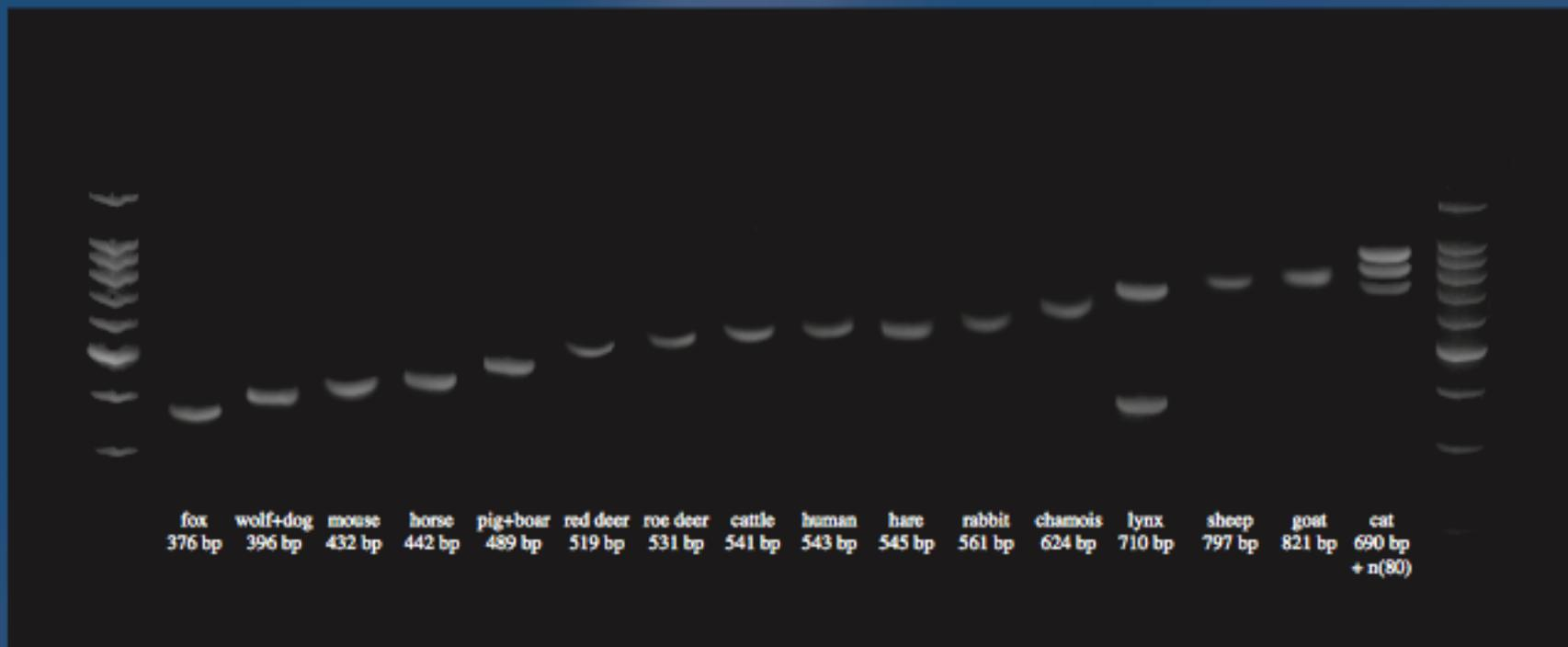
RAPID EVOLUTION, therefore more effective than *cyt b* and other mtDNA coding regions for the identification of closely related species

Presence of tandem repeats: sequence length polymorphism, useful for DNA mixtures



(Fumagalli et al. 1996)

- Species ID



otter 348 bp  
badger 349 bp  
bear 410 bp

(Pun et al. 2009)

# • Species ID



## Animal attacks on animal

- Dead pelicans in a zoo

- Predator from the zoo?



*Dama dama*

*Canis familiaris*

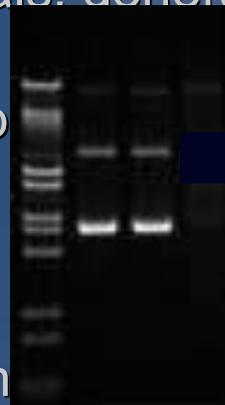
- swabs around wounds
- mtDNA control region: 2 amplicons of 396 bp et de 518 bp, ID by sequencing
- Dog probable predator, fallow deer contamination during storing

# • Species ID



## Animal attacks on animal

- 40+ cases in summer 2005 in NW Switzerland
- Mutilations on domestic animals. generally found dead
- cows, horses, donkeys, sheep
- sexual organs, face, tail, ears
- alleged human mutilator, but no human DNA found
  - Swabs around two dead cows
  - mtDNA control region: 2 amplicons of 541 bp et 376 bp, ID by sequencing
  - Probable *post-mortem* interference



← *Bos taurus*  
← *Vulpes vulpes*

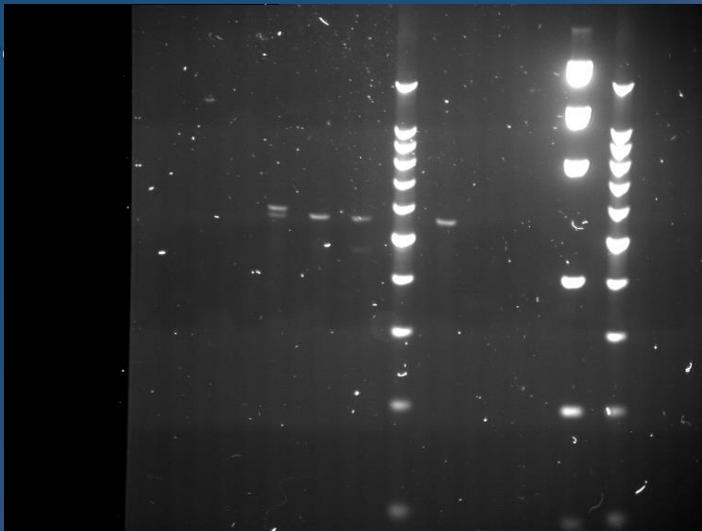
- Species ID



Expertise for exportation permit

- Ancient african mask (ca. 100 years), value ca. 25'000 \$

- Le



- mtDNA control region: amplicons of. 550 bp, ID by sequencing

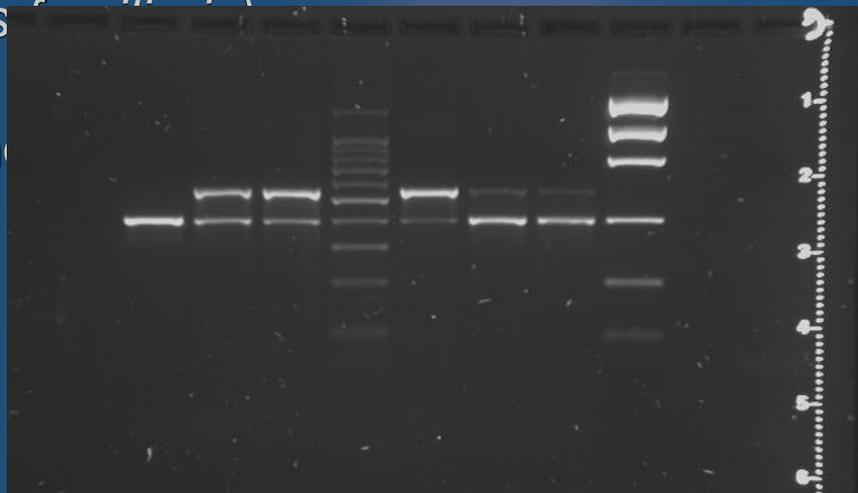
- *Genetta maculata* ou *G. tigrina*

# • Species ID



## Animal attack on human

- dog bite (*Canis familiaris*)
- swabs on clothing



← *Homo sapiens*  
← *Canis familiaris*

- mtDNA control region: 2 amplicons of 543 bp et 396 bp, ID by sequencing
- DNA profiling: 8 STR loci
- Two reference dogs

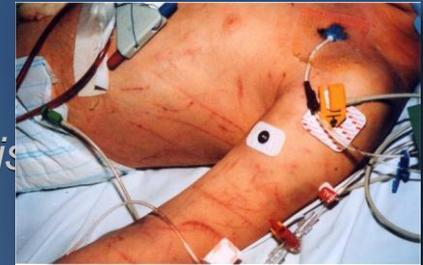
- 6 samples same DNA profile, identical to one of the references,
- allelic  $f$  in a reference population of 63 dogs of various breeds, correction factor theta of 3%
- $LR > 1 \times 10^9$

# • Species ID



## Alleged human attack on human

- Child laying unconscious in the snow, most of his clothes removed or torn, and several wounds on his body. Presence of the family dog (German sheperd).



- mtDNA control region: 30 bands of expected size for *Canis familiaris*
- No human DNA on child except its own
  - Confirmation *C. familiaris* by DNA sequencing on 13 bands
- 5 DNA profiles by 7 samples (our sampling, 8 S (R) dog and clothes), and by us 2
  - Saliva from the family dog (reference)
    - DNA profiles 7 samples identical, and identical to reference
    - allelic  $f$  in a reference population of 95 dogs of various breeds,
      - $P_{(ID)random} < 8/10^{13}$
- Wounds compatible with dog attack (controversial)
  - $P_{(ID)sib} < 1/910^4$

years later (clothes)

- Species ID

Accident caused by animals



- bird strike
- swabs over spots on engine



- mtDNA cyt *b* bird specific universal primers  
→ 3 swabs, species identified: Common kestrel (*Falco tinnunculus*)



- Species ID

Illegal animal traffic (poaching)

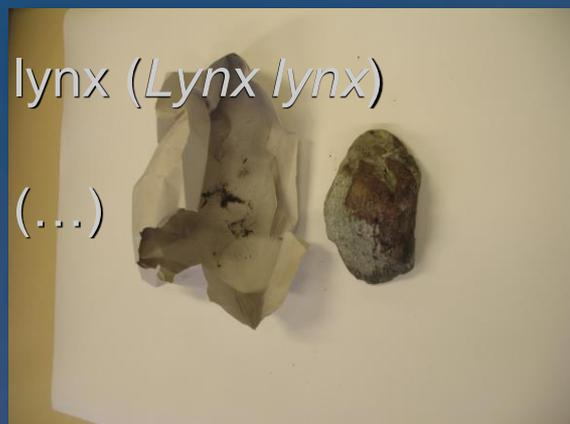


roe deer (*Capreolus capreolus*)

red deer (*Cervus elaphus*)

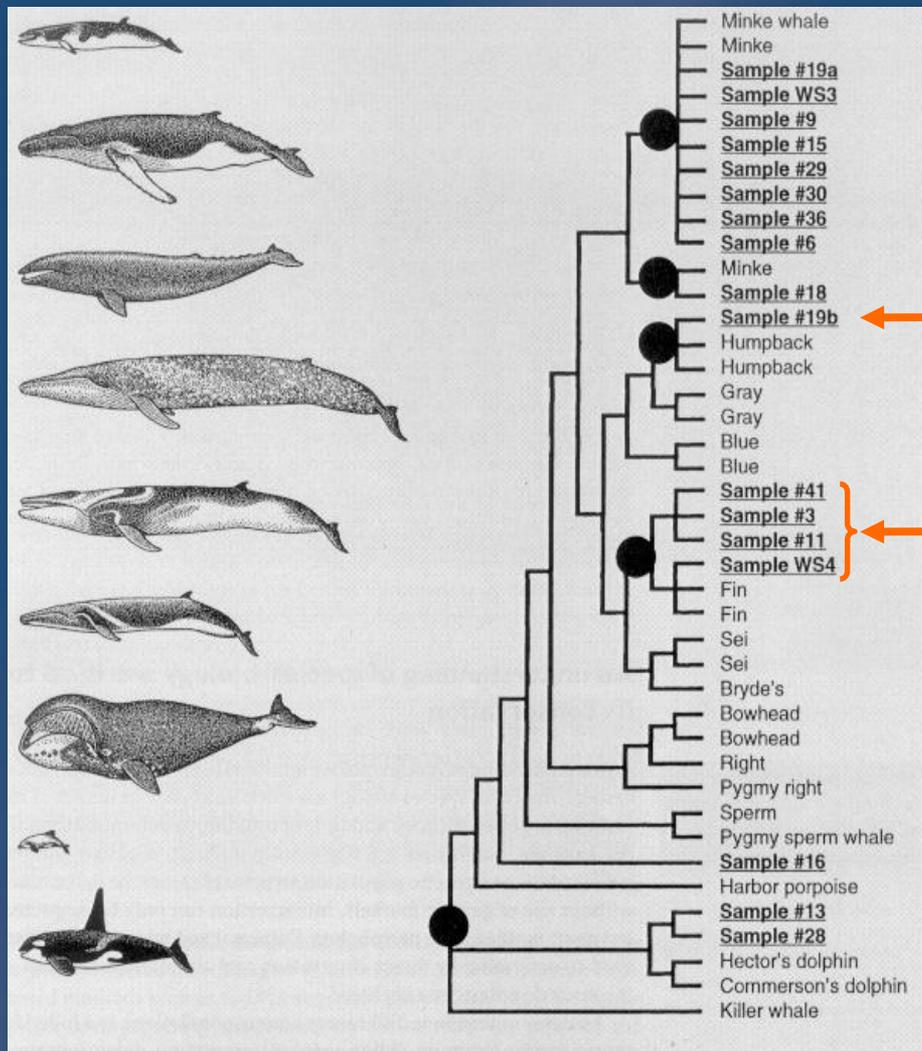
wild boar (*Sus scrofa*)

wolf (*Canis lupus*)



mtDNA: Non-invasive genetic sampling and aDNA

# Species ID



(after Baker & Palumbi 1996)

# • Species ID



## DNA microarrays



### DNA mixtures and multiple species

*Journal of Applied Ecology* 2008, **45**, 967–975

doi: 10.1111/j.1365-2664.2007.01415.x

## Molecular identification of vertebrate species by oligonucleotide microarray in food and forensic samples

Fabrice Teletchea<sup>1,†</sup>, Jacques Bernillon<sup>2</sup>, Marilynne Duffraisie<sup>1,3</sup>, Vincent Laudet<sup>3</sup> and Catherine Hänni<sup>1,3\*</sup>

mtDNA *cyt b* gene  
373 bp (268+124 bp)  
71/77 vertebrate spp.  
unambig. identified  
(mammals, birds, fish)

OPEN ACCESS Freely available online



## Identifying Fishes through DNA Barcodes and Microarrays

Marc Kochzius<sup>1,2\*</sup>, Christian Seidel<sup>1a</sup>, Aglaia Antoniou<sup>3</sup>, Sandeep Kumar Botla<sup>1b</sup>, Daniel Campo<sup>4c</sup>, Alessia Cariani<sup>5</sup>, Eva Garcia Vazquez<sup>4</sup>, Janet Hauschild<sup>1d</sup>, Caroline Hervet<sup>6e</sup>, Sigridur Hjörleifsdottir<sup>7</sup>, Gudmundur Hreggvidsson<sup>7,8</sup>, Kristina Kappel<sup>1</sup>, Monica Landi<sup>5f</sup>, Antonios Magoulas<sup>3</sup>, Viggo Marteinson<sup>7</sup>, Manfred Nölte<sup>9</sup>, Serge Planes<sup>6g</sup>, Fausto Tinti<sup>5</sup>, Cemal Turan<sup>10</sup>, Moleyur N. Venugopal<sup>11</sup>, Hannes Weber<sup>1</sup>, Dietmar Blohm<sup>1</sup>

mtDNA *cyt b*, 16S,  
COI  
30 out of 50 fish spp.  
identified

RESTRICTIONS: development time, costs, no flexibility (species

determined)

- Species ID



Next generation sequencing



A Universal Method for Species Identification of Mammals Utilizing Next Generation Sequencing for the Analysis of DNA Mixtures

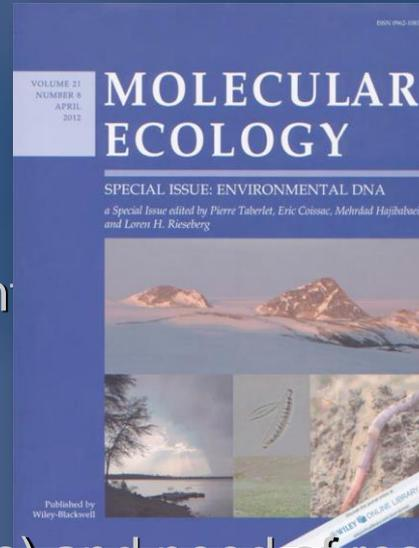
Andreas O. Tillmar, Barbara Dell'Amico, Jenny Welander, Gunilla Holmlund

2013. PLoS ONE 8(12).

eDNA

Biodiversity assessment

Diet analysis



RESTRICTIONS: Slow (few samples) and need of results on a week-by-

x basis

(...)

- Species ID



Next generation sequencing

## Population assignment (SNPs)

### ARTICLE

Received 23 Dec 2011 | Accepted 12 Apr 2012 | Published 22 May 2012

DOI:10.1038/ncomms1845

## Gene-associated markers provide tools for tackling illegal fishing and false eco-certification

Einar E. Nielsen<sup>1</sup>, Alessia Cariani<sup>2,3</sup>, Eoin Mac Aoidh<sup>4</sup>, Gregory E. Maes<sup>3</sup>, Ilaria Milano<sup>2,5</sup>, Rob Ogden<sup>6</sup>, Martin Taylor<sup>7</sup>, Jakob Hemmer-Hansen<sup>1</sup>, Massimiliano Babbucci<sup>5</sup>, Luca Bargelloni<sup>5</sup>, Dorte Bekkevold<sup>1</sup>, Eveline Diopere<sup>3</sup>, Leonie Grenfell<sup>6</sup>, Sarah Helyar<sup>8</sup>, Morten T. Limborg<sup>1</sup>, Jann T. Martinsohn<sup>4</sup>, Ross McEwing<sup>6</sup>, Frank Panitz<sup>9</sup>, Tomaso Patarnello<sup>5</sup>, Fausto Tinti<sup>2</sup>, Jeroen K.J. Van Houdt<sup>3</sup>, Filip A.M. Volckaert<sup>3</sup>, Robin S. Waples<sup>10</sup>, FishPopTrace Consortium\* & Gary R. Carvalho<sup>7</sup>

NATURE COMMUNICATIONS | 3:851 | DOI: 10.1038/ncomms1845 | www.nature.com/naturecommunications

- Difficulty and constraints



- Very high number of different species

### Development of methods

- no commercial kits

- research and development only exclusively performed in « non-

forensic » research laboratories

- From the university to the forensic laboratory

