

A decorative graphic on the left side of the slide consists of a white hexagonal grid pattern. Several hexagons contain scientific images: a DNA double helix, purple rod-shaped bacteria, a colorful fluorescence micrograph of cells, and a person in a lab coat working with petri dishes.

Reclassification of the
Mycobacterium tuberculosis
Complex (MTBC) Species as
Mycobacterium tuberculosis

Marco A. Riojas, Ph.D.
Scientist, ATCC/BEI Resources



Outline

INTERNATIONAL
JOURNAL OF SYSTEMATIC
AND EVOLUTIONARY
MICROBIOLOGY

RESEARCH ARTICLE
Riojas et al., *Int J Syst Evol Microbiol* 2018;68:324–332
DOI 10.1099/ijs.0.002507



Phylogenomic analysis of the species of the *Mycobacterium tuberculosis* complex demonstrates that *Mycobacterium africanum*, *Mycobacterium bovis*, *Mycobacterium caprae*, *Mycobacterium microti* and *Mycobacterium pinnipedii* are later heterotypic synonyms of *Mycobacterium tuberculosis*

Marco A. Riojas,^{1*} Katya J. McGough,^{1,2} Cristin J. Rider-Riojas,³ Nalin Rastogi⁴ and Manzour Hernando Hazbón¹

Abstract

The species within the *Mycobacterium tuberculosis* Complex (MTBC) have undergone numerous taxonomic and nomenclatural changes, leaving the true structure of the MTBC in doubt. We used next-generation sequencing (NGS), digital DNA-DNA hybridization (dDDH), and average nucleotide identity (ANI) to investigate the relationship between these species. The type strains of *Mycobacterium africanum*, *Mycobacterium bovis*, *Mycobacterium caprae*, *Mycobacterium microti* and *Mycobacterium pinnipedii* were sequenced via NGS. Pairwise dDDH and ANI comparisons between these, previously sequenced MTBC type strain genomes (including *Mycobacterium canettii*, *Mycobacterium mungii* and *Mycobacterium orygis*) and *M. tuberculosis* H37Rv were performed. Further, all available genome sequences in GenBank for species in or putatively in the MTBC were compared to H37Rv. Pairwise results indicated that all of the type strains of the species are extremely closely related to each other (dDDH: 91.2–99.2%, ANI: 99.21–99.92%), greatly exceeding the respective species delineation thresholds, thus indicating that they belong to the same species. Results from the GenBank genomes indicate that all the strains examined are within the circumscription of H37Rv (dDDH: 83.5–100%). We, therefore, formally propose a union of the species of the MTBC as *M. tuberculosis*: *M. africanum*, *M. bovis*, *M. caprae*, *M. microti* and *M. pinnipedii* are reclassified as later heterotypic synonyms of *M. tuberculosis*. *M. canettii*, *M. mungii*, and *M. orygis* are classified as strains of the species *M. tuberculosis*. We further recommend use of the infrasubspecific term 'variant' (var.) and infrasubspecific designations that generally retain the historical nomenclature associated with the groups or otherwise convey such characteristics, e.g. *M. tuberculosis* var. *bovis*.

INTRODUCTION

The species within the *Mycobacterium tuberculosis* Complex (MTBC) have undergone numerous taxonomic and nomenclatural changes, leaving the true structure of the MTBC in doubt.

At the time of writing, the species within the MTBC with validly published names are *Mycobacterium tuberculosis* (also the type species of the genus), *Mycobacterium africanum*, *Mycobacterium bovis*, *Mycobacterium caprae*,

Mycobacterium microti and *Mycobacterium pinnipedii*, which are all very closely related [1, 2]. Even these species have undergone some taxonomic and nomenclatural changes. For example, the species most recently known as *M. caprae* was first proposed as *M. tuberculosis* subsp. *caprae* [3], then was later renamed as *M. bovis* subsp. *caprae* [4], and finally elevated to the rank of species [5]. These nomenclatural changes have resulted in equally valid basonyms for the same organism. Numerous other 'species' with similar properties have been identified but have not been

Author affiliations: ¹ATCC[®], Manassas, VA 20110, USA; ²George Mason University, Fairfax, VA 22030, USA; ³Unaffiliated; ⁴Institut Pasteur de la Guadeloupe, Les Abymes 97139, Guadeloupe, France.
*Correspondence: Marco A. Riojas, mriojas@atcc.org
Keywords: *Mycobacterium tuberculosis* complex; *Mycobacterium* reclassification.

Abbreviations: ANI, average nucleotide identity; dDDH, digital DNA-DNA hybridization; MTBC, *Mycobacterium tuberculosis* complex; NGS, next-generation sequencing. The GenBank accession numbers for the whole genome of MWXF01.1, *Mycobacterium tuberculosis* var. *bovis* ATCC[®] 11762, MWXD01.1, *Mycobacterium tuberculosis* var. *microti* ATCC[®] 35061, and MWXB01.1, *Mycobacterium tuberculosis* var. *pinnipedii* ATCC[®] 35062 are available with the online version of this article.

Eight supplementary tables are available with the online version of this article.

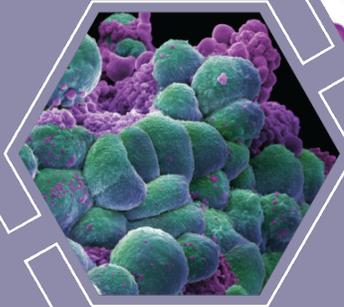
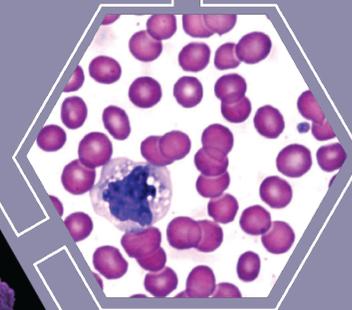
00507 © 2018 by American Type Culture Collection (ATCC). Downloaded from https://academic.oup.com/ijse/advance-article-abstract/doi/10.1093/ijse/ijse00507/5070000 by American Type Culture Collection (ATCC) user on 12 October 2018

Riojas et al., *Int J Syst Evol Microbiol* 68(1): 324–332, 2018

- About ATCC
- Determination of species
 - What is a species?
 - How are species assigned?
 - DNA-DNA hybridization (DDH)
 - Digital DDH (dDDH)
- Research results and conclusions
 - MTBC results
 - Nomenclatural proposal



About ATCC



About ATCC

- Founded in 1925, ATCC is a non-profit organization with HQ in Manassas, VA, and an R&D and Services center in Gaithersburg, MD
- World's premier biological materials resource and standards development organization
 - 5,000 cell lines
 - 80,000 microorganisms
 - Genomic & synthetic nucleic acids
 - Media/reagents
- ATCC collaborates with and supports the scientific community with industry-standard biological products and innovative solutions
 - Growing portfolio of products and services
 - Sales and distribution in 150 countries, 15 international distributors
- Talented team of 450+ employees, over one-third with advanced degrees

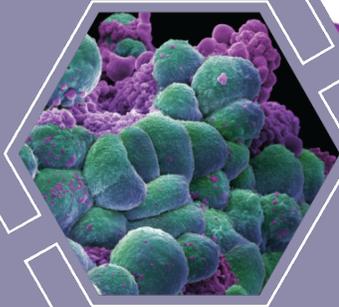
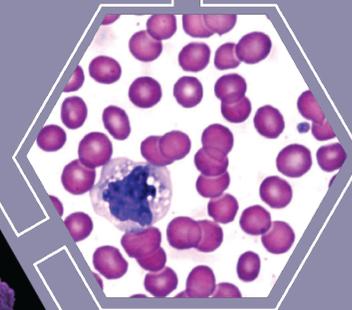


An innovative
global partner for
authentic
biomaterials,
standards and
services





Determination of Species



How are Species Defined?

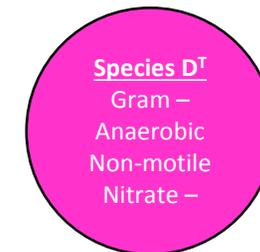
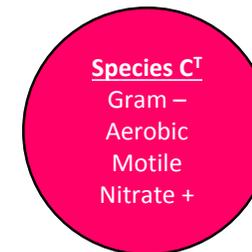
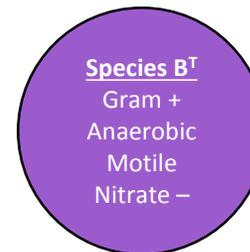
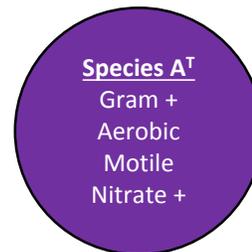


- Each species is represented by a **type strain** and a description of that strain
 - Usually the first strain identified
 - Not necessarily the most typical or the representative of the species
 - Designated with a superscripted, uppercase T after the strain name (e.g., *Mycobacterium tuberculosis* H37Rv^T)
 - **The type strain is essentially the “definition” of a species**
- Serves as the comparative reference point for the identification of new strains and the assignment to existing species

How is a New Strain Assigned to a Species?



- The characteristics of the new strain are compared to the characteristics of species type strains
 - Historically: phenotypic characteristics
- A strain that shares enough of the essential characteristics of a type strain is said to be within the **circumscription** of that species/type strain
 - Therefore, it belongs to that species

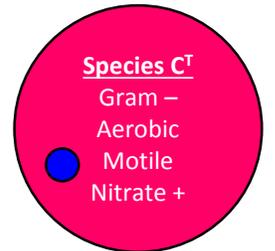


● Gram –
Aerobic
Motile
Nitrate +

How is a New Strain Assigned to a Species?

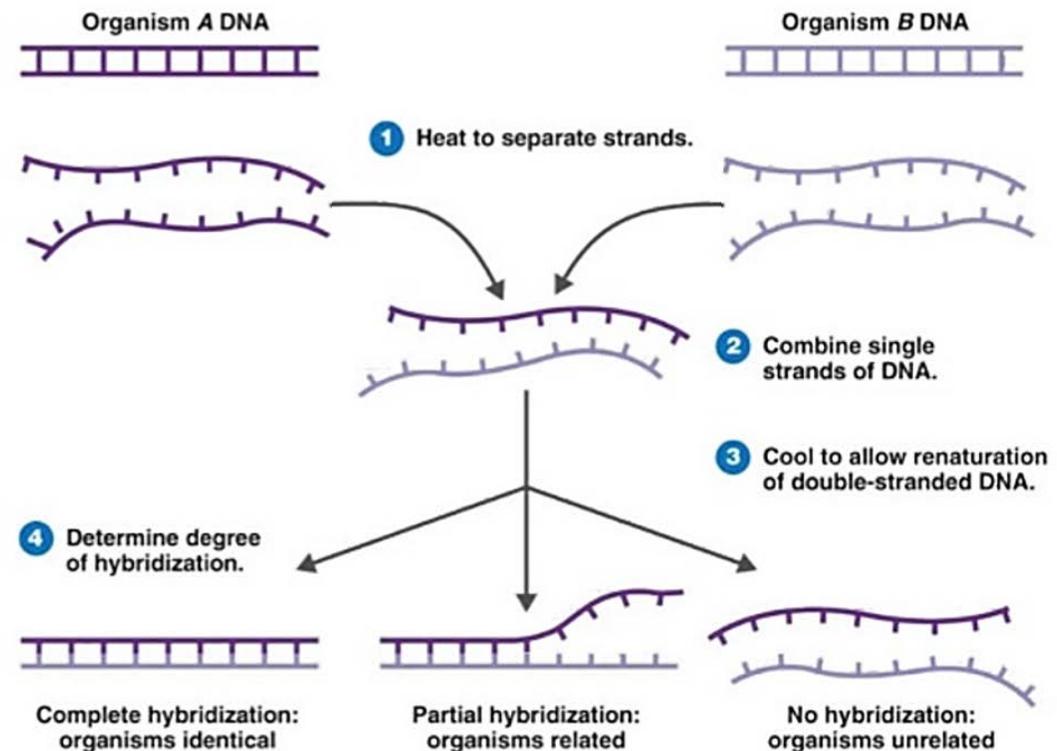


- Recognizing that phenotypes can be quite unreliable, today we rely more heavily on genotypic comparisons
 - 16S rRNA genes, *hsp65*, *rpoB*
 - Multilocus Sequence Typing (MLST)
- Single or small numbers of genes can still provide misleading results
 - Vulnerable to selection bias
- The most objective and accurate comparison would be between whole genomes

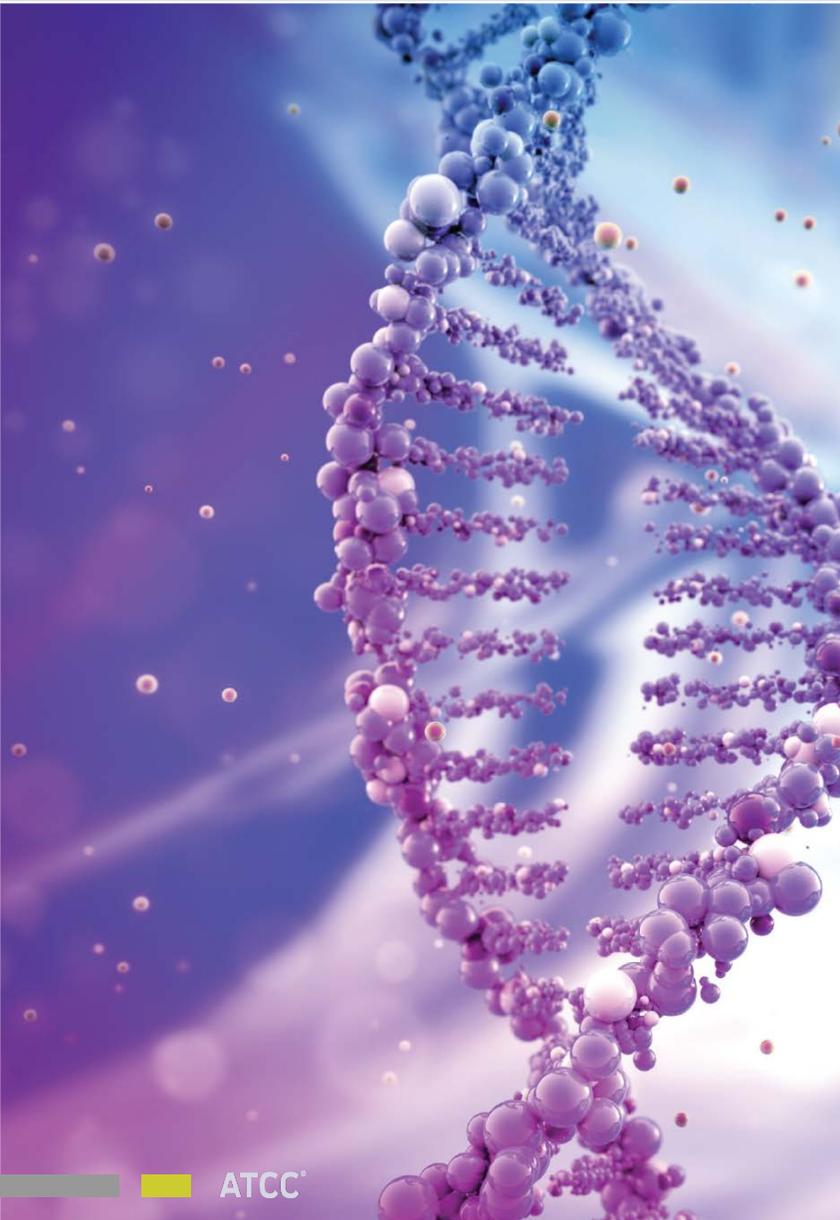


DNA-DNA Hybridization (DDH)

- Denature genomic DNA (gDNA) mixture for organisms A and B
- Allow gDNA to anneal; hybrids result
- Reassociation of gDNA \approx sequence similarity



DNA-DNA Hybridization (DDH)



- Denature genomic DNA (gDNA) mixture for organisms A and B
- Allow gDNA to anneal; hybrids result
- Reassociation of gDNA \approx sequence similarity

- **DDH similarity >70%: same species**
- **DDH similarity <70%: different species**

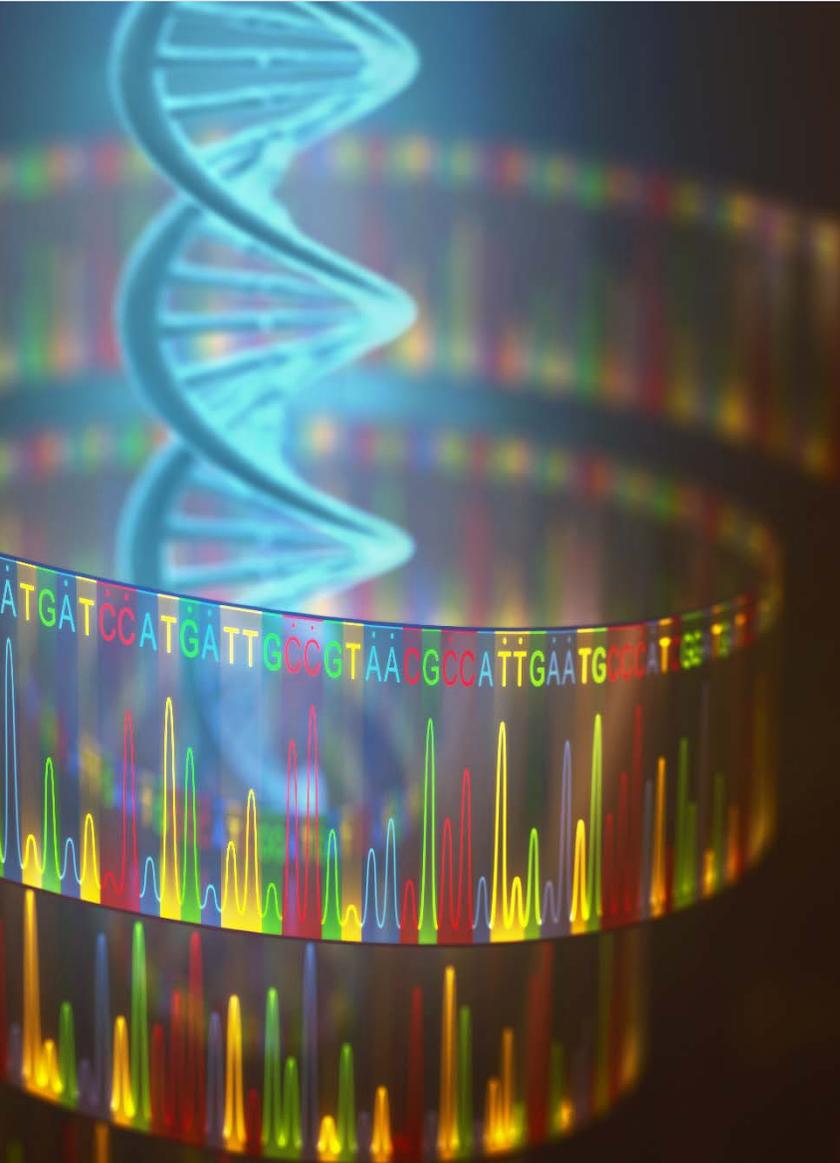
- Considered the gold standard for species definition
 - A putatively new species must be shown to have <70% DDH similarity to existing species (*i.e.*, type strains) to be recognized as a legitimately new species

DNA-DNA Hybridization (DDH)



- DDH poses significant challenges
 - Highly labor intensive
 - Time consuming
 - Error prone
 - Poorly reproducible
- To overcome these challenges, an analogous sequence-based method has been developed
 - Digital DDH (dDDH)
- This method uses the whole-genome sequence of two strains to compare the strains bioinformatically and infer their distance from each other

Digital DNA-DNA Hybridization (dDDH)



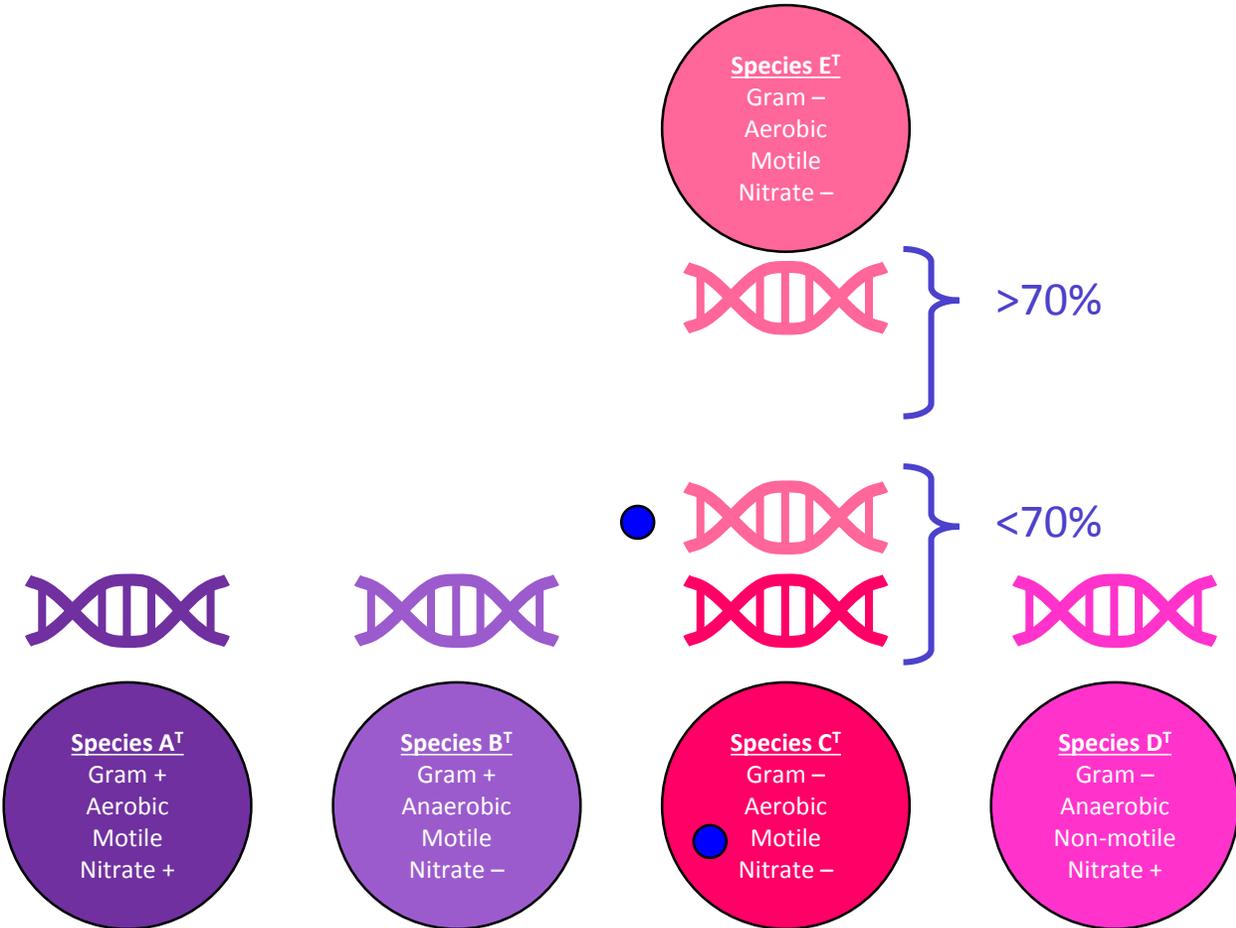
- Available via web server
 - GGDC: Genome-to-Genome Distance Calculator
ggdc.dsmz.de
- Calibrated to the DDH 70% threshold
 - No difference in scale, *e.g.*, 95% average nucleotide identity (ANI)
- Data supports the use of dDDH for assignment of subspecies
 - Delineation threshold proposed at 79–80%

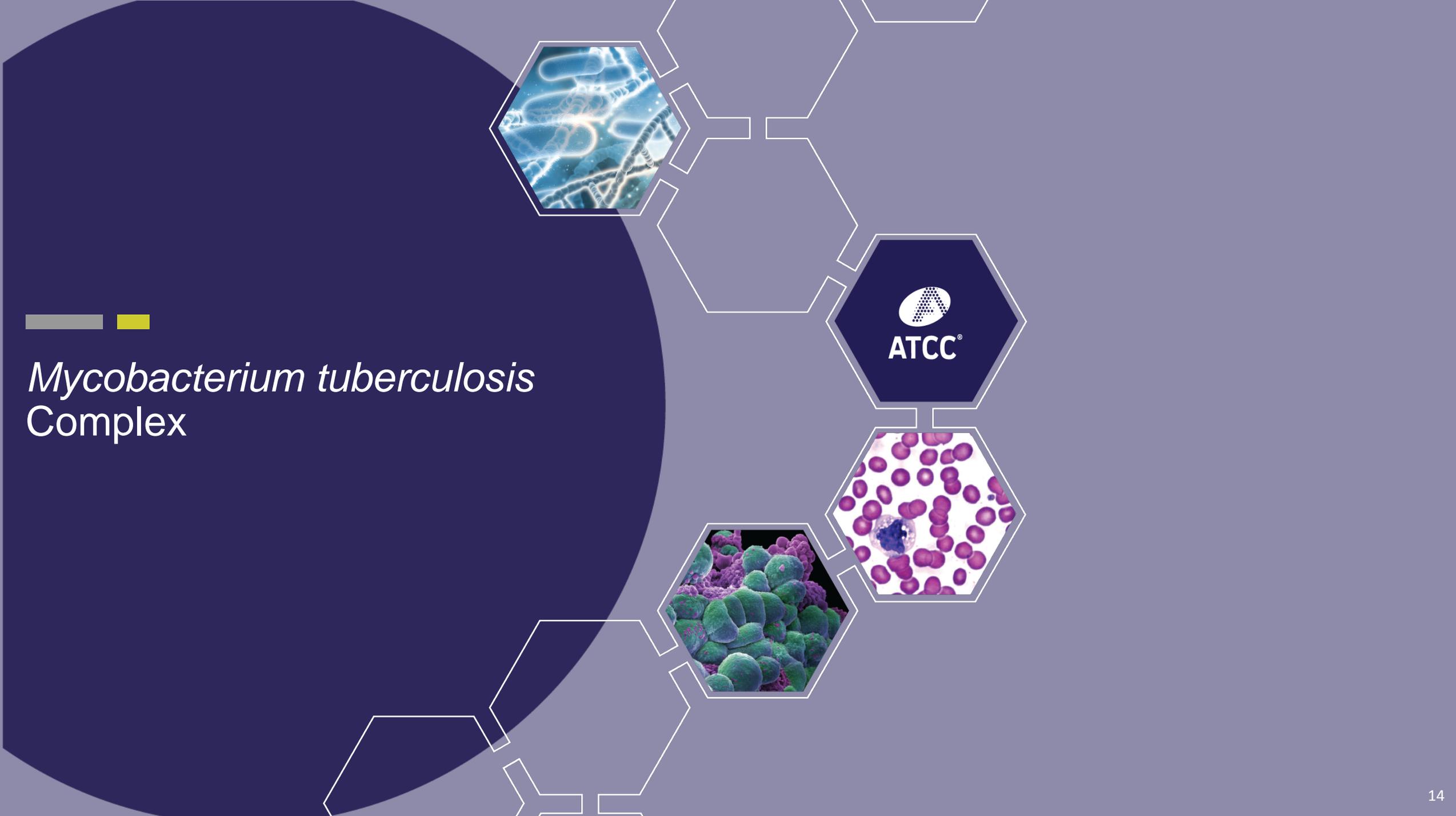
dDDH Range	Interpretation	
≥ 80%	Same species	Same subspecies
70 – 80%	Same species	Different subspecies
< 70%	Different species	

Transfer of Species Based on Genomic Data

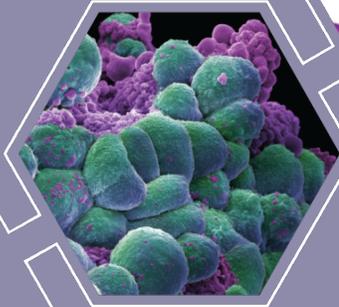
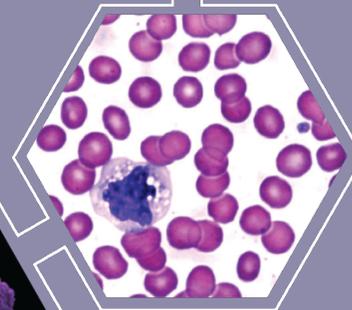


- Genomic analysis can provide updated taxonomy





Mycobacterium tuberculosis
Complex



Mycobacterium tuberculosis Complex (MTBC)



- Consists of six officially recognized (validly published) species, all very closely related
 - *Mycobacterium tuberculosis*
 - *Mycobacterium africanum*
 - *Mycobacterium bovis*
 - *Mycobacterium caprae*
 - *Mycobacterium microti*
 - *Mycobacterium pinnipedii*
 - Not validly published
 - “*Mycobacterium canettii*”
 - “*Mycobacterium mungi*”
 - “*Mycobacterium orygis*”
- It has been hypothesized for several decades that these may actually represent a single species

Mycobacterium tuberculosis Complex (MTBC)



- Recent work has shown that some strains of the related species fall within the circumscription of *M. tuberculosis* H37Rv^T

Species	Strain
<i>Mycobacterium tuberculosis</i>	H37Rv ^T
<i>Mycobacterium tuberculosis</i>	H37Ra
<i>Mycobacterium tuberculosis</i>	CDC15151
<i>Mycobacterium tuberculosis</i>	F11
<i>Mycobacterium tuberculosis</i>	KZN 1435
<i>Mycobacterium bovis</i>	AF 2122/97
<i>Mycobacterium bovis</i>	Pasteur 1173P2
<i>Mycobacterium bovis</i>	Tokio 172
<i>Mycobacterium africanum</i>	GM041182
" <i>Mycobacterium canettii</i> "	CIPT 140010059

Mycobacterium tuberculosis Complex (MTBC)



- Recent work has shown that some strains of the related species fall within the circumscription of *M. tuberculosis* H37Rv^T
 - However, all strains compared were **non-type strains**
 - Conclusion: “mycobacterial scientists should agree an accord that designates MTBC as a single species in the official taxonomic rules of nomenclature.”
 - Only applicable to those strains, not the species as a whole

Species	Strain
<i>Mycobacterium tuberculosis</i>	H37Rv ^T
<i>Mycobacterium tuberculosis</i>	H37Ra
<i>Mycobacterium tuberculosis</i>	CDC15151
<i>Mycobacterium tuberculosis</i>	F11
<i>Mycobacterium tuberculosis</i>	KZN 1435
<i>Mycobacterium bovis</i>	AF 2122/97
<i>Mycobacterium bovis</i>	Pasteur 1173P2
<i>Mycobacterium bovis</i>	Tokio 172
<i>Mycobacterium africanum</i>	GM041182
“ <i>Mycobacterium canettii</i> ”	CIPT 140010059



Sequencing of MTBC Type Strains



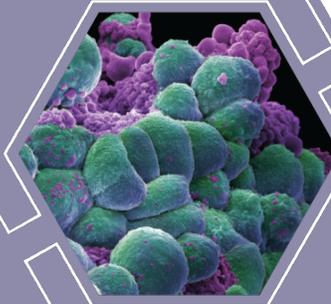
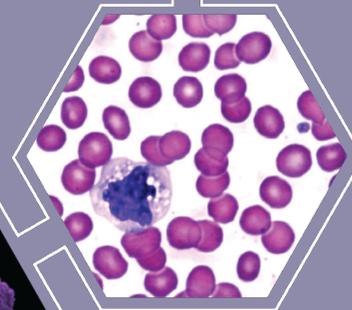
- Extracted DNA from the type strains of the missing five species
- Phylogenomically compared the WGS sequences of these 5 species, *M. tuberculosis* H37Rv^T, and 3 non-validly published “species”, plus 2 outgroups
 - Digital DNA-DNA hybridization (dDDH)
 - Average nucleotide identity (ANI)

Table 1. Genomes of type (or treated as type) strains of the MTBC

Current organism name	Strain	Genome	Genome source
<i>M. tuberculosis</i>	H37Rv ^T	NC_000962.3	GenBank
<i>M. africanum</i>	ATCC 25420 ^T	MWXF01.1	This Work
<i>M. bovis</i>	ATCC 19210 ^T	MWXE01.1	This Work
<i>M. caprae</i>	ATCC BAA-824 ^T	MWXD01.1	This Work
<i>M. microti</i>	ATCC 19422 ^T	MWXC01.1	This Work
<i>M. pinnipedii</i>	ATCC BAA-688 ^T	MWXB01.1	This Work
' <i>M. canettii</i> '	CIPT 140010059	NC_015848.1	GenBank
' <i>M. mungi</i> '	BM22813	LXTB01.1	GenBank
' <i>M. orygis</i> '	112400015	APKD01.1	GenBank
<i>M. pseudoshottsii</i>	L15 ^T	BCND01.1	GenBank
<i>Nocardia asteroides</i>	NBRC 15531 ^T	BAFO02.1	GenBank



Research Results and
Conclusions



Genomic Comparison of MTBC Type Strains



Table 2. Genomic Distance Analysis of type strains of members of the MTBC

Pairwise genome-to-genome distances (GGDs) for the type strains of the species within the *Mycobacterium tuberculosis* complex ('MTBC') and two outgroups ('OUT'). Results from digital DNA–DNA hybridization (dDDH) are shown above the self-comparison diagonal, and average nucleotide identity (ANI) results are shown below the diagonal. Species delineation thresholds used were 80 % for dDDH and 96 % for ANI.

	Species/ Strain	dDDH	MTBC1	MTBC2	MTBC3	MTBC4	MTBC5	MTBC6	MTBC7	MTBC8	MTBC9	OUT1	OUT2
MTBC1	<i>M. tuberculosis</i> H37Rv ^T	ANI	100	97.7	97.5	97.9	98.7	97.3	91.2	97.9	97.6	21.1	18.8
MTBC2	<i>M. africanum</i> ATCC® 25420 ^T		99.88	100	98	98.5	98.9	98.2	92.2	98.9	98.4	22.2	18.8
MTBC3	<i>M. bovis</i> ATCC® 19210 ^T		99.78	99.82	100	98.7	98.2	97.8	91.6	98.1	97.8	23.3	19
MTBC4	<i>M. caprae</i> ATCC® BAA-824 ^T		99.86	99.87	99.83	100	98.8	97.7	91.7	98.6	98.2	21.1	18.8
MTBC5	<i>M. microti</i> ATCC® 19422 ^T		99.89	99.8	99.79	99.91	100	99.2	93.3	99	98.9	23.3	18.9
MTBC6	<i>M. pinnipedii</i> ATCC® BAA-688 ^T		99.84	99.86	99.78	99.84	99.92	100	91.6	98.3	97.6	21.1	18.8
MTBC7	" <i>M. canettii</i> " CIPT 140010059		99.25	99.28	99.21	99.26	99.36	99.26	100	92.5	91.8	22.2	19
MTBC8	" <i>M. mungi</i> " BM22813		99.88	99.91	99.81	99.88	99.92	99.89	99.31	100	98.6	22.2	18.8
MTBC9	" <i>M. orygis</i> " 112400015		99.85	99.87	99.80	99.87	99.91	99.86	99.28	99.88	100	21.1	18.8
OUT1	<i>M. pseudoshottsii</i> L15 ^T		79.33	79.27	79.21	79.34	79.34	79.25	79.29	79.25	79.37	100	19.2
OUT2	<i>Nocardia asteroides</i> NBRC 15531 ^T		70.92	70.76	70.96	70.94	70.88	71.02	70.99	71.00	70.84	70.75	100

Color Key

ANI	dDDH
Self-Comparison	Self-Comparison
96 – 100	80 – 100
< 96	< 80

Genomic Comparison of MTBC Type Strains

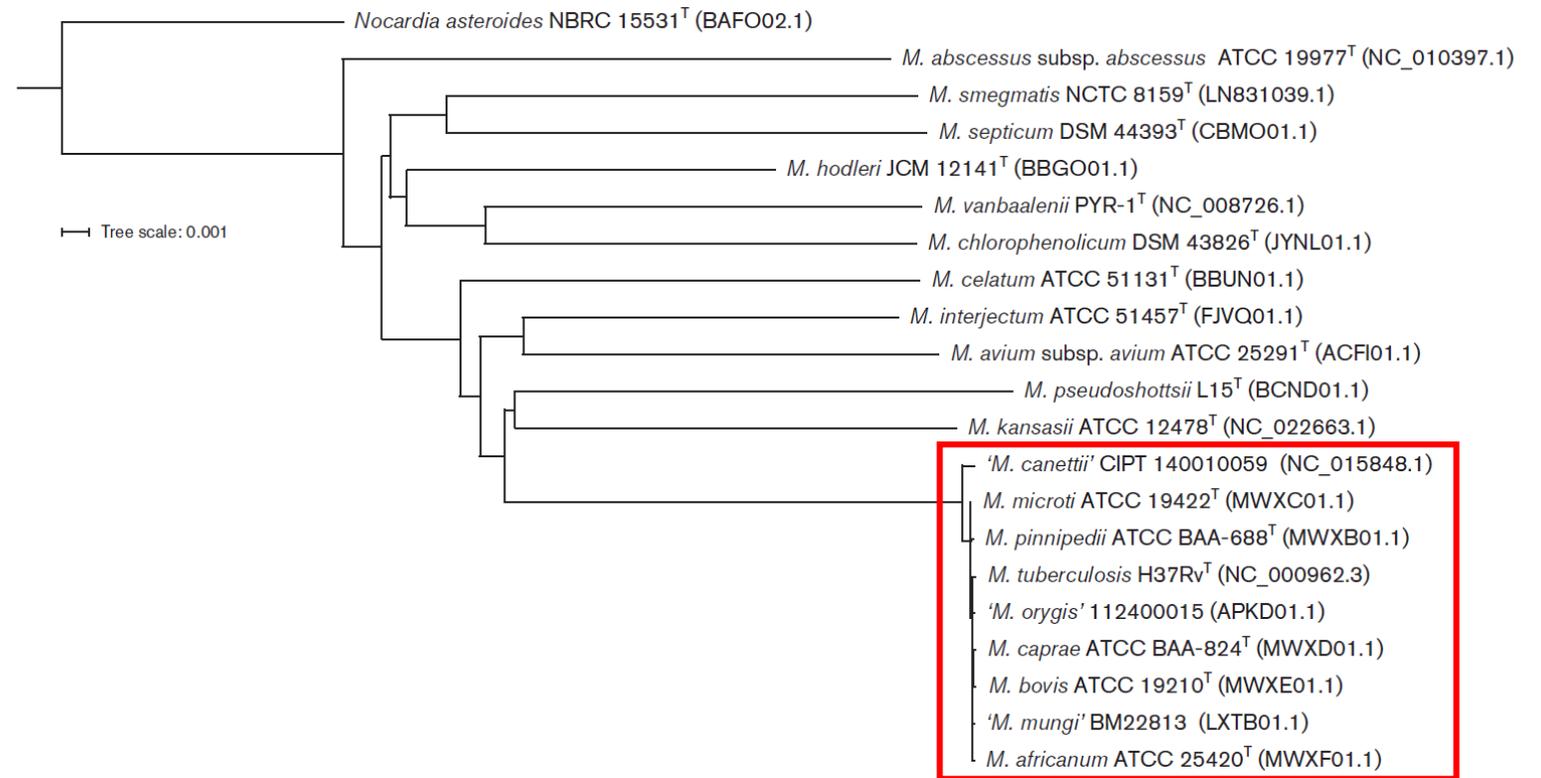


Fig. 1. Phylogenomic tree showing the relationship between the whole genomes of type strains of species of MTBC and the type strains of various other species of the genus *Mycobacterium* with *Nocardia asteroides* as an outgroup. Bar, 0.001 substitutions per site.

Genomic Comparison of MTBC Type Strains

dDDH GGD values are between the GenBank genomes identified as respective species and the type strain of *M. tuberculosis* (H37Rv^T).

GenBank organism identifier	<i>n</i>	dDDH GGD to <i>M. tuberculosis</i> , H37Rv ^T (%)		
		Minimum	Mean	Maximum
<i>M. africanum</i>	30	96.7	97.5	98.3
<i>M. bovis</i>	69	95.7	97.8	99.1
' <i>M. canettii</i> '	9	80.1	89.8	94.1
<i>M. caprae</i>	2	97.4	97.7	97.9
<i>M. microti</i>	1	97.1	97.1	97.1
<i>M. tuberculosis</i>	3631*	83.5	98.6	100.0

*Excludes one single outlier which was examined separately (see text and Table S8).

↳ dDDH: 22.2%, ANI: 79.39%

Genomic Comparison of MTBC Type Strains



Table S8. dDDH GGDs Between *M. tuberculosis* TTK-01-0051 in GenBank vs. *M. tuberculosis* H37Rv[†]

Query genome	Reference genome	Formula 2			
		DDH	Model C.I.	Distance	Prob. DDH >= 70%
JLXW01.1_Mtuberculosis	TS1_NC_010397.1_M_abscessus_abscessus	21.2	[18.9 - 23.6%]	0.2072	0
JLXW01.1_Mtuberculosis	TS2_AHAS01.1_M_abscessus_bollettii	19.5	[17.3 - 21.9%]	0.2252	0
JLXW01.1_Mtuberculosis	TS3_GH1_Mafricanum	22.1	[19.9 - 24.6%]	0.1982	0
JLXW01.1_Mtuberculosis	TS6_GH1_Malgericum	20.5	[18.3 - 23%]	0.2138	0
JLXW01.1_Mtuberculosis	TS11_JALN02.1_M_aromaticivorans	20.1	[17.9 - 22.6%]	0.2182	0
JLXW01.1_Mtuberculosis	TS14_CCB01.1_M_asiaticum	21.8	[19.5 - 24.2%]	0.2014	0
JLXW01.1_Mtuberculosis	TS15_GH1_Maubagnense	19.4	[17.2 - 21.8%]	0.2263	0
JLXW01.1_Mtuberculosis	TS16_CVQ01.1_M_aurum	19.7	[17.5 - 22.1%]	0.223	0
JLXW01.1_Mtuberculosis	TS17_CCAW01.1_M_austroafricanum	21.2	[18.9 - 23.6%]	0.2074	0
JLXW01.1_Mtuberculosis	TS18_ACFI01.1_M_avium_avium	30.3	[27.9 - 32.8%]	0.1406	0.12
JLXW01.1_Mtuberculosis	TS19_AGAR01.1_M_avium_paratuberculosis	30.7	[28.3 - 33.2%]	0.1385	0.14
JLXW01.1_Mtuberculosis	TS20_AYOC01.1_M_avium_silvaticum	30.7	[28.4 - 33.3%]	0.1381	0.14
JLXW01.1_Mtuberculosis	TS23_CSTD01.1_M_bohemicum	24.8	[22.5 - 27.3%]	0.1755	0.01
JLXW01.1_Mtuberculosis	TS27_GH1_Mbovis	22.1	[19.9 - 24.6%]	0.198	0
JLXW01.1_Mtuberculosis	TS30_FJNX01.1_M_brumae	20.4	[18.2 - 22.9%]	0.2149	0
JLXW01.1_Mtuberculosis	TS31_BCSY01.1_M_canariense	20.5	[18.3 - 22.9%]	0.2145	0
JLXW01.1_Mtuberculosis	TS32_GH1_Mcaprae	22	[19.8 - 24.5%]	0.1992	0
JLXW01.1_Mtuberculosis	TS33_BBUN01.1_M_celatum	22.5	[20.2 - 24.9%]	0.195	0
JLXW01.1_Mtuberculosis	TS35_CP010946.1_M_chelonae	19.4	[17.3 - 21.8%]	0.226	0
JLXW01.1_Mtuberculosis	TS36_MRBR01.1_M_chimaera	31.5	[29.1 - 34.1%]	0.134	0.2
JLXW01.1_Mtuberculosis	TS38_JYNL01.1_M_chlorophenolicum	20.8	[18.5 - 23.2%]	0.2116	0
JLXW01.1_Mtuberculosis	TS39_NC_018027.1_M_chubuense	20	[17.8 - 22.4%]	0.2194	0
JLXW01.1_Mtuberculosis	TS40_AFWW02.1_M_colombiense	81.3	[78.4 - 83.9%]	0.0219	91.52
JLXW01.1_Mtuberculosis	TS41_CTEF01.1_M_conceptionense	21.1	[18.8 - 23.5%]	0.2085	0
JLXW01.1_Mtuberculosis	TS45_CCB01.1_M_cosmeticum	20.9	[18.7 - 23.3%]	0.21	0
JLXW01.1_Mtuberculosis	TS46_BBHD01.1_M_crocinum	22	[19.7 - 24.4%]	0.1996	0
JLXW01.1_Mtuberculosis	TS50_GH1_Melephantis	22.7	[20.4 - 25.2%]	0.1929	0
JLXW01.1_Mtuberculosis	TS54_CCA01.1_M_farcinogenes	20	[17.8 - 22.4%]	0.2193	0
JLXW01.1_Mtuberculosis	TS57_BBFT01.1_M_fluoranthenvivorans	21.6	[19.3 - 24%]	0.2033	0
JLXW01.1_Mtuberculosis	TS58_CP014258.1_M_fortuitum_fortuitum	19.8	[17.6 - 22.2%]	0.2217	0
JLXW01.1_Mtuberculosis	TS59_BCSZ01.1_M_fortuitum_acetamidolyticum	20.7	[18.5 - 23.1%]	0.212	0
JLXW01.1_Mtuberculosis	TS62_GH1_Mfrederiksbergense	20.9	[18.6 - 23.3%]	0.2103	0
JLXW01.1_Mtuberculosis	TS64_AZYN01.1_M_gastri	22.5	[20.2 - 24.9%]	0.1948	0
JLXW01.1_Mtuberculosis	TS65_JAGZ01.1_M_genavense	23.7	[21.4 - 26.2%]	0.1843	0
JLXW01.1_Mtuberculosis	TS69_CP011883.2_M_haemophilum	22.4	[20.1 - 24.8%]	0.1957	0

Color Key	
dDDH Range	Interpretation [†]
80.0 - 100	Same species and subspecies
70.0 - 79.9	Same species, different subspecies
50.0 - 69.9	Different species
30.0 - 49.9	
0.0 - 29.9	

[†]Meier-Kolthoff *et al.*, 2014

(ANI: 97.73%)

Unification of MTBC Species as *M. tuberculosis*

- Unites the MTBC as *Mycobacterium tuberculosis*
- Submitted emended species description in *IJSEM*
 - Type strain: *M. tuberculosis* H37Rv^T

Emended description of *Mycobacterium tuberculosis* (Zopf 1883) Lehmann and Neumann 1896 (Approved Lists 1980)

M. africanum, *M. bovis*, *M. caprae*, *M. microti* and *M. pinnipedii* are reclassified as later heterotypic synonyms of *M. tuberculosis*. The strains described with the effectively but not validly published names '*M. canettii*' [6], '*M. mungi*' [7], and '*M. orygis*' [8] are likewise reclassified as belonging to the species *M. tuberculosis*. The phylogenetic groups that correspond to the previously named species should be considered infrasubspecific subdivisions, i.e. variants. Thus, the characteristics described for the above previously named species are now included in the description of *M. tuberculosis*. Although infrasubspecific subdivisions are not governed by the rules of prokaryotic nomenclature, it is recommended that these infrasubspecific subdivisions retain the previous specific epithet as the infrasubspecific designations, e.g. *M. tuberculosis* var. *bovis* (see Table 4).

Unification of MTBC Species as *M. tuberculosis*

- Unites the MTBC as *Mycobacterium tuberculosis*
- Submitted emended species description in *IJSEM*
 - Type strain: *M. tuberculosis* H37Rv^T
 - All other species type strains downgraded to reference strains
 - Recommendation: use former species name as intrasubspecific designation

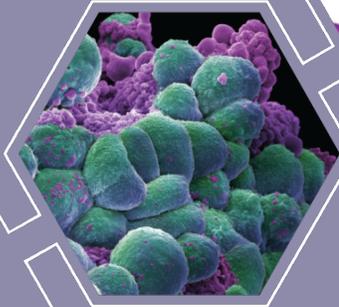
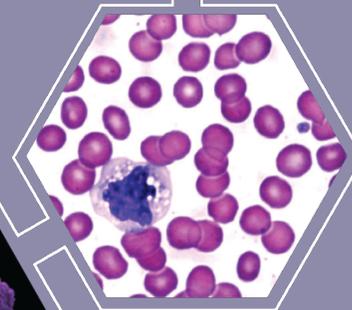
Table 4. Recommended infrasubspecific designations and reference strains

Current Name	Recommended Name	Reference Strain
<i>M. tuberculosis</i>	<i>M. tuberculosis</i> var. <i>tuberculosis</i>	H37Rv ^T (ATCC 27294 ^T) (type strain of species)
<i>M. africanum</i>	<i>M. tuberculosis</i> var. <i>africanum</i>	ATCC 25420
<i>M. bovis</i>	<i>M. tuberculosis</i> var. <i>bovis</i>	ATCC 19210
<i>M. bovis</i> BCG	<i>M. tuberculosis</i> var. BCG	–
<i>M. caprae</i>	<i>M. tuberculosis</i> var. <i>caprae</i>	ATCC BAA-824
<i>M. microti</i>	<i>M. tuberculosis</i> var. <i>microti</i>	ATCC 19422
<i>M. pinnipedii</i>	<i>M. tuberculosis</i> var. <i>pinnipedii</i>	ATCC BAA-688
' <i>M. canettii</i> '	<i>M. tuberculosis</i> var. <i>canettii</i>	CIPT 140010059
' <i>M. mungi</i> '	<i>M. tuberculosis</i> var. <i>mungi</i>	BM22813
' <i>M. orygis</i> '	<i>M. tuberculosis</i> var. <i>orygis</i>	112400015
' <i>M. suricattae</i> '*	<i>M. tuberculosis</i> var. <i>suricattae</i>	–
Dassie bacillus*	<i>M. tuberculosis</i> var. <i>dassie</i>	–
Chimpanzee bacillus*	<i>M. tuberculosis</i> var. <i>chimpanzee</i>	–

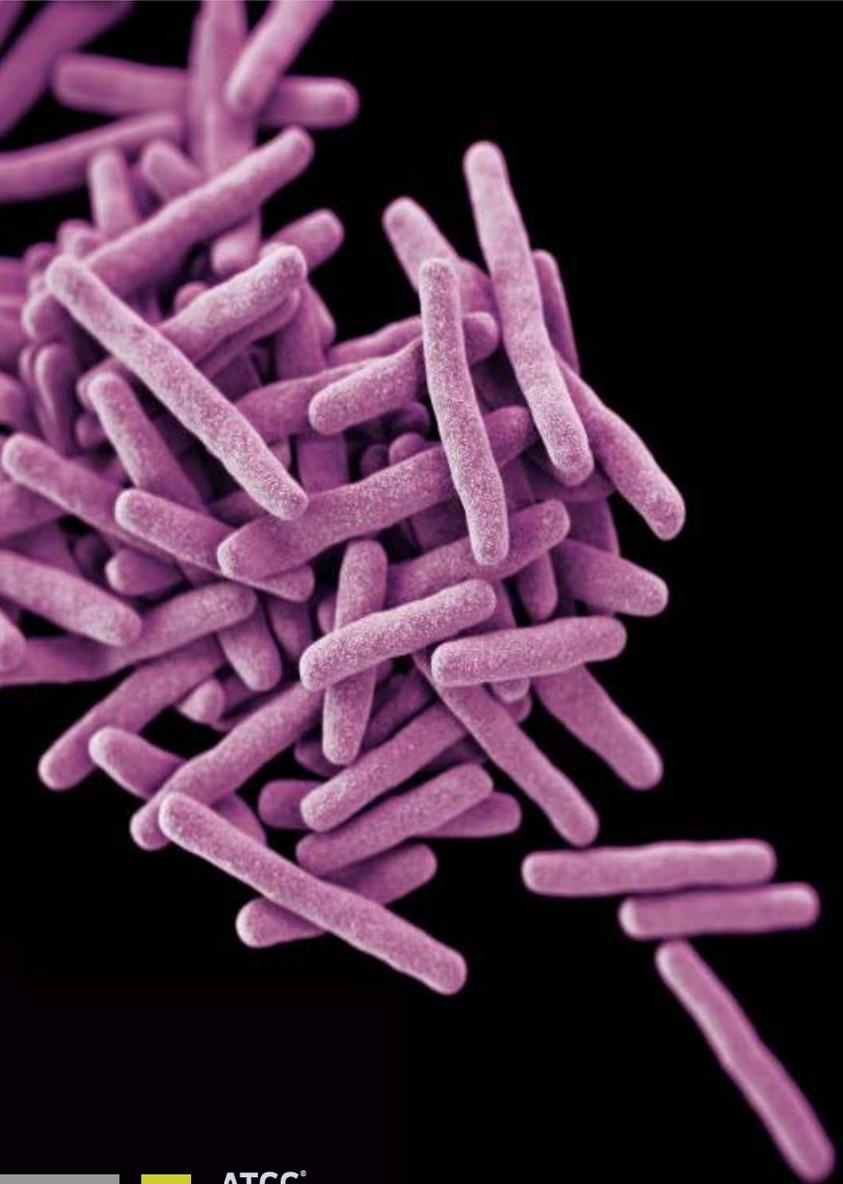
*These organisms were not included in this study, but recommended infrasubspecific designations (based upon the typically used nomenclature for these organisms) are included for completeness. Note that the strains that were previously listed as the species type strains are now considered reference strains of the respective varieties.



Conclusions



Conclusions



- Phylogenomic analysis comparing the type strains of the “species” within the MTBC shows that they are, in fact, a single species
- dDDH comparison against a larger set of sequenced type strains enabled discovery of the identity of a strain misclassified as *M. tuberculosis*
- Whole-genome sequencing of bacterial type strains enables the elucidation of the true relationships between species and strains
 - Modernize taxonomic relationships using complete genetic information
 - Develop detailed phylogenomic trees
 - Determine the identity of unknown sequenced samples

References

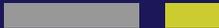
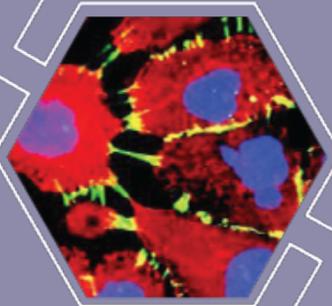
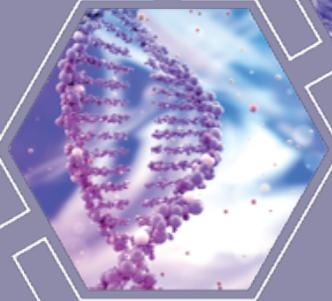
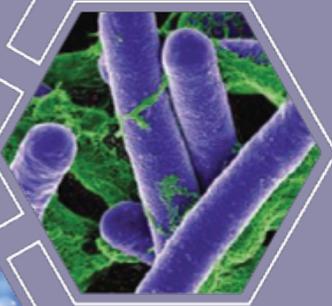


- **Riojas MA, et al.** Phylogenomic analysis of the species of the *Mycobacterium tuberculosis* complex demonstrates that *Mycobacterium africanum*, *Mycobacterium bovis*, *Mycobacterium caprae*, *Mycobacterium microti* and *Mycobacterium pinnipedii* are later heterotypic synonyms of *Mycobacterium tuberculosis*. *Int J Syst Evol Microbiol* 68: 324–332, 2018.
- **Auch AF, von Jan M, Klenk HP & Göker M.** Digital DNA-DNA hybridization for microbial species delineation by means of genome-to-genome sequence comparison. *Stand Genomic Sci* 2: 117-134, 2010.
- **Auch AF, Klenk HP & Göker M.** Standard operating procedure for calculating genome-to-genome distances based on high-scoring segment pairs. *Stand Genomic Sci* 2: 142-148, 2010.
- **Garcia-Betancur JC, et al.** Alignment of multiple complete genomes suggests that gene rearrangements may contribute towards the speciation of Mycobacteria. *Infect Genet Evol* 12: 819–826, 2012.
- **Meier-Kolthoff JP, et al.** Complete genome sequence of DSM 30083^T, the type strain (U5/41^T) of *Escherichia coli*, and a proposal for delineating subspecies in microbial taxonomy. *Stand Genomic Sci* 9:2, 2014.
- **Meier-Kolthoff JP, et al.** Genome sequence-based species delimitation with confidence intervals and improved distance functions. *BMC Bioinformatics* 14:60, 2013.

Disclaimers



© American Type Culture Collection. The ATCC trademark and trade name, and any other trademarks listed in this publication are trademarks owned by the American Type Culture Collection unless indicated otherwise.



Thank You

Register for more ATCC webinars at
www.atcc.org/Webinars

