

## Q&A ATCC® *Excellence in Research* Webinar “Enhancing Vector-borne Research with Biological and Molecular Standards”

### General Questions

1. Will we be able to download the presentation?  
This presentation will be available to watch on demand on the ATCC website.
2. What primer sets can I use with the Synthetic Dengue RNA?  
The ATCC Synthetic Dengue RNA was designed and developed to contain key target regions of the viral genome as well as sequences complementary to primers used in the CDC DENV 1-4 Real-Time PCR Assay for the detection and serotype identification of dengue virus as well as the PLoS Neglected Tropical Diseases 2013 publication by Waggoner *et al.*, entitled Single-Reaction, Multiplex, Real-Time RT-PCR for the Detection, Quantitation, and Serotyping of Dengue Viruses.
3. ATCC recommends that the synthetic West Nile virus RNA be stored at 8°C or colder, but ships them on dry ice. Are there any stability problems if these are stored at 8°C rather than -80°C?  
Accelerated stability studies at 45°C have been performed on this sample, and no degradation was observed.
4. I’m trying to culture *Trypanosoma*. Is ATCC medium 431 suitable for culturing procyclic forms and mammal-infective forms?  
The various media, including ATCC medium 431, listed on the ATCC website for cultivation of the different strains of *Trypanosoma* are all suitable for cultivation of the insect-infective forms (hence the 25°C recommended incubation temperature).  
  
ATCC does not culture the blood-dwelling trypomastigote forms. You may be able to find useful information on the cultivation of those forms in the publication: Taylor AE and Baker JR, *In Vitro Methods for Parasite Cultivation*. Academic Press, 1987. Also, please download a free copy of the ATCC Protistology Culture Guide from our website at [www.atcc.org/guides](http://www.atcc.org/guides) for more information on how to culture protists.
5. Does ATCC distribute mosquitoes, antibodies, or other resources for malaria research?  
ATCC may distribute some items which may be useful for malaria research, such as various *Plasmodium* species and associated DNA products.

For more resources and information regarding malaria research, please visit the Malaria Research and Reference Reagent Resource Center (MR4) online at [www.mr4.org](http://www.mr4.org). The Malaria Research and Reference Reagent Resource Center (MR4) was established to provide a central source of quality controlled malaria-related reagents and information to the international malaria research

community. Materials available to qualified, registered users include parasites, mosquitoes, vectors, antibodies, antigens, gene libraries, molecular probes, and constructs.

6. Where can I find other arthropod vectors for vector-borne research?

The Vector Resource center within BEI Resources currently supplies mosquitoes, ticks, reduviids, sand flies, and snails to qualified researchers. Please visit BEI Resources online at [www.beiresources.org](http://www.beiresources.org).