

Carbapenem-resistant Enterobacteriaceae (CRE) – A Growing Superbug Population

Cara Wilder, Ph.D.
Technical Writer, ATCC
May 5, 2016



About ATCC

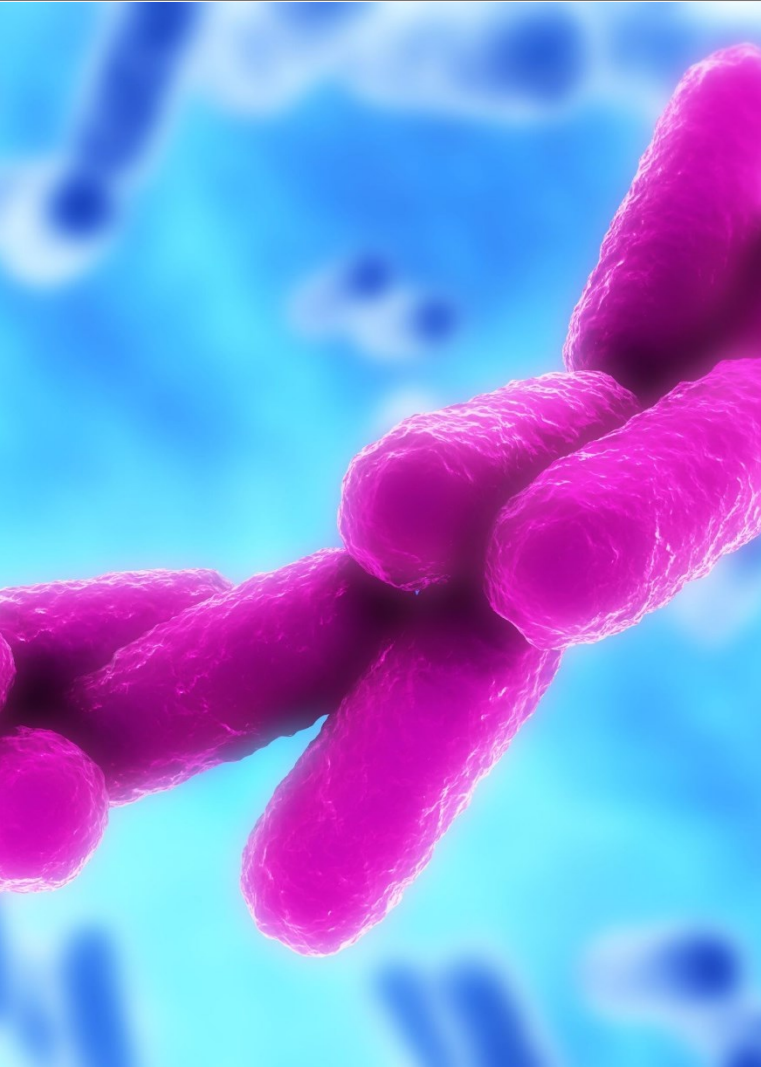
- Founded in 1925, ATCC is a non-profit organization with headquarters in Manassas, VA
- World's premiere biological materials resource and standards development organization
- ATCC collaborates with and supports the scientific community with industry-standard biological products and innovative solutions
- Strong team of 400+ employees; over one third with advanced degrees



Established partner to global researchers and scientists



Outline



1. Background on antibiotic resistance
2. Emergence and spread of carbapenem-resistant Enterobacteriaceae (CRE)
3. CRE strains available from ATCC
 - KPC
 - NDM
 - OXA-48

Multidrug Resistance (MDR) is an Emerging Threat

- Antimicrobial resistance is present in all parts of the world
- The CDC estimates that every year in the United States:
 - **2 million** people become infected with antibiotic-resistant bacteria
 - **23,000** people die as a direct result of these infections
 - **\$20 billion** in excess direct healthcare costs
 - **\$35 billion** cost associated with lost productivity



Bad Bugs, No Drugs: No ESKAPE!



Enterococcus faecium

Staphylococcus aureus

Klebsiella sp.

Acinetobacter baumannii

Pseudomonas aeruginosa

Enterobacter sp.



These pathogens cause the majority of US hospital infections and can effectively “escape” the effects of antimicrobial therapeutics

Multidrug Resistance is Widespread

Minimum Estimates of Morbidity and Mortality from Antibiotic-Resistant Infections

Antibiotic-Resistant Microorganism	Estimated Annual Number of Cases	Estimated Annual Number of Deaths
Carbapenem-resistant Enterobacteriaceae (CRE)	9,300	610
Extended-spectrum β -lactamase producing Enterobacteriaceae (ESBL)	26,000	1,700
Vancomycin-resistant <i>Enterococcus</i> (VRE)	20,000	1,300
Methicillin-resistant <i>Staphylococcus aureus</i> (MRSA)	80,000	11,000
Multidrug-resistant <i>Acinetobacter</i> (≥ 3 drug classes)	7,300	500
Multidrug-resistant <i>Pseudomonas aeruginosa</i> (≥ 3 drug classes)	6,700	440

National Action Plan for Combating Antibiotic-Resistant Bacteria

The White House – March 2015

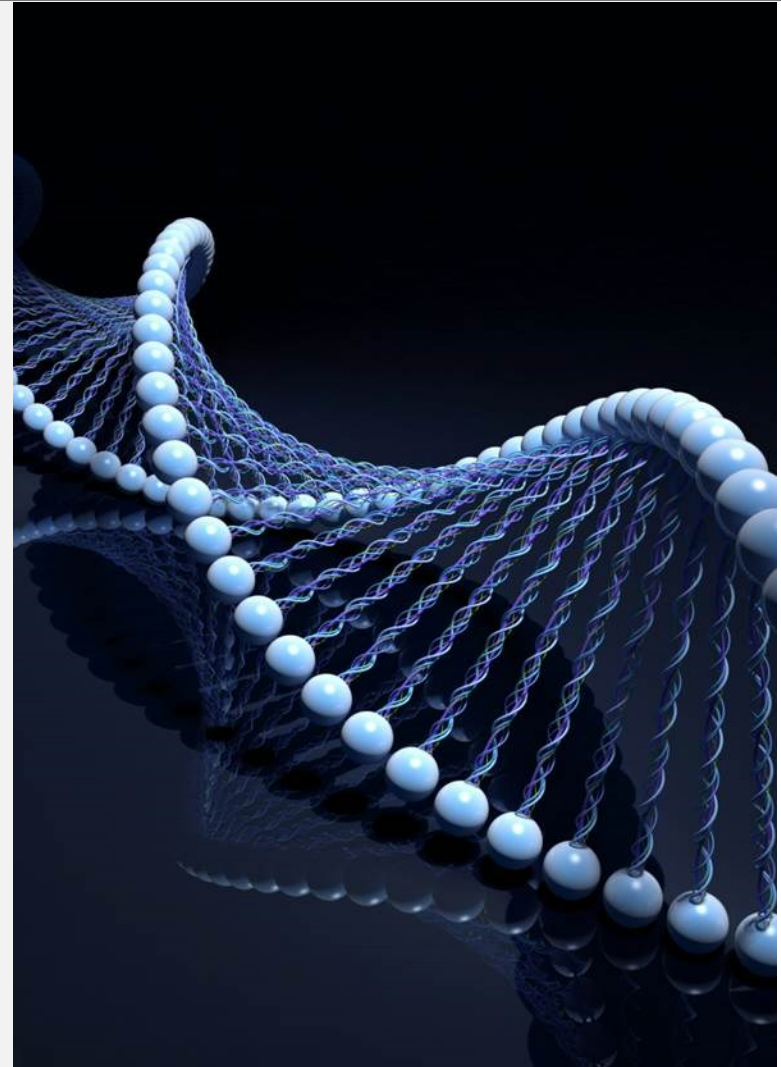
Goals:

- Slow the emergence of resistant bacteria and prevent the spread of resistant infections
- Strengthen national One Health surveillance efforts to combat resistance
- Advance development and use of rapid and innovative diagnostic tests for identification and characterization of resistant bacteria
- Accelerate basic and applied research and development for new antibiotics, other therapeutics, and vaccines
- Improve international collaboration and capacities for antibiotic-resistance prevention, surveillance, control, and antibiotic research and development



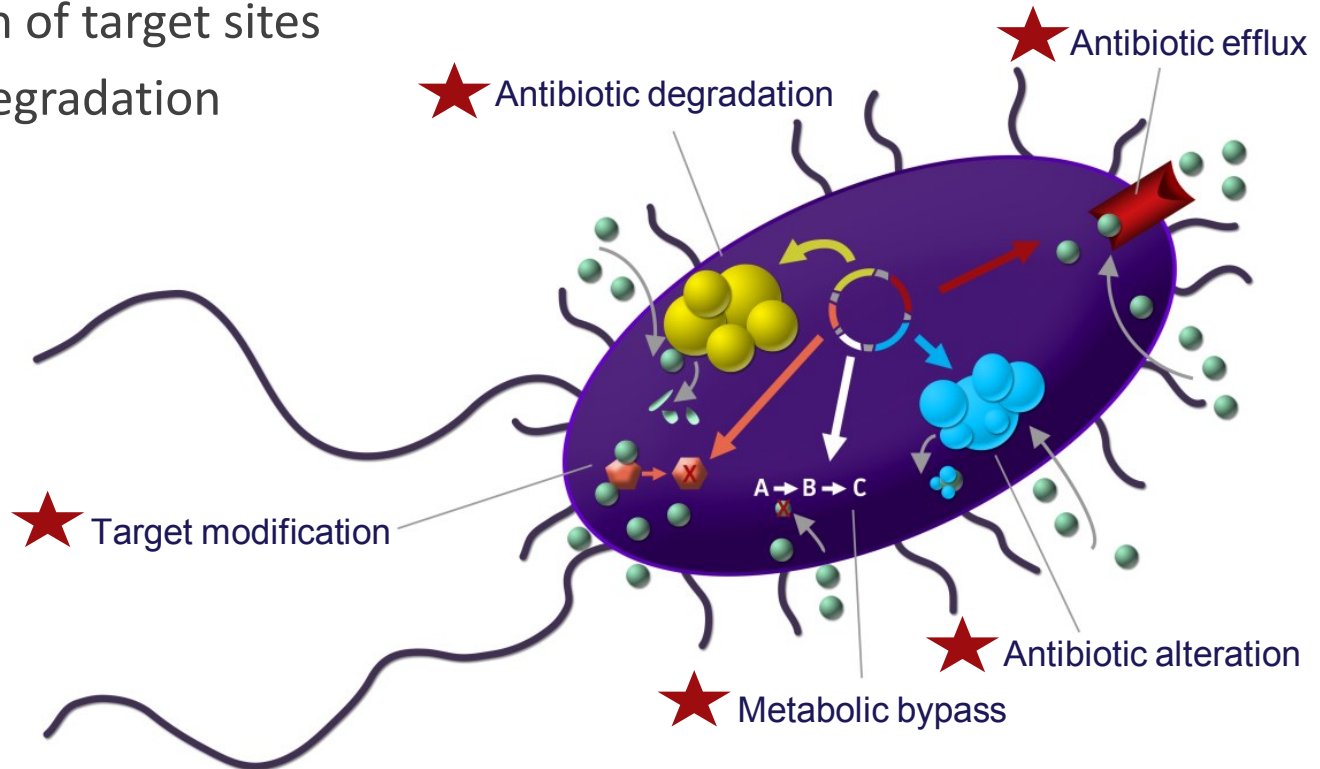
Antibiotic Resistance – Evolution & Dissemination

- Inherent resistance
- Genetic mutation
- Horizontal gene transfer
 - Transformation
 - Transduction
 - Conjugation

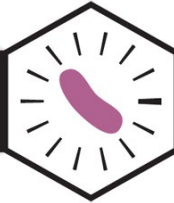


Antibiotic Resistance – Evolution & Dissemination

- Reduced drug accumulation
- Antibiotic alteration
- Metabolic bypass
- Modification of target sites
- Antibiotic degradation



Antibiotic Resistance – Evolution & Dissemination



How Antibiotic Resistance Happens

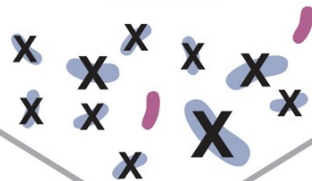
1.

Lots of germs.
A few are drug resistant.



2.

Antibiotics kill
bacteria causing the illness,
as well as good bacteria
protecting the body from
infection.



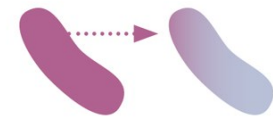
3.

The drug-resistant
bacteria are now allowed to
grow and take over.

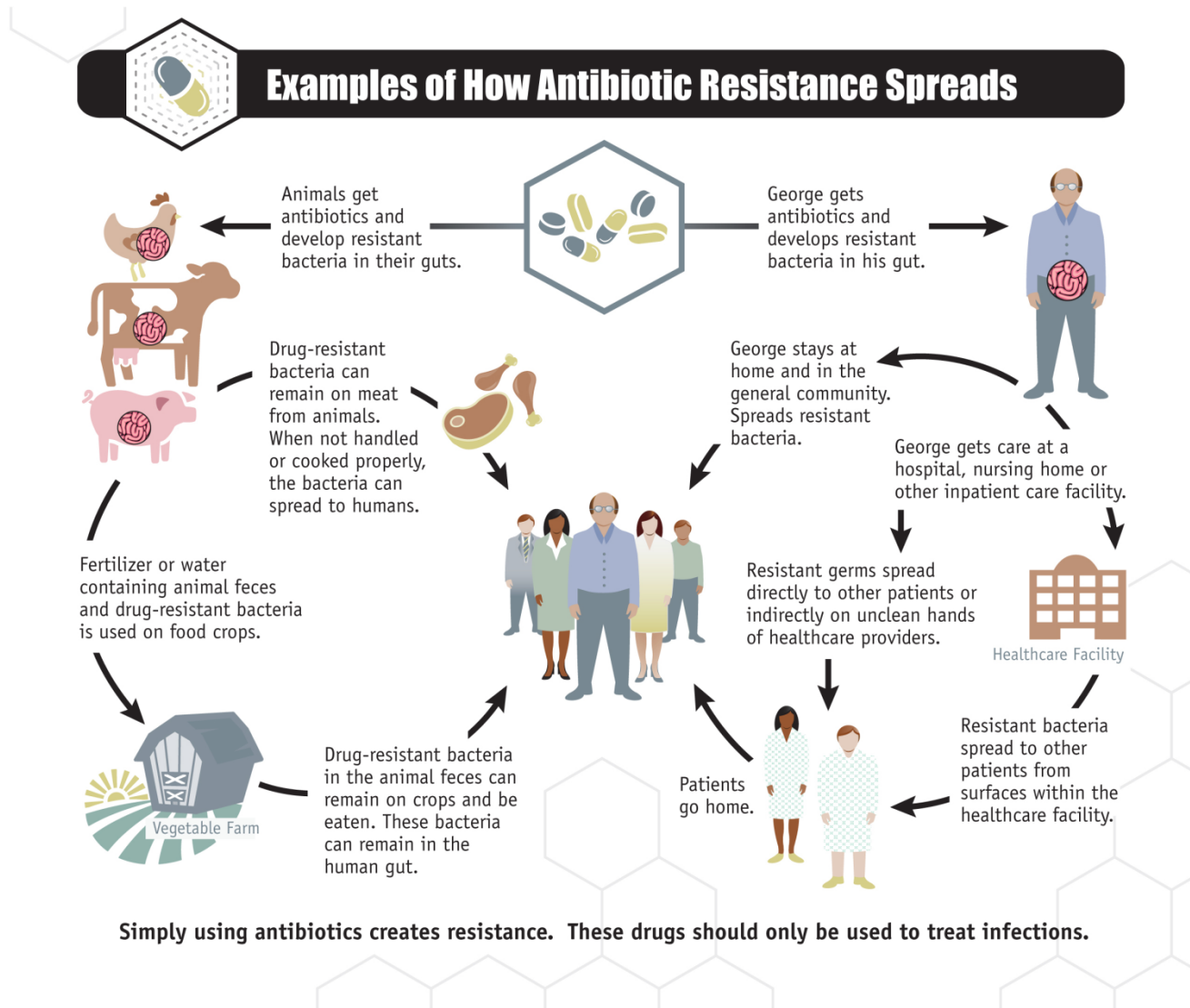


4.

Some bacteria give
their drug-resistance to
other bacteria, causing
more problems.



Antibiotic Resistance – Evolution & Dissemination



CRE – Hazard Level Urgent

HAZARD LEVEL
URGENT



These are high-consequence antibiotic-resistant threats because of significant risks identified across several criteria. These threats may not be currently widespread but have the potential to become so and require urgent public health attention to identify infections and to limit transmission.



THREAT LEVEL
URGENT



This bacteria is an immediate public health threat that requires urgent and aggressive action.

CARBAPENEM-RESISTANT ENTEROBACTERIACEAE



9,000

DRUG-RESISTANT
INFECTIONS
PER YEAR



600

DEATHS

CARBAPENEM-
RESISTANT
KLEBSIELLA SPP.

7,900



1,400

CARBAPENEM-
RESISTANT
E. COLI



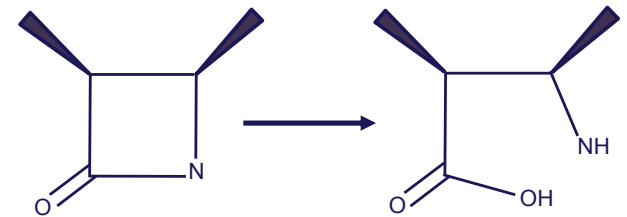
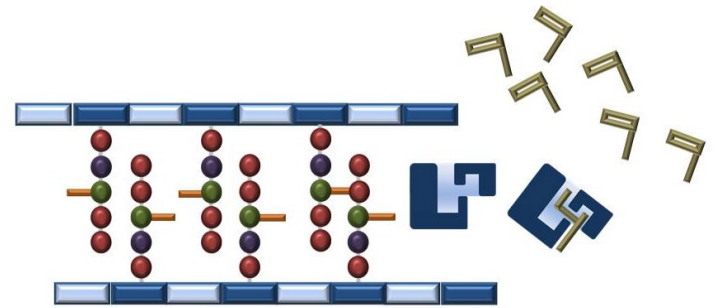
**CRE HAVE BECOME RESISTANT TO ALL
OR NEARLY ALL AVAILABLE ANTIBIOTICS**



Carbapenem Resistance

- Carbapenem antibiotics
 - β -lactam antibiotic
 - Inhibit peptidoglycan synthesis

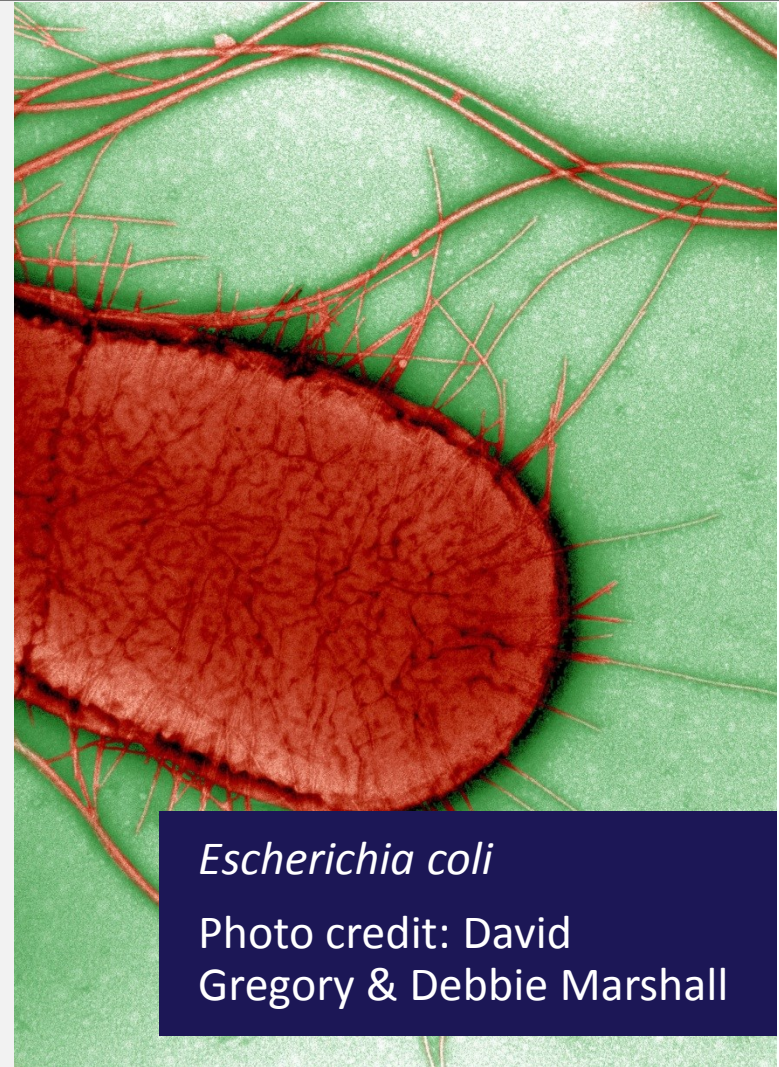
- Mechanisms of carbapenem-resistance
 - β -lactamase production combined with porin mutations
 - Carbapenemase production



CRE Definitions and Recommendations

CDC has defined CRE as Enterobacteriaceae that are:

- Resistant to any carbapenem antimicrobial (i.e. MIC of $\geq 4 \mu\text{g/mL}$ for doripenem, meropenem, or imipenem OR $\geq 2 \mu\text{g/mL}$ for ertapenem)
- Documented to produce carbapenemase



CRE Definitions and Recommendations

- Lower CLSI break points allow easier detection
- More information is available:
 - CLSI M100-S25 2015
 - CDC 2015 CRE Toolkit
 - AHRQ Carbapenem-Resistant *Enterobacteriaceae* (CRE) Control and Prevention Toolkit

Agent	Previous Breakpoints (M100-S19) MIC (µg/mL)			Current Breakpoints (M100-S25) MIC (µg/mL)		
	Susceptible	Intermediate	Resistant	Susceptible	Intermediate	Resistant
Doripenem	--	--	--	≤ 1	2	≥ 4
Ertapenem	≤ 2	4	≥ 8	≤ 0.5	1	≥ 2
Imipenem	≤ 4	8	≥ 16	≤ 1	2	≥ 4
Meropenem	≤ 4	8	≥ 16	≤ 1	2	≥ 4

Concerns About CRE as an Emerging Threat



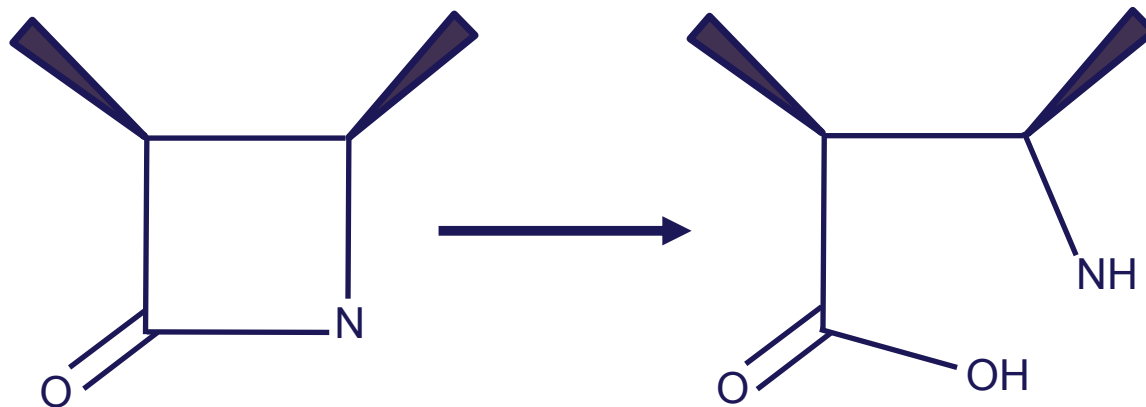
Klebsiella pneumoniae

Photo credit: NIAID

- Multidrug-resistant
- High mortality rates for invasive infections
 - Up to 50% in some studies
- Rapid spread in healthcare settings
- Potential to become widespread in the community
- Carbapenemase genes can be transmitted from one Enterobacteriaceae to another
- Increase in CRE strains
 - 1.2% in 2001 to 4.2% in 2011

CRE strains

Carbapenemase	Ambler Class	Known Bacterial Carriers
KPC	A	<i>K. pneumoniae</i> , <i>E. coli</i> , and <i>Enterobacter</i> spp.
NDM	B	<i>E. coli</i> , <i>K. pneumoniae</i> , and <i>E. cloacae</i>
VIM	B	<i>P. aeruginosa</i> and <i>P. putida</i>
IMP	B	<i>Pseudomonas</i> and <i>Acinetobacter</i> spp.
CMY	C	<i>E. aerogenes</i>
OXA	D	<i>K. pneumoniae</i> , <i>E. coli</i> , <i>Acinetobacter</i> spp.



CRE strains

Carbapenemase	Ambler Class	Known Bacterial Carriers
KPC	A	<i>K. pneumoniae</i> , <i>E. coli</i> , and <i>Enterobacter</i> spp.
NDM	B	<i>E. coli</i> , <i>K. pneumoniae</i> , and <i>E. cloacae</i>
VIM	B	<i>P. aeruginosa</i> and <i>P. putida</i>
IMP	B	<i>Pseudomonas</i> and <i>Acinetobacter</i> spp.
CMY	C	<i>E. aerogenes</i>
OXA	D	<i>K. pneumoniae</i> , <i>E. coli</i> , <i>Acinetobacter</i> spp.



DEADLY DRUG-RESISTANT BUG
C.R.E. HAS A DEATH RATE OF ABOUT 50%

By Peter Coates (H)
A NEW "superbug" has been found in two patients here, one of whom died. The bug is a type of carbapenem-resistant enterobacter (CRE) that is highly resistant to antibiotics. It is a type of bacteria that is often found in hospitals and is a common cause of infection in patients who have been in a hospital. The bug is a type of bacteria that is often found in hospitals and is a common cause of infection in patients who have been in a hospital.

New superbug found in two patients here

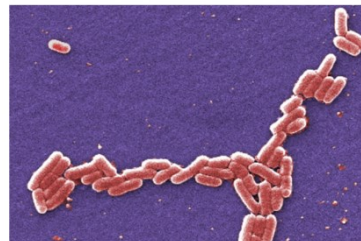
Both cases successfully contained an expert, large health officials to track deadly bacteria

The state health department confirmed the presence of a new strain of carbapenem-resistant enterobacter (CRE) in two patients here. The bug is a type of bacteria that is often found in hospitals and is a common cause of infection in patients who have been in a hospital.

The bug is a type of bacteria that is often found in hospitals and is a common cause of infection in patients who have been in a hospital. The bug is a type of bacteria that is often found in hospitals and is a common cause of infection in patients who have been in a hospital.

Last-Ditch Resistance: More Countries, More Dire Results

POSTED THU, 01/17/2015



An E. coli bacterium. PHOTOGRAPH BY THE PUBLIC HEALTH SERVICE, CDC/DOH

The frantic international hunt triggered by the discovery of genetically mobile resistance to colistin, a last-resort antibiotic, is producing many more findings this evening. The resistance factor is showing up in more countries, but, much

'Phantom Menace' Superbug Has Alarming Ability

Antibiotic resistance travels between bacteria

By Rob Quinn, Newser Staff
Posted Dec 4, 2013 5:50 AM CST



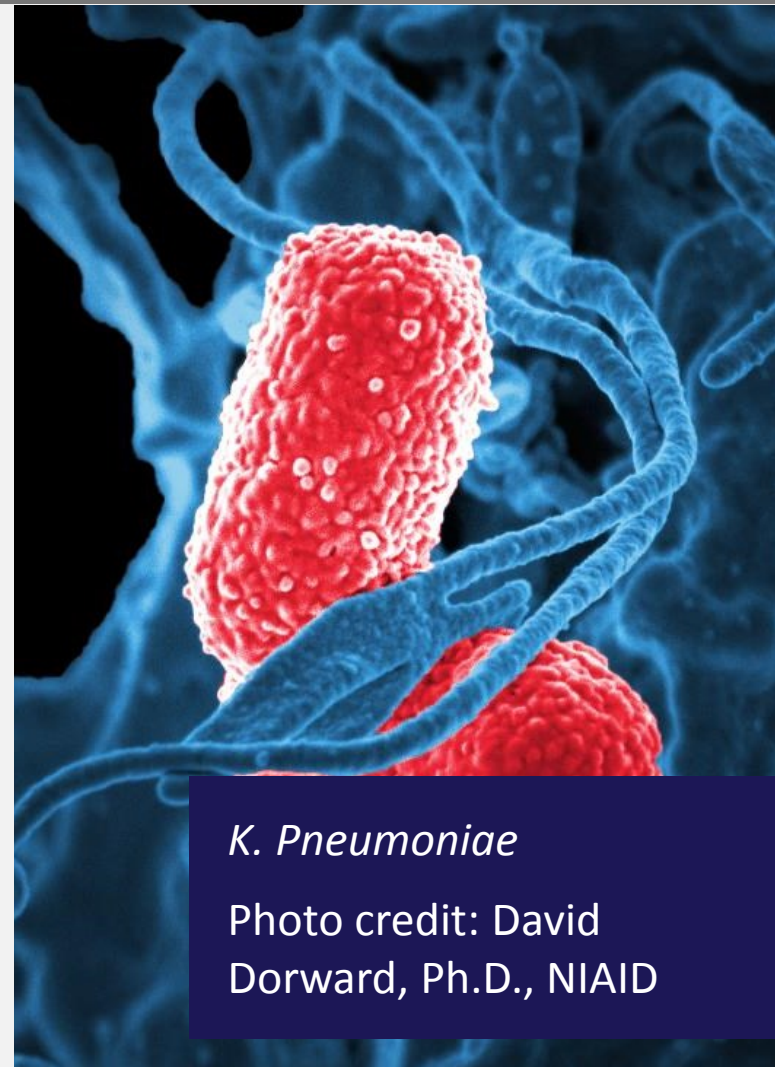
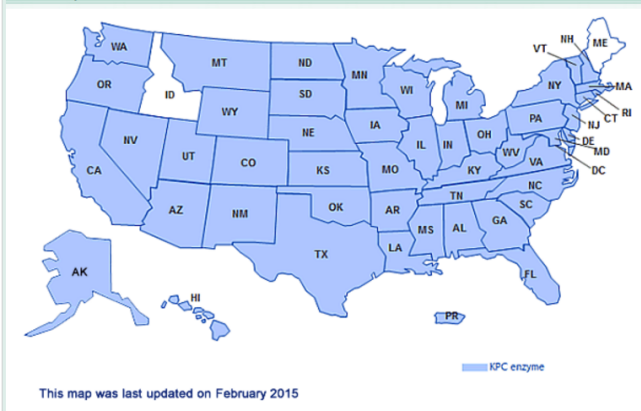
This illustration released by the CDC depicts a group of carbapenem-resistant Enterobacteriaceae bacteria. (AP Photo/Centers for Disease Control)

(NEWSER) – The latest strains of antibiotic-resistant bacteria to give researchers sleepless nights aren't the most common—or even the most antibiotic-resistant—but they have an ability that could make them a serious danger to public health. They contain enzymes known as "OXA-48-like carbapenemases" that can break down antibiotics and transfer that ability to normal bacteria in the body through mobile pieces of DNA, reports the *Washington Post* and *Live Science*. The enzymes have been nicknamed the "phantom menace" by researchers because they can be tough to detect. In a report issued this week, the CDC says it identified 43 cases in the US involving the superbug between June 2010 and August of this year, mainly involving people who had

Klebsiella pneumoniae Carbapenemase (KPC)

- KPC hydrolyzes all β -lactam agents
- Encoded by the plasmid-associated gene *bla*_{KPC}
- May be difficult to detect using higher (older) breakpoints
- The CDC has confirmed the presence of KPC throughout most of the United States

States with KPC-producing Carbapenem-resistant Enterobacteriaceae (CRE) reported to the Centers for Disease Control and Prevention (CDC) as of February 2015



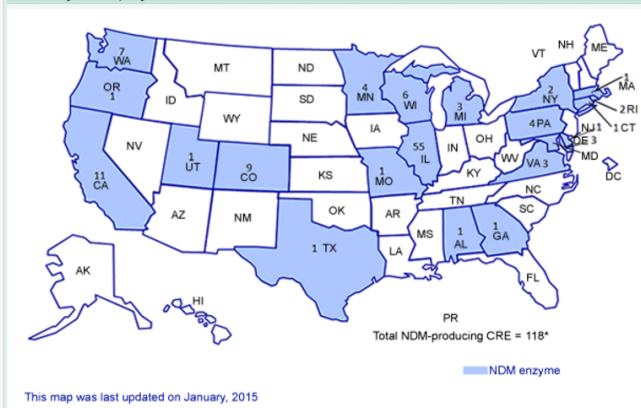
K. Pneumoniae

Photo credit: David Dorward, Ph.D., NIAID

New Delhi Metallo- β -lactamase (NDM)

- Encoded by the plasmid-associated gene *bla*_{NDM}
- First identified in 2008
- It has since been detected worldwide

NDM-producing Carbapenem-resistant Enterobacteriaceae (CRE) isolates reported to the Centers for Disease Control and Prevention (CDC) as of January 2015, by state

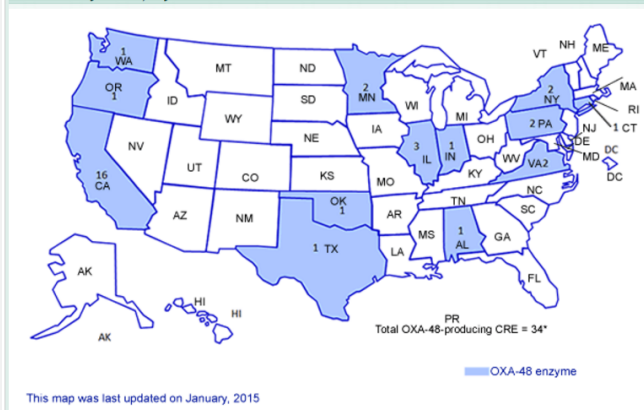


Enterobacteriaceae
Photo credit: CDC

Oxacillinase-48 (OXA-48)

- Penicillinases that can hydrolyze penicillins and imipenem
- Encoded by the plasmid-associated gene *bla*_{OXA-48}
- First isolated in 2001 in Istanbul, Turkey, has since spread in Europe, the Middle East, India, North Africa, and the US

OXA-48-Type-producing Carbapenem-resistant Enterobacteriaceae (CRE) isolates reported to the Centers for Disease Control and Prevention (CDC) as of January 2015, by state



Acinetobacter sp.

Photo credit: CDC

CRE – Identification

Modified Hodge Test

Photo credit: James
Gathany



- CHROM agar
- MacConkey agar plates supplemented with meropenem
- Modified Hodge Test
- Carba NP
- Antibiotic susceptibility disc testing
- PCR
- Check-Direct CPE assay (Check-Points)
- Next-generation sequencing

CRE – Prevention



Photo credit:
Amanda Mills

Improve sanitation procedures and barrier precautions

- Hand hygiene
- Contact precautions
- Disposable equipment
- Environmental cleaning
- Chlorhexidine bathing
- Limit indwelling devices

CRE – Prevention

Implement a surveillance program and communicate outbreaks

- CRE screening of patients that meet pre-specified criteria
- Screen contacts of CRE patients
- Laboratory notification
- Inter-facility communication/identification of CRE patients at admission



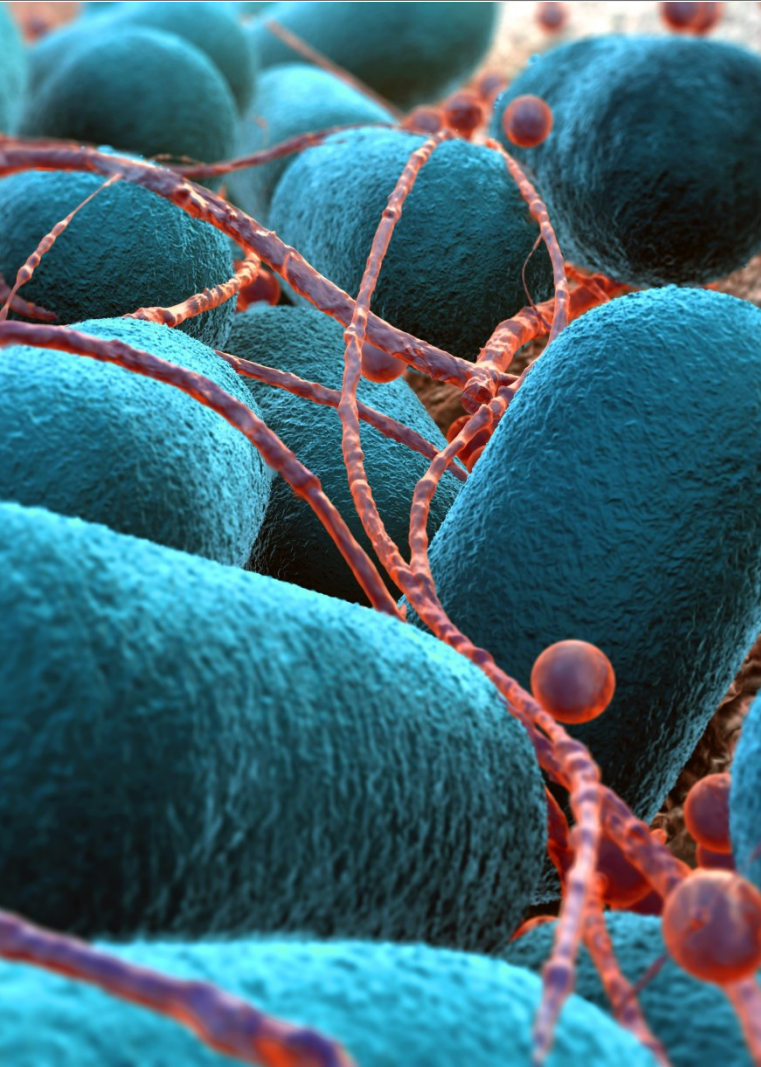
CRE – Prevention



Antimicrobial stewardship

- Promote the appropriate use of antimicrobials
- Appoint a drug expert
- Monitor and report antibiotic prescriptions and resistance patterns
- Educate clinicians about resistance and optimal prescribing

ATCC – Aiding the Scientific Community



ATCC provides top-quality, authenticated reference strains and associated molecular materials

- Enhance diagnostics
- Analyze novel therapeutics
- Improve sterility protocols

ATCC KPC Strains

ATCC® No.	Species	Strain Designation	Presence of Select Virulence Genes
BAA-1705™	<i>Klebsiella pneumoniae</i>	ART 2008133 [D-05, 1338]	<i>bla</i> _{KPC} ⁺ / <i>bla</i> _{NDM} ⁻
BAA-1898™	<i>Klebsiella pneumoniae</i>	-	<i>bla</i> _{KPC} ⁺
BAA-1899™	<i>Klebsiella pneumoniae</i>	-	<i>bla</i> _{KPC} ⁺
BAA-1900™	<i>Klebsiella pneumoniae</i>	-	<i>bla</i> _{KPC} ⁺
BAA-1902™	<i>Klebsiella pneumoniae</i>	-	<i>bla</i> _{KPC} ⁺
BAA-1903™	<i>Klebsiella pneumoniae</i>	-	<i>bla</i> _{KPC} ⁺
BAA-1904™	<i>Klebsiella pneumoniae</i>	-	<i>bla</i> _{KPC} ⁺
BAA-1905™	<i>Klebsiella pneumoniae</i>	-	<i>bla</i> _{KPC} ⁺
BAA-2078™	<i>Klebsiella pneumoniae</i>	-	<i>bla</i> _{KPC} ⁺
BAA-2082™	<i>Enterobacter hormaechei</i>	-	<i>bla</i> _{KPC} ⁺
BAA-2340™	<i>Escherichia coli</i>	1101362	<i>bla</i> _{KPC} ⁺ / <i>bla</i> _{NDM} ⁻
BAA-2341™	<i>Enterobacter cloacae</i>	1101152	<i>bla</i> _{KPC} ⁺ / <i>bla</i> _{NDM} ⁻
BAA-2342™	<i>Klebsiella pneumoniae</i>	1101160	<i>bla</i> _{KPC} ⁺ / <i>bla</i> _{NDM} ⁻
BAA-2343™	<i>Klebsiella pneumoniae</i>	1101172	<i>bla</i> _{KPC} ⁺ / <i>bla</i> _{NDM} ⁻
BAA-2344™	<i>Klebsiella pneumoniae</i>	1101200	<i>bla</i> _{KPC} ⁺ / <i>bla</i> _{NDM} ⁻

KPC Strains Panel
(ATCC® MP-24™)

ATCC NDM Strains

ATCC® No.	Species	Strain Designation	Presence of Select Virulence Genes
BAA-2146™	<i>Klebsiella pneumoniae</i>	1000527, 7561	<i>bla</i> _{NDM} ⁺ / <i>bla</i> _{KPC} ⁻
BAA-2452™	<i>Escherichia coli</i>	NDM-1	<i>bla</i> _{NDM} ⁺ / <i>bla</i> _{KPC} ⁻
BAA-2468™	<i>Enterobacter cloacae</i>	1000654	<i>bla</i> _{NDM} ⁺ / <i>bla</i> _{KPC} ⁻
BAA-2469™	<i>Escherichia coli</i>	1001728	<i>bla</i> _{NDM} ⁺ / <i>bla</i> _{KPC} ⁻
BAA-2470™	<i>Klebsiella pneumoniae</i> subsp. <i>pneumoniae</i>	1002565	<i>bla</i> _{NDM} ⁺ / <i>bla</i> _{KPC} ⁻
BAA-2471™	<i>Escherichia coli</i>	1100101	<i>bla</i> _{NDM} ⁺ / <i>bla</i> _{KPC} ⁻
BAA-2472™	<i>Klebsiella pneumoniae</i> subsp. <i>pneumoniae</i>	1100975	<i>bla</i> _{NDM} ⁺ / <i>bla</i> _{KPC} ⁻
BAA-2473™	<i>Klebsiella pneumoniae</i>	1100770	<i>bla</i> _{NDM} ⁺ / <i>bla</i> _{KPC} ⁻
BAA-2566™	<i>Escherichia coli</i>	--	<i>bla</i> _{NDM} ⁺ / <i>bla</i> _{KPC} ⁻
BAA-2578™	<i>Klebsiella pneumoniae</i>	--	<i>bla</i> _{NDM} ⁺ / <i>bla</i> _{KPC} ⁻

NDM-1 Strains Panel
(ATCC® MP-18™)

ATCC OXA-48 Strains

ATCC® No.	Species	Strain Designation	Relevant Phenotype*
BAA-2523™	<i>Escherichia coli</i>	bMx# 1109131	Produces OXA-48
BAA-2524™	<i>Klebsiella pneumoniae</i> subsp. <i>pneumoniae</i>	bMx# 1103199	Produces OXA-48
BAA-2525™	<i>Providencia rettgeri</i>	bMX# 1103204	Produces OXA-48

*Depositor statement

Antibiotic*	BAA-2523™	BAA-2524™	BAA-2525™
Meropenem	R	R	S
Ertapenem	R	R	I
Imipenem	R	R	R

R = Resistant, S = Susceptible, I = Intermediate susceptibility

*Antibiotic susceptibility determined using E-Test Strips; results may vary depending on the assay and susceptibility cut-offs used

Microbial Strain Authentication



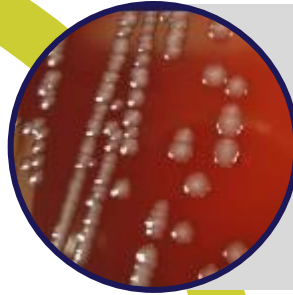
ATCC utilizes both classical and modern techniques

- Phenotypic analysis
- Genotypic analysis
- Functional analysis

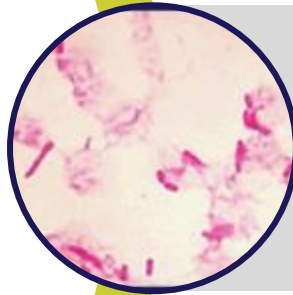
No single method of identification is sufficient

Phenotypic Testing

Culture Purity and
Biochemical
Properties



Colony Morphology



Cell Attributes



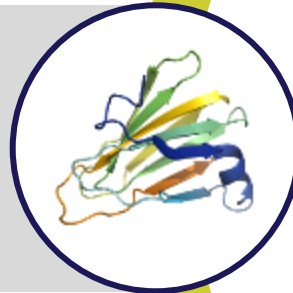
Biochemical Analysis

Genotypic & Protein Testing

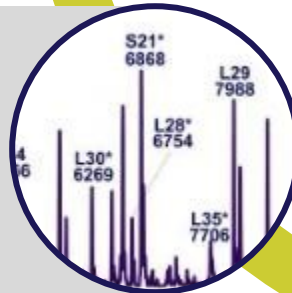
Sequencing



Toxinotyping



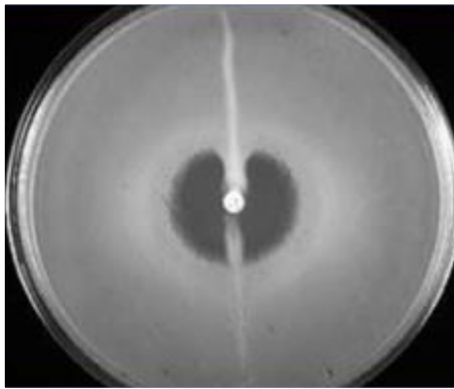
MALDI-TOF



Targeted Gene Sequencing

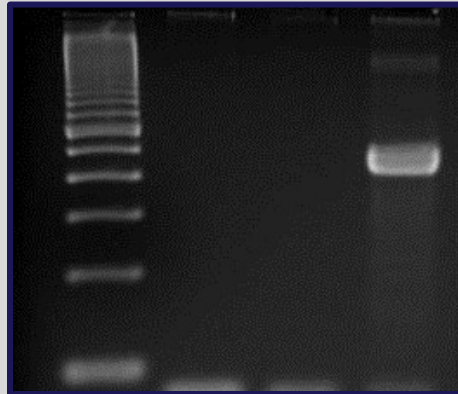
Verification of Drug Resistance

Modified Hodge Test



- Recommended by CLSI and the CDC for the detection of carbapenemase production

Endpoint PCR



- Endpoint PCR used to detect the presence or absence of genes required for antibiotic production

Antibiotic Profiling



- VITEK[®] 2 used to analyze resistance to various antibiotic classes
 - Penicillins
 - Cephalosporins
 - Carbapenems
 - Quinolones
 - Aminoglycosides

Enhance Your Research with ATCC Strains



ATCC – Leading the fight
against superbugs!

- Rapid detection methods
- Innovative therapeutic techniques
- Novel antibiotics
- Updated sterility protocols

Conclusion

- Multidrug-resistant strains are an emerging problem throughout the world
- ATCC acquires, authenticates, and distributes clinically relevant strains that are essential to the scientific community
 - Phenotypic, genotypic, and functional testing
- KPC, NDM, and OXA strains are now available at ATCC
 - Individual strains
 - Microbial panels

Identity
Purity
Authenticity
Homogeneity
Stability
Functionality



Comparability
Quality
Reproducibility
Standardization
Development
Verification



Enterobacteriaceae

Photo credit: CDC

Thank you for joining today!

Register for more ATCC “*Excellence in Research*” webinars, or watch recorded webinars, at www.atcc.org/webinars.



WEBINAR
ONLINE
LEARNING

Please email additional questions to: tech@atcc.org