DEL RESOURCES

SUPPORTING INFECTIOUS DISEASE RESEARCH

D3403

ABSTRACT

The Biological and Emerging Infections Research Resources Program (BEI Resources) developed by the National Institute of Allergy and Infectious Diseases (NIAID) functions as a centralized bioresource for the accessibility of reference strains of arthropod vectors to the global public health scientific community (<u>www.beiresources.org</u>). The BEI Vector Resources program currently provides over 60 colonies of live mosquitoes as well as Ixodid and soft ticks, sandflies, black flies, and triatomines. These vectors are competent hosts for transmission of blood-borne arboviruses, bacterial, and parasitic pathogens including Dengue, Zika, Powassan viruses, Rocky Mountain Spotted Fever rickettsia, Lyme borreliosis bacteria, Chagas, Leishmania, Babesia, and malaria parasites. Catalog holdings cover different stages of live insects, transgenic strains expressing reporter genes, nucleic acids, plasmids, and cell banks. Methods and protocol manuals for mosquito, tick, and sandfly insectary management and research are available for free download from the BEI Resources website. Part of our community outreach includes webinars to discuss current procedures for the care and maintenance of insect colonies in the laboratory and examples of investigator-led research projects enabled by BEI Vector Resources. This presentation will provide an overview of the biological resources available to the researcher, the benefits of depositing and registering with BEI Resources, and an assessment of publications from investigators utilizing arthropod vectors and associated reagents from the repository. The support from BEI's Vector Resources remains critical in research focused on the elucidation of the mechanisms used by highly adapted vector-borne pathogens to block or subvert host processes and the development of therapies and vaccines.

MISSION OF BEI RESOURCES

- BEI Resources provides NIAID and registered researchers worldwide with a central repository for the acquisition, authentication and distribution of a broad range of unique and quality assured pathogen research materials that will aid in the development and evaluation of vaccines, therapeutics, and diagnostics (Fig 1).
- BEI Resources covers NIAID Category A, B and C priority pathogens, including emerging infectious disease (EID) pathogens, arthropod vectors and reagents.
- Available arthropod Vector Resources materials include mosquitoes, sand flies, ticks, and triatomines, species known to be competent for transmission of viral, parasitic, and bacterial EID pathogens.
- Investigators can register at <u>www.beiresources.org</u> for free access to research materials at no cost. Institutional review and approval is required for MTA and for required material biosafety level classification.

Figure 1. BEI Resources procures, expands, authenticates and distributes research strains and reagents for investigators conducting pathogen and vector research.



BEI RESOURCES

10801 University Boulevard, Manassas, Virginia 20110-2209 **Phone:** 800.359.7370

BEI Resources is funded by the National Institute of Allergy and Infectious Diseases (NIAID) and managed by ATCC. Materials are distributed only to established qualified research laboratories with facilities and safety programs appropriate for the level of material requested. © 2019 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.

The Role of BEI Resources in the Support of Arthropod **Vector Research**

Robert E. Molestina, Rebecca Bradford, and Sujatha Rashid

BEI Resources, American Type Culture Collection, Manassas, VA 20110

ARTHROPOD VECTOR RESOURCES

Vector	Species Represented
Mosquitoes	 Aedes aegypti Aedes albopictus Anopheles albimanus Anopheles arabiensis Anopheles arabiensis Anopheles atroparvus Anopheles coluzzii Ngousso Anopheles dirus Anopheles farauti Anopheles freeborni Anopheles funestus Anopheles gambiae Anopheles merus Anopheles merus Anopheles quadriannulatus Anopheles stephensi Culex quinquefasciatus Culex tarsalis
Sand Flies	 Lutzomyia longipalpis Phlebotomus papatasi Phlebotomus duboscqi Phlebotomus sergenti Phlebotomus arabicus
Ticks	 Amblyomma americanum Dermacentor variabilis Haemaphysalis longicornis Ixodes pacificus Ixodes ricinus Ixodes scapularis Rhipicephalus sanguineus
Reduviids	 Rhodnius prolixus
Black flies	Simulium vittatum

Research Resources for Arthropod Vector Biology

• BEI Resources provides free of charge insectary protocol manuals for mosquitoes, ticks, fleas and sand flies, containing an array of insectary management and experimental approaches for Arthropod Vector Biology. www.beiresources.org/Catalog/

VectorResources.aspx





• Larvae (L3) / Pupae Adult



- Larvae
- Nymph
- Adult





- Adult
- Eggs
- Larvae
- Adult

EXAMPLES OF ARTHROPOD VECTOR COMPETENCE

BEI Vector	
 Culex quinquefasciatus (Southern house mosquito) Culex tarsalis YOLO (Meetern encenhalitie) 	St. I <i>Wud</i> WE
 (Western encephalitis mosquito) <i>Rhodnius prolixus</i> (Assassin bug) 	Тгур
 Phlebotomus papatasi PPNS (Sand fly; Egypt) 	<i>Leis</i> Nap
 Lutzomyia longipalpis LLJB (Sand fly, Brazil) 	Leis
 Ixodes scapularis (deer / black legged tick) 	<i>Bori</i> age

VECTOR RESOURCES CITATIONS 2013-2023



SUMMARY

- resource availability.
- depositor recognition.

ACKNOWLEDGEMENTS

nor by the U.S. Government.

© ATCC 2019. The ATCC trademark, trade name, any, and all ATCC catalog numbers listed in this publication are trademarks of the American Type Culture Collection unless indicated otherwise.



Pathogen Transmission Competence

Louis encephalitis (SLEV), West Nile (WNV), Rift Valley Fever (RVFV), chereria bancrofti (filariasis).

EV, SLEV, WNV, RVFV, CEV, avian *Plasmodium* spp.

panosoma cruzi (Chagas' Disease)



relia burgdorferi, Anaplasma phagocytophilum, Ehrlichia muris-like agent, Powassan virus, Babesia spp.

Search terms performed in Google Scholar

Mosquitos: Aedes OR Anopheles OR Culex AND BEI Resources Ticks: Amblyomma OR Dermacentor OR Ixodes OR Rhipicephalus **AND BEI Resources Reduviids:** Rhodnius OR Triatomine OR Reduviid AND BEI Resources Sand flies: Lutzomyia OR Phlebotomus AND BEI Resources Black flies: Simulium AND BEI Resources Reduviids Sandflies Black flies

• BEI Resources provides vector and pathogen research materials for investigators worldwide. Metrics on program usage are critical for assessment and continued

• As a program requirement, literature citations to BEI Resources materials define arthropod vector strain origin and provide

• Program usage based on research publications indicate that materials are used for vector biology and genetics research, pathogen transmission models, and vector control strategies.



BEI Resources thanks Ellen Dotson and Michael Levin of the Centers for Disease Control and Prevention, Tobin Roland of Walter Reed Army Institute for Research, and Elmer W. Gray of University of Georgia for providing resources and expertise for deposited materials. Vector images courtesy of James Gathany, Public Health Image Library (PHIL), CDC.

BEI Resources is funded under contract 75N93024C00001 by the National Institute of Allergy and Infectious Diseases, National Institutes of Health, Department of Health and Human Services. The views expressed herein neither imply review nor endorsement by HHS