

# Validation data for A549-RepTor™ cells

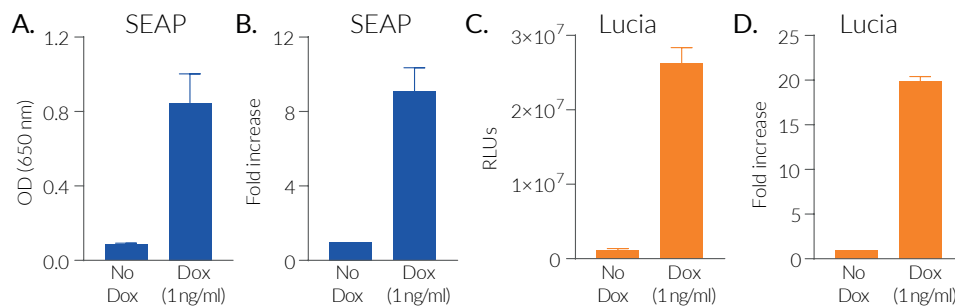
<https://www.invivogen.com/tet-on-a549-reptor-cells>

For research use only

Version 24A25-NJ

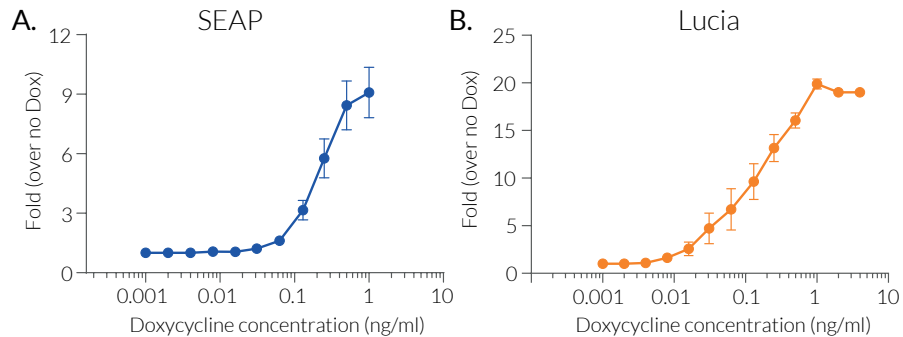
A549-RepTor™ cells are designed for the Tet-on inducible expression of a protein. The absence of expression leakage has been verified in A549-RepTor™ cells transfected with pTiGer plasmids coding for the SEAP (secreted embryonic alkaline phosphatase) or Lucia luciferase reporter proteins (Figure 1). SEAP and Lucia are conditionally expressed by A549-RepTor™ SEAP and A549-RepTor™ Lucia cells upon incubation with Doxycycline (Dox), in a dose-dependent manner (Figure 2).

## No expression leakage of protein of interest in A549-RepTor™ cells



**Figure 1: Expression of SEAP or Lucia reporter proteins in transfected A549-RepTor™ cells.** A549-RepTor™ cells were transfected with the gene coding for SEAP (secreted embryonic alkaline phosphatase) or Lucia luciferase cloned into a pTiGer plasmid. The cells were then treated or not with Doxycycline (Dox) at 1 ng/ml for 24 hours. (A, B) The SEAP activity in the supernatant of A549-RepTor™ SEAP cells was assessed using QUANTI-Blue™ detection reagent. The data is shown as (A) OD at 650 nm and (B) fold increase (mean + SEM). (C, D) The Lucia activity in the supernatant of A549-RepTor™ Lucia cells was assessed using QUANTI-Luc™ 4 Lucia/Gaussia detection reagent. The data is shown as (C) relative light units (RLUs) and (D) fold increase (mean + SEM).

## Inducible SEAP or Lucia expression in A549-RepTor™ cells



**Figure 2: Dose-dependent Doxycycline-mediated expression of SEAP or Lucia in A549-RepTor™ cells.** A549-RepTor™ cells were transfected or not with the gene coding for SEAP (secreted embryonic alkaline phosphatase) or Lucia luciferase cloned into a pTiGer plasmid. A549-RepTor™ SEAP cells and A549-RepTor™ Lucia cells were incubated with increasing concentrations of Doxycycline. After 24 hours, the SEAP or Lucia activity in the supernatant was assessed using (A) QUANTI-Blue™ or (B) QUANTI-Luc™ 4 Lucia/Gaussia, respectively. The data is shown as fold induction over no Dox treatment (mean + SEM).

### TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873  
InvivoGen USA (International): +1 (858) 457-5873  
InvivoGen Europe: +33 (0) 5-62-71-69-39  
InvivoGen Asia: +852 3622-3480  
E-mail: [info@invivogen.com](mailto:info@invivogen.com)



Any questions about our cell lines?  
Visit our FAQ page.

 **InvivoGen**  
[www.invivogen.com](http://www.invivogen.com)