



## Q&A ATCC® *Excellence in Research* Webinar “Seeing is Believing – Reporter-Labeled Microbial Control Strains”

### General Questions

1. Will we be able to download the presentation?

This presentation will be available to watch on demand on the ATCC website, or [click here](#).

2. Are there licensing requirements to consider when choosing between GFP- and NanoLuc®-labeled strains?

GFP products are distributed under the terms of the ATCC MTA and a Limited Use License from Life Technologies. The commercial use of GFP and research use of GFP by for-profit institutions both require a license from Life Technologies. Many large companies already have a license in place. Academic and government institutions are generally exempt from this requirement. The Limited Use License is available on the product pages of our website.

NanoLuc® and Nano-Glo® products are distributed under the terms of the ATCC MTA and a Limited Use License from Promega. Promega does not require for-profit institutions and academic and government institutions to have a license for the research use of NanoLuc® and Nano-Glo®. The commercial use of NanoLuc® and Nano-Glo® requires a license from Promega. The Limited Use License is available on the product pages of our website.

3. What are the storage conditions for the NanoLuc® substrate?

The labeled strains will be shipped to you on dry ice with the substrate and lysis buffer. Store the lysis buffer and the substrate separately at -20°C. The strains should be stored at -70°C or colder. Make the reaction solution by mixing the substrate and the lysis buffer. The reaction solution will keep for up to a week at 4°C, but will slowly lose activity. For the brightest response, prepare only as much fresh reaction solution as required for each experiment, storing the substrate and lysis buffer separately until needed.

4. How long is the NanoLuc® label visible?

The reporter will produce an intense blue glow-response signal until the substrate runs out or the reaction dries. The light is generally visible for several minutes.

© 2015 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. NanoLuc® and Nano-Glo® are registered trademarks of Promega Corporation.