

Development of a Murine Model of *Pseudomonas aeruginosa* Lung Infection

Presented at, Advancing Animal Models for Antibacterial Drug
Development Workshop

U.S. Food and Drug Administration, White Oak MD
March 5, 2020

Lynn Miesel, Ph.D.
Pharmacology Discovery Services, Ltd.
25, Wugong 6th Road, Wugu District,
New Taipei City, Taiwan 24891

Eurofins Panlabs, Inc.
6 Research Park Drive
St. Charles, MO 63304

Lynn.Miesel@eurofins.com

www.pharmacologydiscoveryservices.com

- Mission: streamline the discovery of therapeutics for *P. aeruginosa*
- *P. aeruginosa* is a leading pathogen that causes VABP and HABP
- Carbapenem resistant *P. aeruginosa* (CRPA) infections have few treatment options
- Related projects
 - Conduct testing services to evaluate therapeutic candidates
 - Developed mouse thigh and lung infection models with MDR isolates
 - Generated PK/PD tutorials with example studies of standard drugs
 - Protocols and example data are available to the drug discovery community

P. aeruginosa Mouse Lung Infection Model

Model features

- Correlate infection mortality with pathogen burden, dissemination and tissue pathology
- Measured endpoints: mortality onset, microbial burden, tissue pathology
- An extended infection period (target ≥ 48 h)
- Compare intranasal (IN) and intratracheal (IT) infection routes
- Conduct with MDR isolates, CDC & FDA AR-BANK

Approach

- Host: persistently neutropenic mice
- Development steps
 - Inoculum optimization
 - Characterize the natural history of infection
 - Benchmark with approved antibiotics



JOURNALS
investing in science

FEMS Pathogens and Disease, 73, 2015, ftv025

doi: 10.1093/femsdp/ftv025

Advance Access Publication Date: 9 April 2015

Research Article

RESEARCH ARTICLE

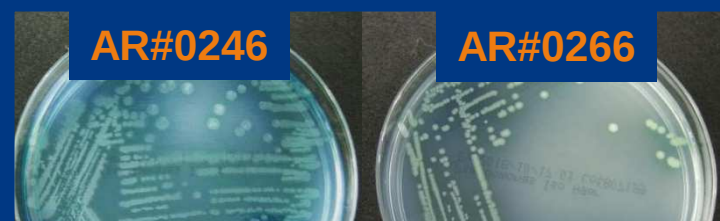
Development and evaluation of murine lung-specific disease models for *Pseudomonas aeruginosa* applicable to therapeutic testing

Matthew B. Lawrenz^{1,2,*}, Ashley E. Biller¹, Daniel E. Cramer¹,
Jennifer L. Kraenzle¹, Julie B. Sotsky¹, Carol D. Vanover¹,
Deborah R. Yoder-Himes³, Angela Pollard⁴ and Jonathan M. Warawa^{1,2,*}

FEMS Pathogens and Disease, 73, 2015 ftv025

P. aeruginosa isolates CDC & FDA AR Isolate Bank

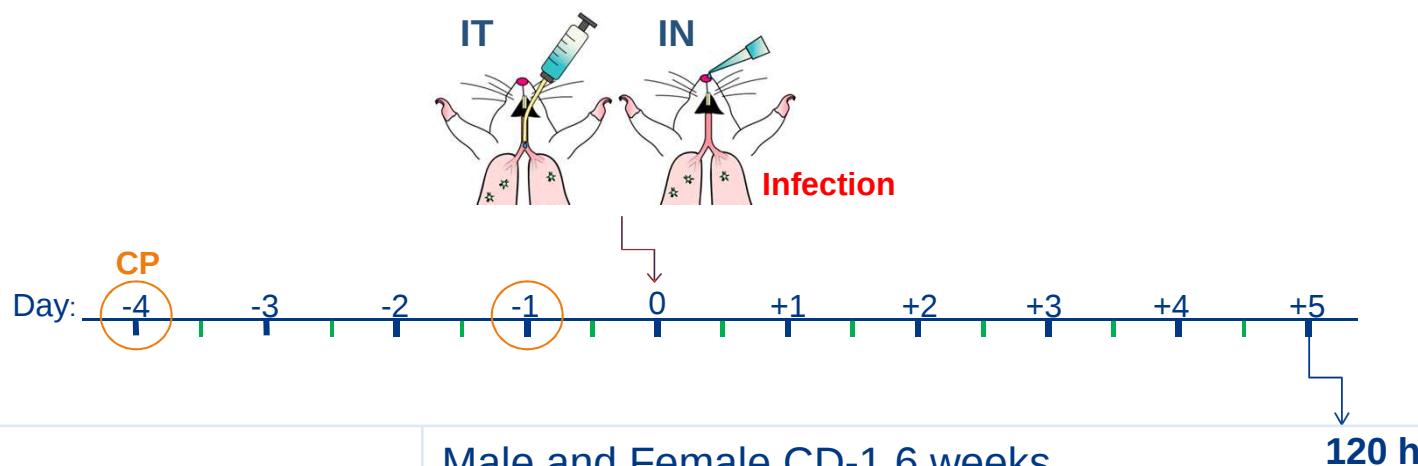
■ Pharmacology
■ ■ Discovery
■ ■ ■ Services



AR-BANK#0246 CRPA, NDM-1		AR-BANK#0266 CRPA
<u>Antibiotic susceptibility</u>		
Notable AMR gene*	NDM-1	None Detected
Carbapenems	R	R
Piperacillin/tazobactam	R	I
Aztreonam	R	S
Ceftazidime/avibactam	CAZ-R, AVYCAZ-R	CAZ-R, AVYCAZ-S
Amikacin	R	S
Ciprofloxacin, Levofloxacin	R	R
Colistin	S	S
<u>Virulence determinants</u>		
Produce pyocyanin and alginate, motile		

*Genomic information from CDC & FDA AR-Bank

Inoculum Optimization



Mice	Male and Female CD-1 6 weeks	120 h
Neutropenia	CP 150 mg/kg D-4, 100 mg/kg D-1	
Inoculum	Titration	
Observations	body temp 6 h intervals, and body weight daily	
Humane endpoints	4°C body temp Δ, 20% weight loss, moribund, and diarrhea	

- Temperature transponders were used for temperature reads

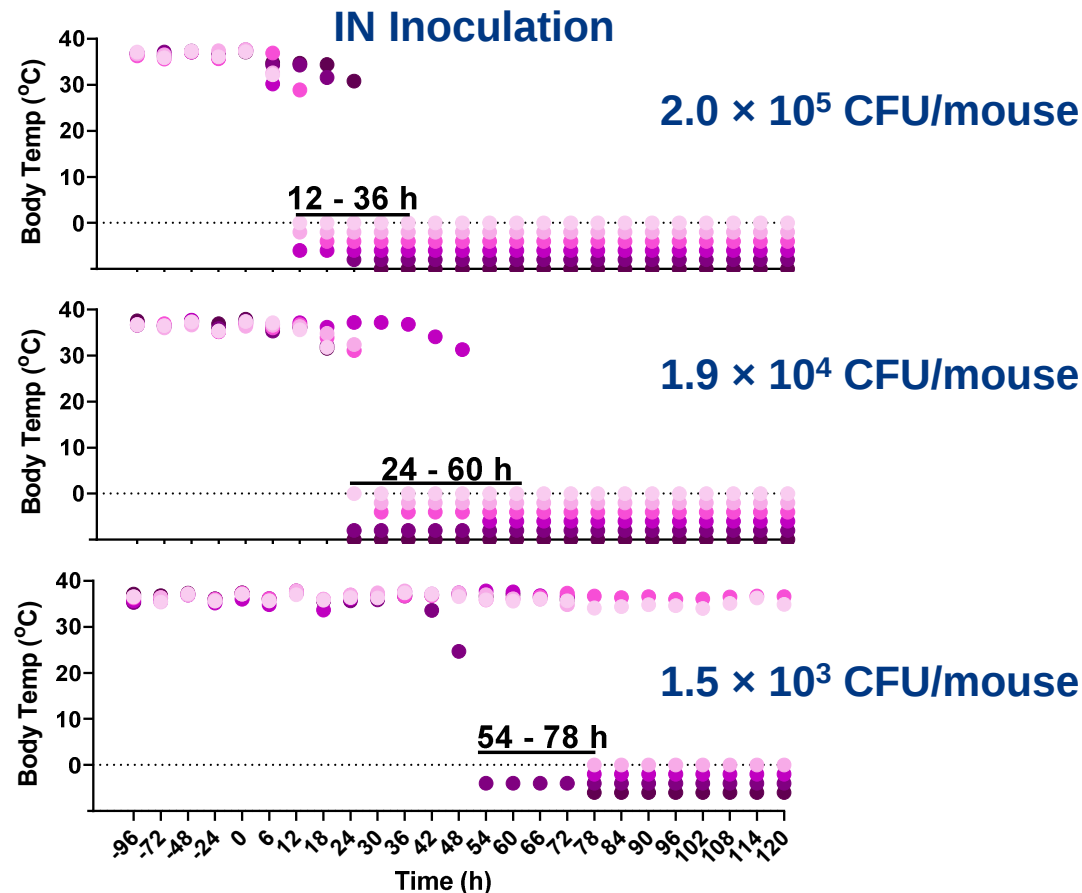


IPTT300 Bio Medic Data Systems

Mouse figure: van Erp EA *et al.* Viruses 2019, 11(6), 508

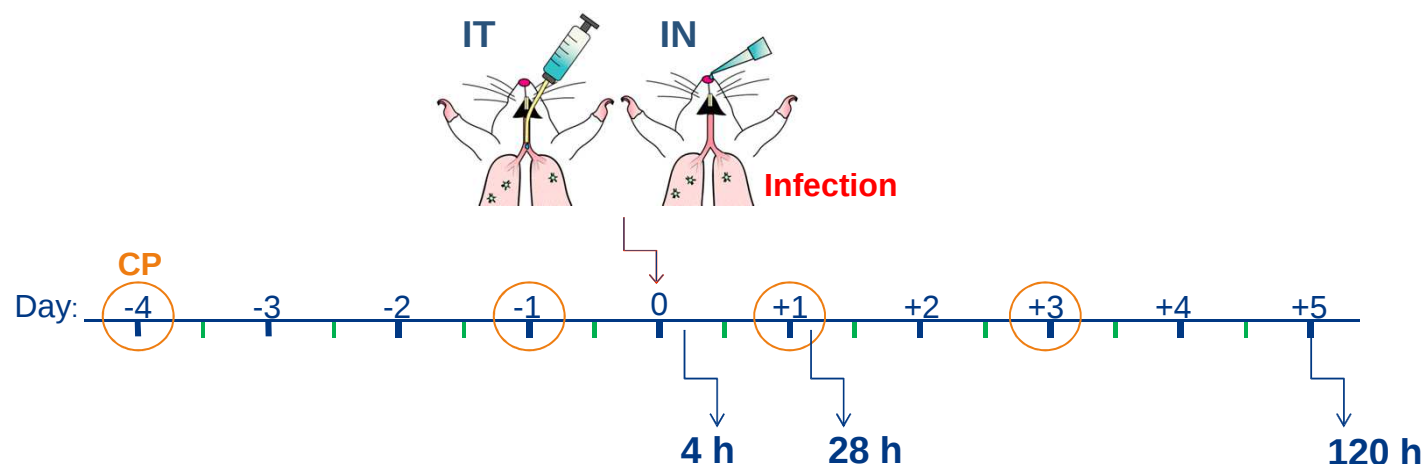
Inoculum Optimization

AR-BANK#0246 NDM-1



- Iterative titration studies were conducted for each strain
- Body temperature telemetry was highly effective for humane terminal endpoints
- Persistent neutropenia was added to subsequent studies to prevent immune recovery
- Male mice only were selected for subsequent studies to minimize variables

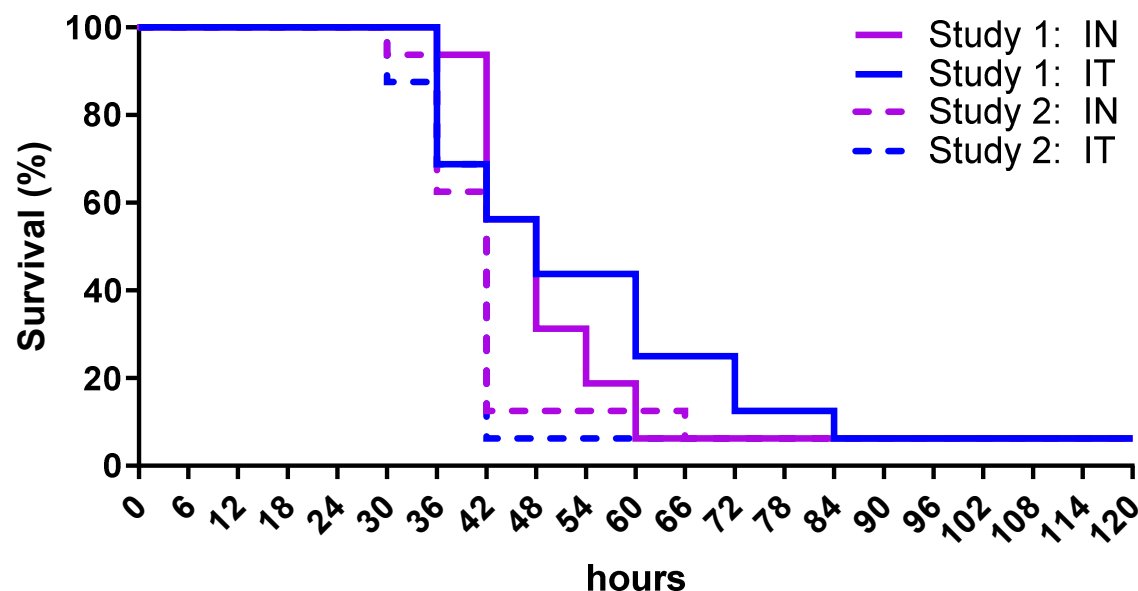
Natural History of Infection



Mice	Male CD-1 6 weeks
Persistent neutropenia	CP 150 mg/kg D-4, 100 mg/kg D-1, D+1, D+3
Target inoculum	#0246 1×10^3 ; #0266 3×10^3
Observations	body temp 6 h, body weight daily
Humane endpoints	4°C body temp Δ , 20% weight loss, moribund, and diarrhea
Scheduled sacrifice	4 h, 28 h, or 5D post infection
Measurements	mortality onset (h) bacterial CFU/g lung and spleen lung gross pathology and histopathology

Natural History of Infection

AR-BANK#0246 NDM-1

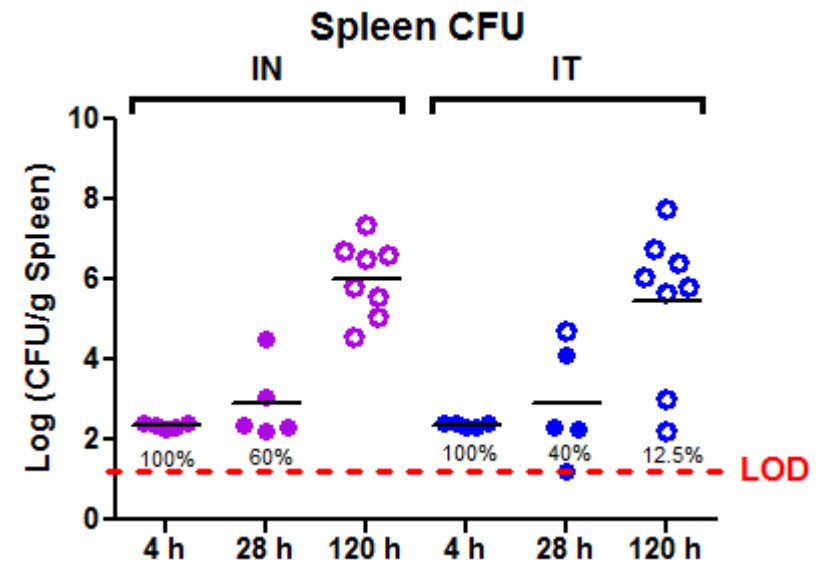
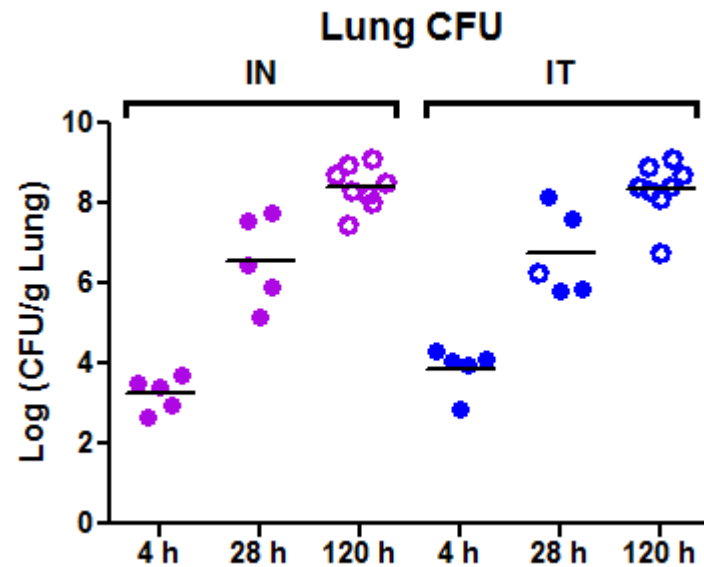


	Study 1	Study 2
Inoculum count (target 1×10^3)	8.8×10^2	1.6×10^3
Mean survival time (h)	IN 46 ± 5 IT 50 ± 6	IN 40 ± 6 IT 39 ± 6

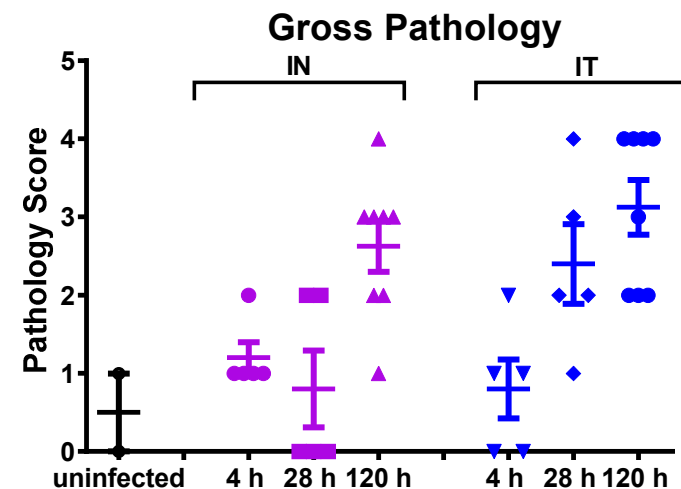
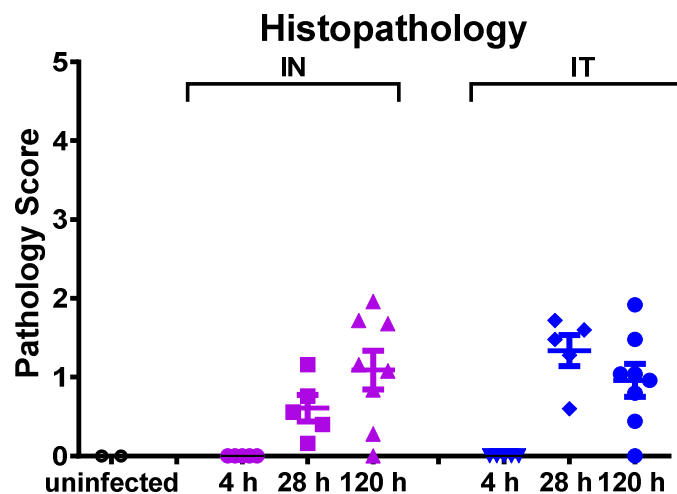
Natural History of Infection

AR-BANK#0246 NDM-1

Pharmacology
Discovery
Services



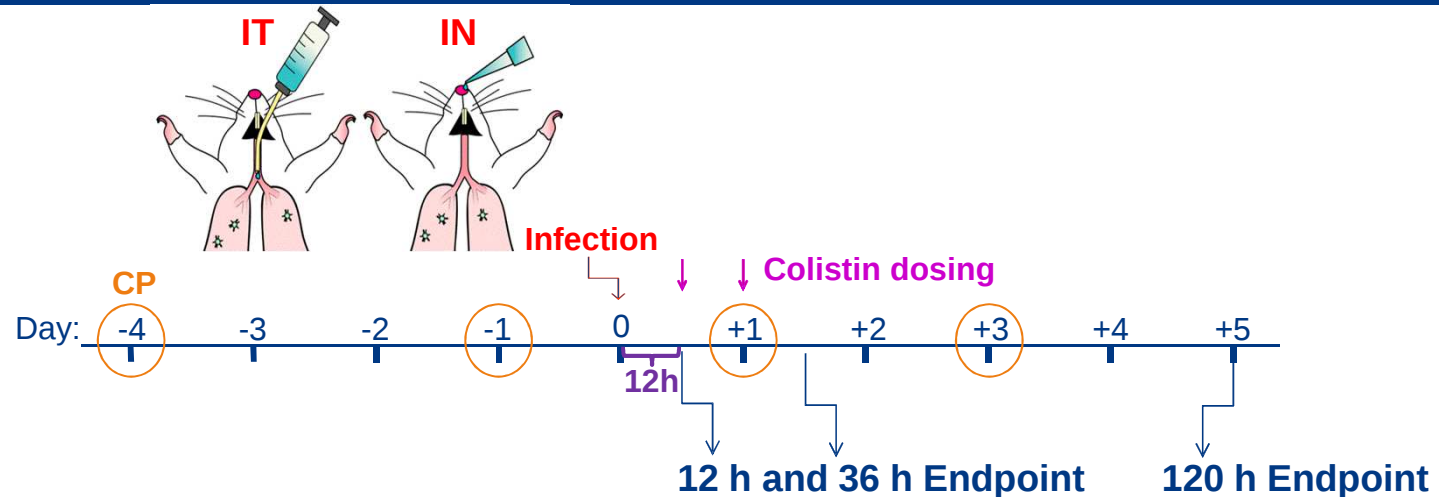
● mice sacrificed at scheduled time points; ○ mice sacrificed at mortality onset



Antibiotic Efficacy

AR-BANK#0246 (NDM-1), Colistin (CST)

Pharmacology
Discovery
Services



Mice:	Male CD-1 6 weeks
Persistent neutropenia	CP 150 mg/kg D-4, 100 mg/kg D-1, D+1, D+3
Target Inoculum	1×10^3 CFU
CST dosing	Start 12 h after infection, 30, 20 and 10 mg/kg BID q12h, one day
Humane endpoints	4°C body temp Δ , 20% weight loss, moribund, diarrhea
Scheduled sacrifice times	12 h, 36 h, or 5 D post infection
Measurements	time to mortality onset bacterial counts in lung and spleen lung gross pathology and histopathology

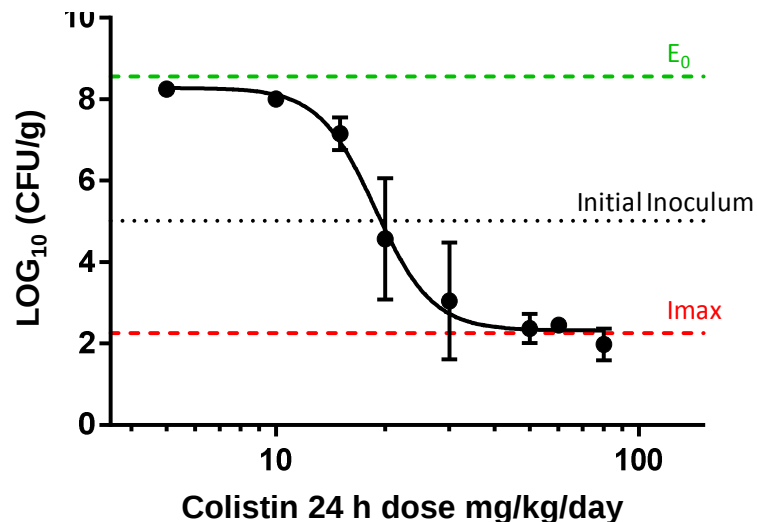
Antibiotic Efficacy

AR-BANK#0246 (NDM-1), Colistin (CST)

Colistin dose selection

- Literature data: Dudhani RF, *et al.* 2010 Antimicrob. Agents Chemother. 54:1117
- Colistin tolerability in persistently neutropenic mice, 48 h 30 mg/kg q12h
- Colistin efficacy, thigh infection model

Colistin q12h BID vs. *P. aeruginosa* AR#0246
Thigh infection model

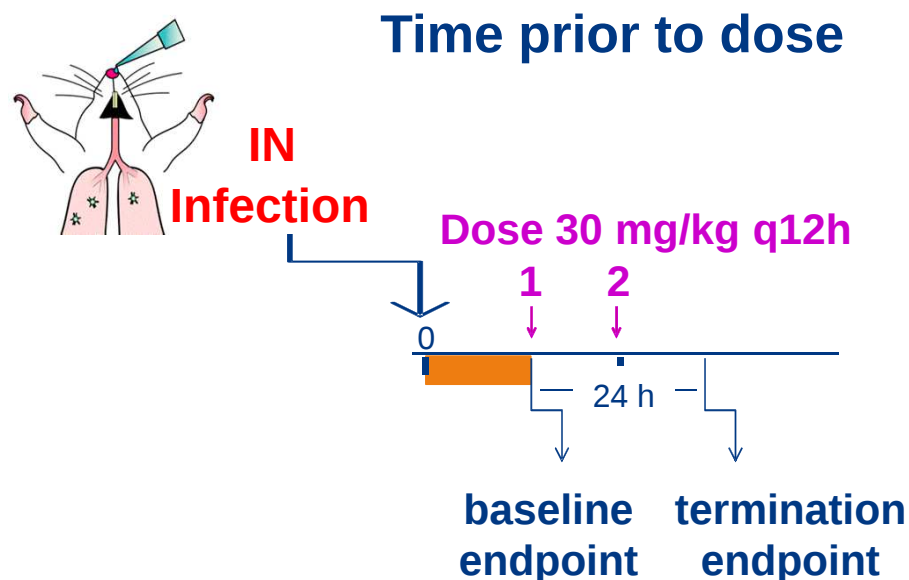


Efficacious dose range:
10 - 40 mg/kg q12h BID

HHSN27200005 NIAID PCMIID A25
Pharmacology Discovery Services
J. Bulitta, University of Florida team

Antibiotic Efficacy

AR-BANK#0246 (NDM-1), Colistin (CST)

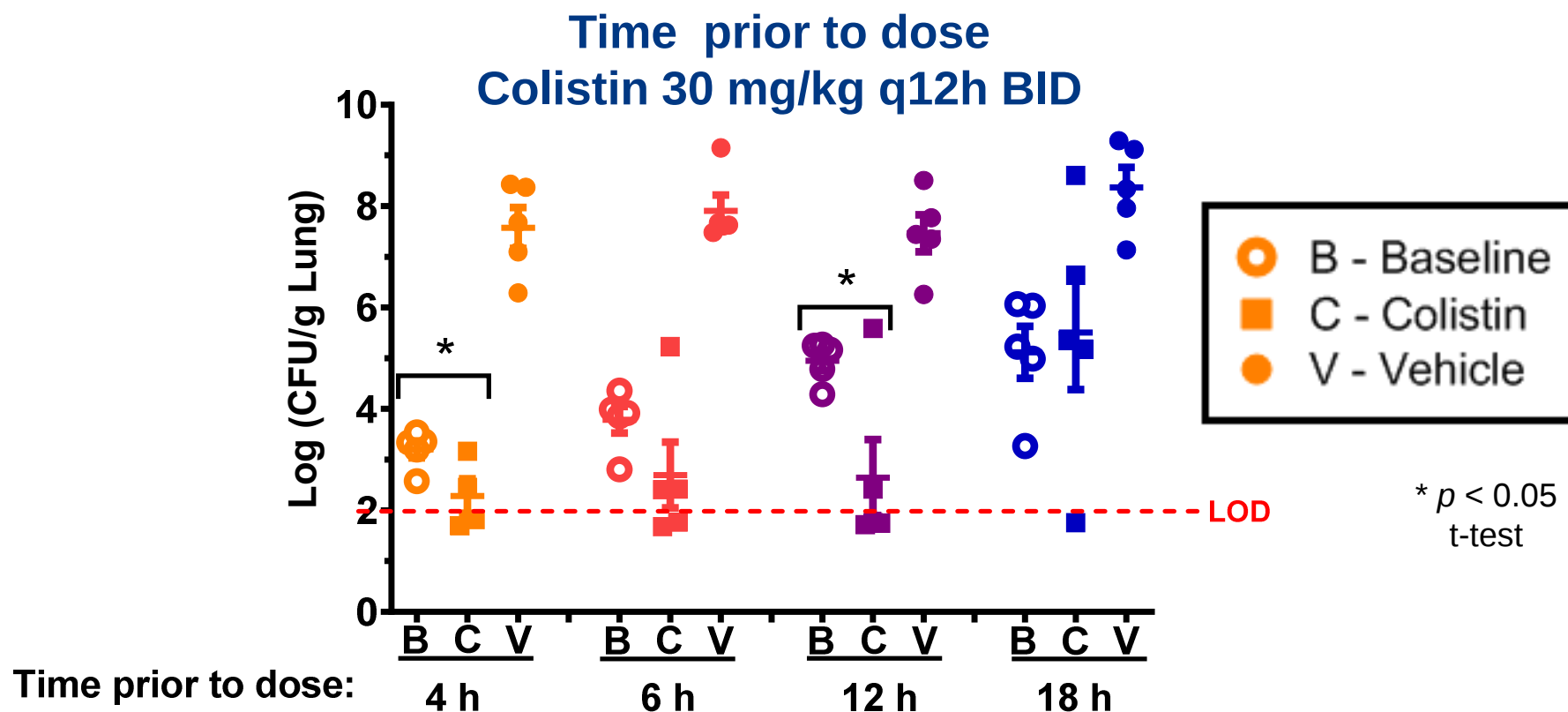


24 h treatment period		
Time of dose 1 ← 12h → Time of dose 2 ← 12h → Termination		
4 h	16 h	28 h
6 h	18 h	30 h
12 h	24 h	36 h
16 h	28 h	40 h

Antibiotic Efficacy

AR-BANK#0246 (NDM-1), Colistin (CST)

Pharmacology
Discovery
Services



The 12 h time to dose was selected:

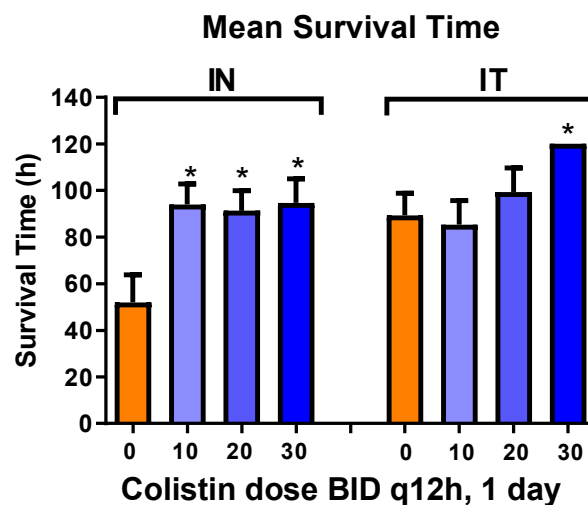
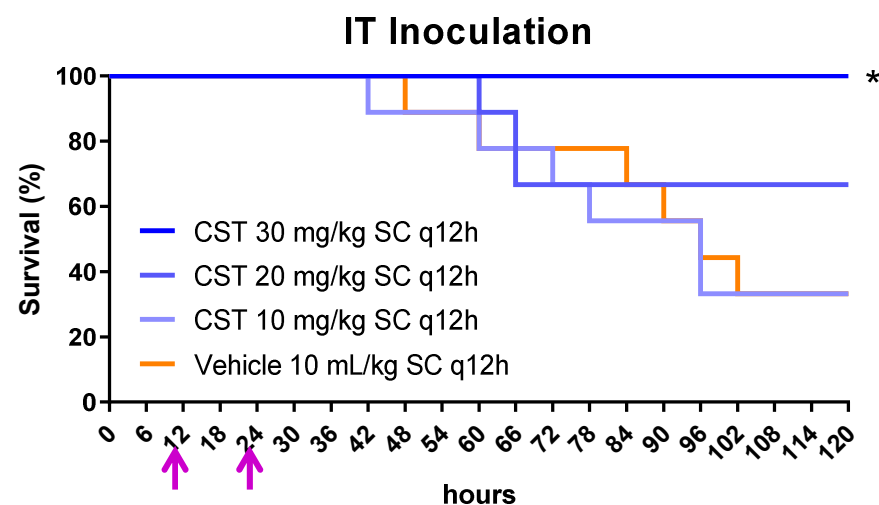
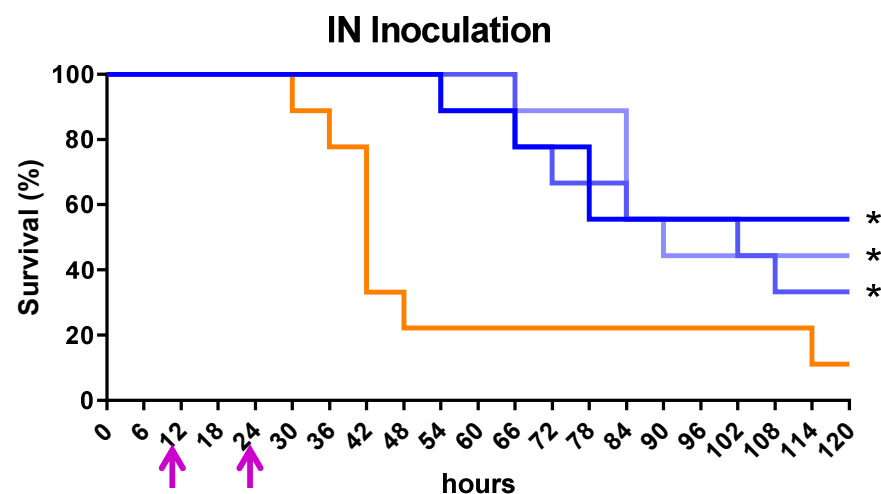
(1) $\sim 10^5$ CFU/g at baseline

(2) Colistin caused a significant reduction in counts relative to baseline

Antibiotic Efficacy

AR-BANK#0246 (NDM-1), Colistin (CST)

Colistin dosage: BID, q12h, 24 h



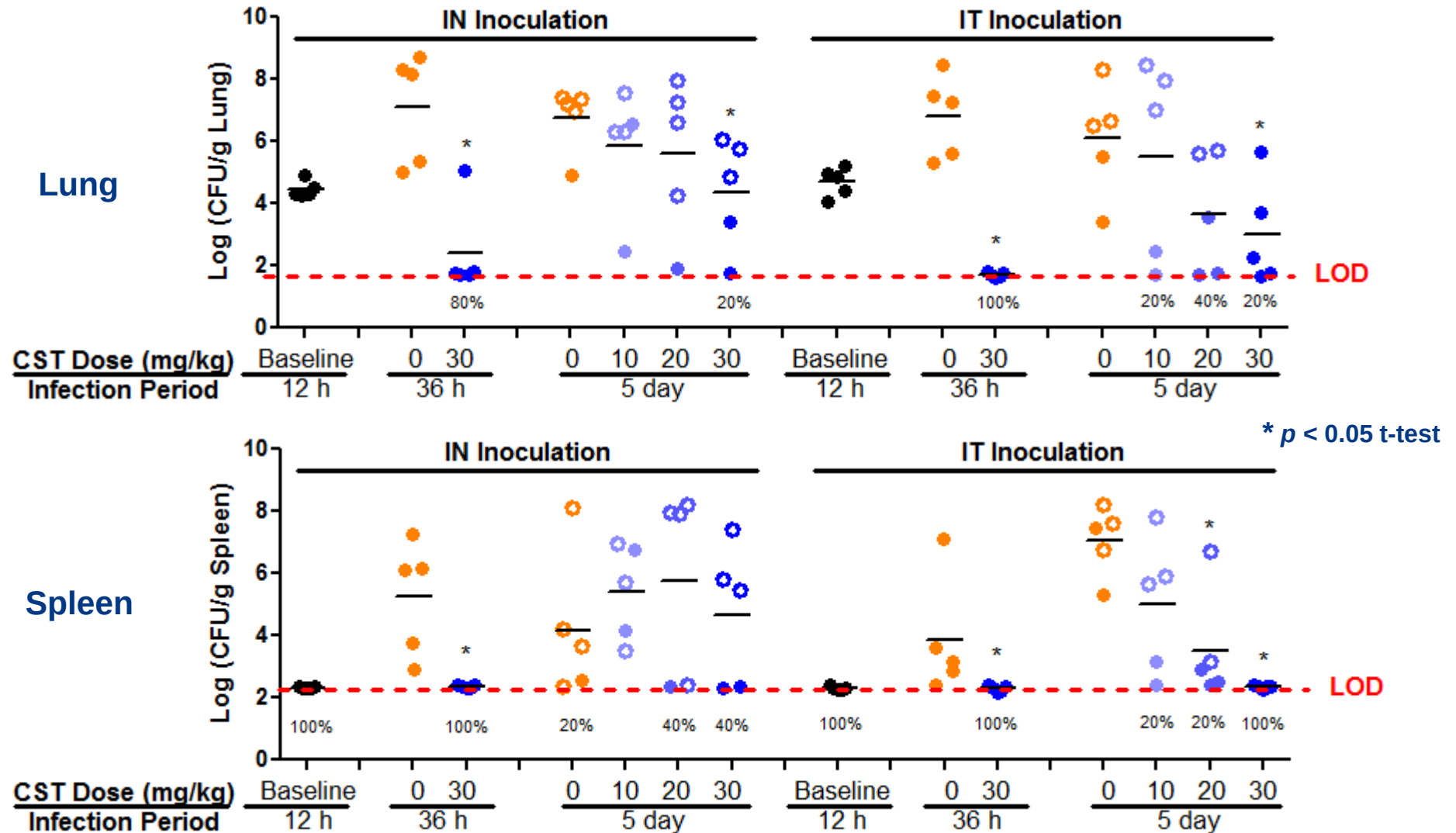
Inoculum count: 1.5×10^3
Kaplan Meier: Fisher's exact test
Mean survival time: On-way ANOVA
* $p < 0.05$

Antibiotic Efficacy

AR-BANK#0246 (NDM-1), Colistin (CST)

Pharmacology
Discovery
Services

Colistin dosage: BID, q12h, 24 h



Summary Across Multiple Studies FDA & CDC AR-BANK#0246 NDM-1

 Pharmacology
  Discovery
   Services

Untreated (vehicle treated) mice at the humane endpoint (or 120 h)

	IN Inoculation	IT Inoculation	IN vs IT
Survival (percent)	3.9%	12%	n.s.
Number of animals	77	41	N/A
Survival time (mean, SEM)	42 ± 3	54 ± 5	*
Lung CFU (Log mean, SEM)	7.7 ± 0.2	7.3 ± 0.4	n.s.
Spleen CFU (Log, mean, SEM)	5.5 ± 0.2	5.8 ± 0.4	n.s.
Gross pathology score (SD)	2.5 ± 1.3	3.2 ± 1.3	n.s.
Histopathology score (SD)	1.0 ± 1.1	1.2 ± 1.3	*
Recommended group size	13		

Pooled data from four studies

* $p < 0.05$

Survival comparison: Fisher's exact test

Pairwise comparisons: t-test

Summary

- Body temperature monitoring facilitates correlation of disease severity with pathogen burden, dissemination, and tissue pathology
- The approach facilitates selection of time points for dosing and sacrifice
- The isolates were highly virulent in neutropenic mice
- Approved antibiotics (colistin and amikacin) were efficacious
- IN and IT infection yielded similar study results
- Limitation, IT infection yielded more variability: survival time and % mortality

Next Steps and Future Directions

- PK/PD to correlate drug exposure with multiple treatment outcomes
- Establish lung infection models with other pathogens
- Establish lung infection models with immune competent mice

Other Resources, NIAID and PDS

- Drug discovery testing services, 75N93019F00131 (A-32)
- Protocols for murine thigh infection models with MDR isolates, HHSN27200003 (A-25)
- PK/PD tutorial PK/PD tutorials with example studies of standard drugs, HHSN27200003 (A-25)
- Protocols and reports are available to the drug discovery community

Acknowledgments

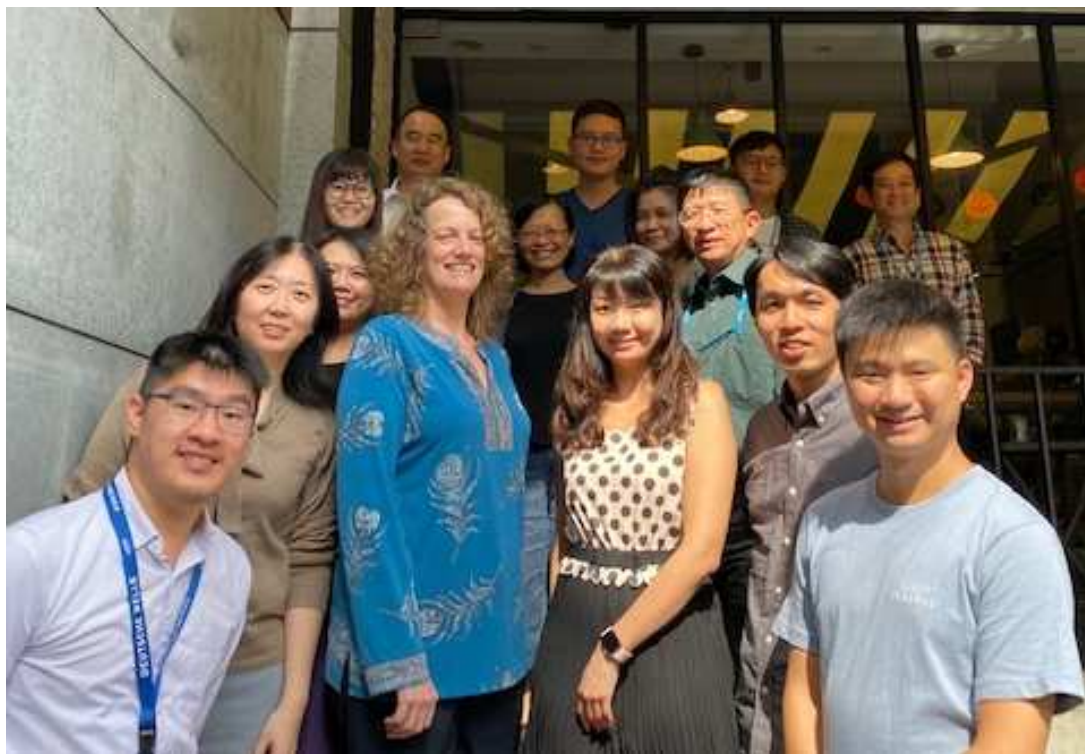
- We thank the CDC & FDA AR Isolate Bank for supplying strains
- This project was funded in whole with Federal funds from the HHS/NIH/NIAID, under Contracts No. HHSN272201700020I / Task Order HHSN27200003 task order A10

Acknowledgments

Name	Organization	Name	Organization
Kun-Yuan Lin	PDST	Chin-Wei Chiang	PDST
Lucy Chia	PDST	David Lo	PDST
Pony Lee	PDST	Hsiao-Lung Chen	PDST
Tim Yeh	PDST	Jiunn-Wang Liao	NCHU
Nien-Chen Lee	PDST	Ray Slay	NIAID
Wei-Ting Liao	PDST	Erin Zeituni	NIAID
Gin-Feng Lai	PDST	Anna Chiang	NIAID
Jui-Che Chien	PDST	Charles Jackson	NIAID
Mei-Lee Hsieh	PDST	Ann Eakin	NIAID
Candance Lee	PDST	Erica Raterman	NIAID
Timmy Huang	PDST	Jane Knisely	NIAID

Come visit us!

■ Pharmacology
■ ■ Discovery
■ ■ ■ Services



www.pharmacologydiscoveryservices.com